

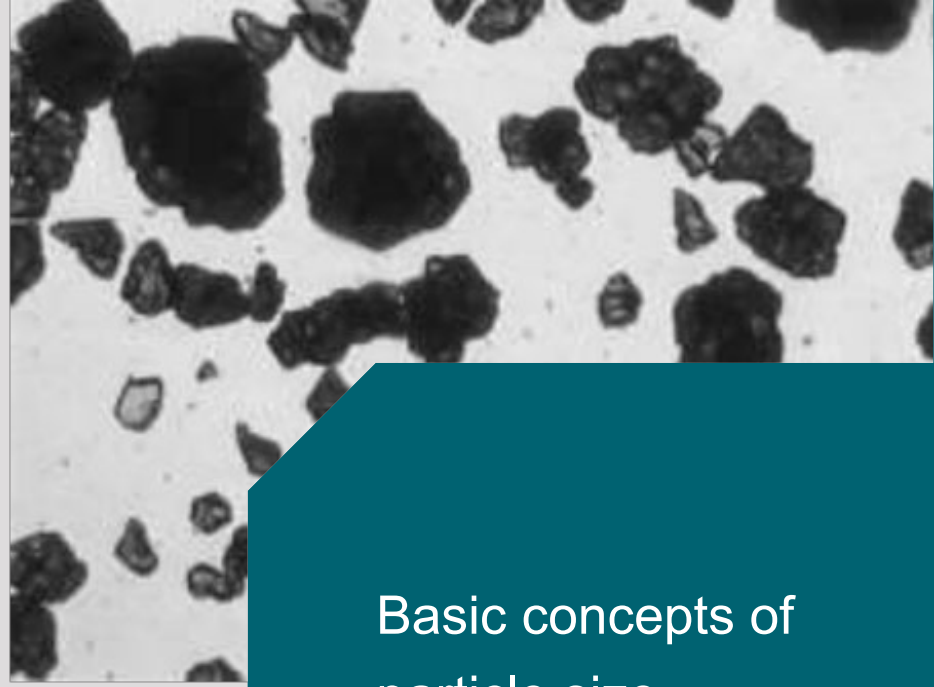
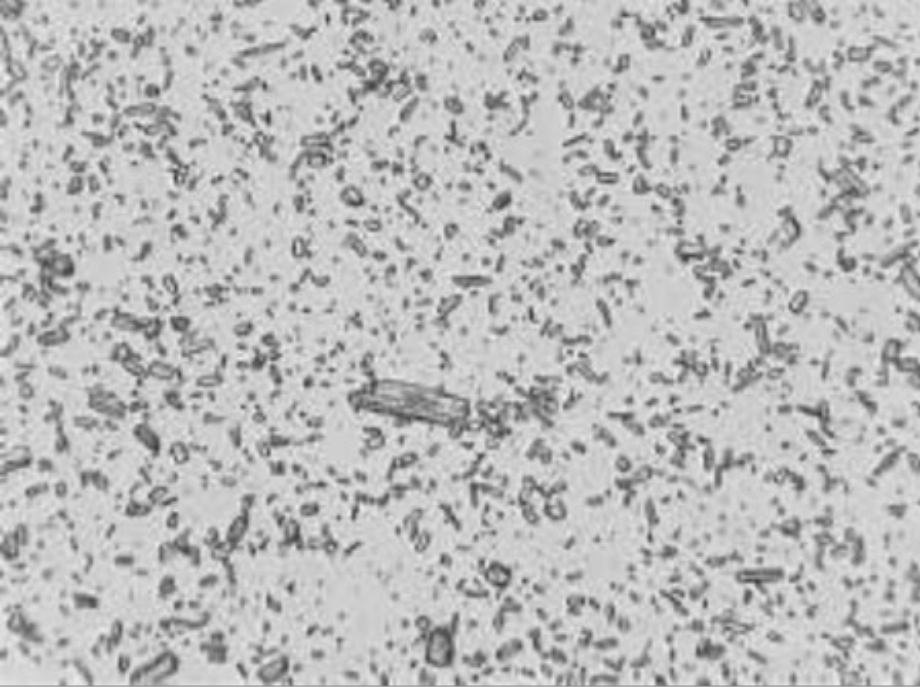


**Malvern
Panalytical**

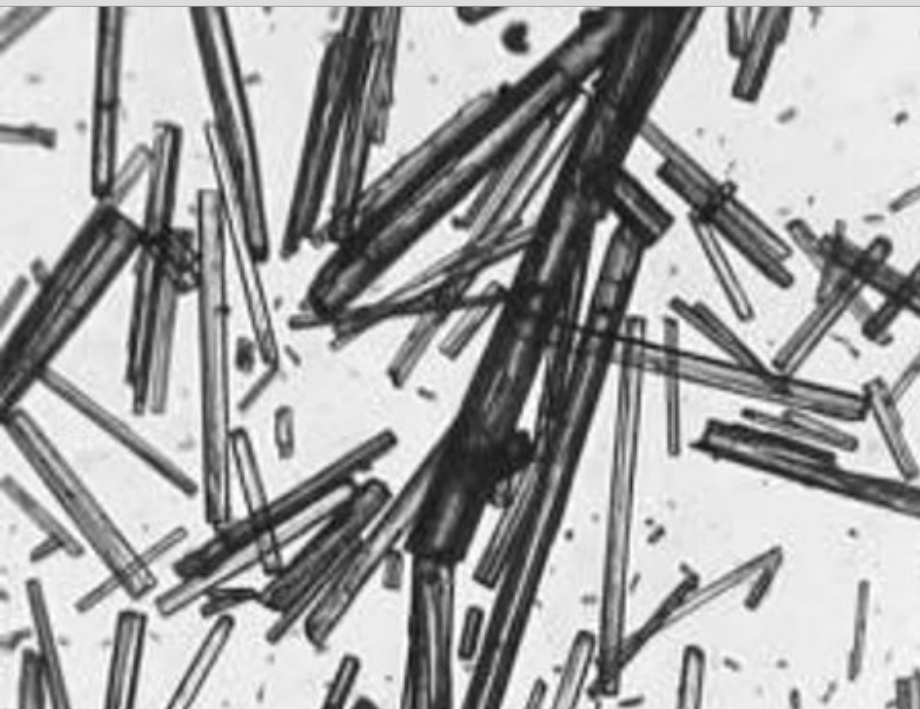
Mastersizer 3000 Customer Training Course

Part 1: Basic Principles and Data Quality



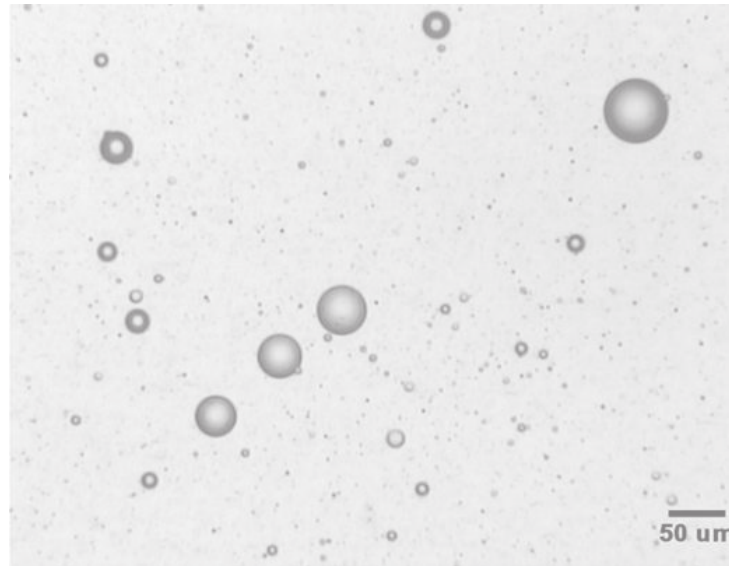


Basic concepts of
particle size



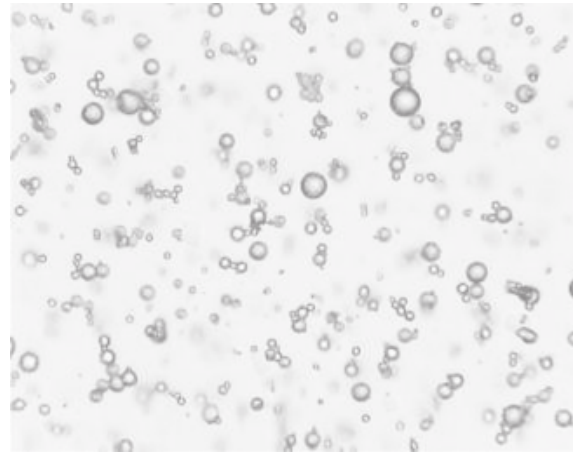
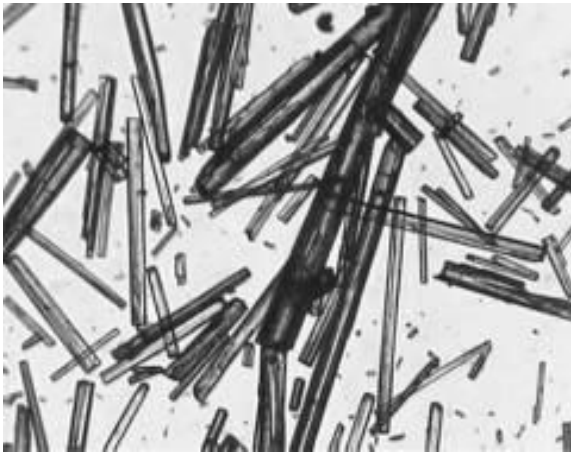
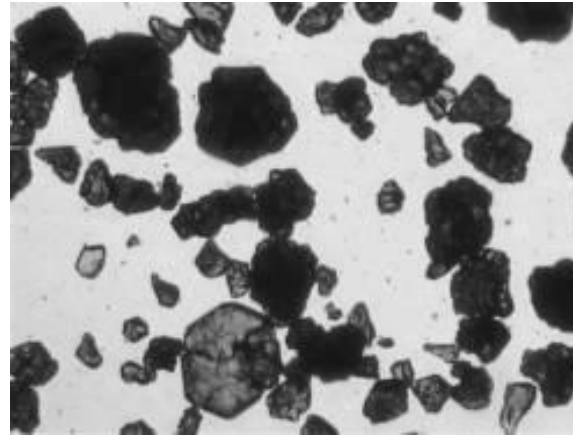
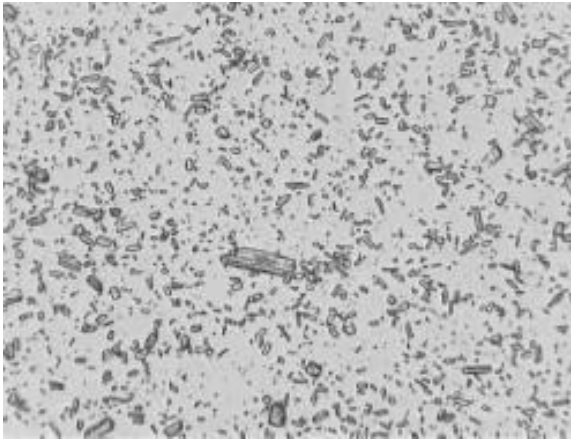
What do we mean by a particle?

- A particle can be described as a discrete sub-portion of a substance, e.g.
 - solid particles
 - gas bubbles
 - or liquid droplets



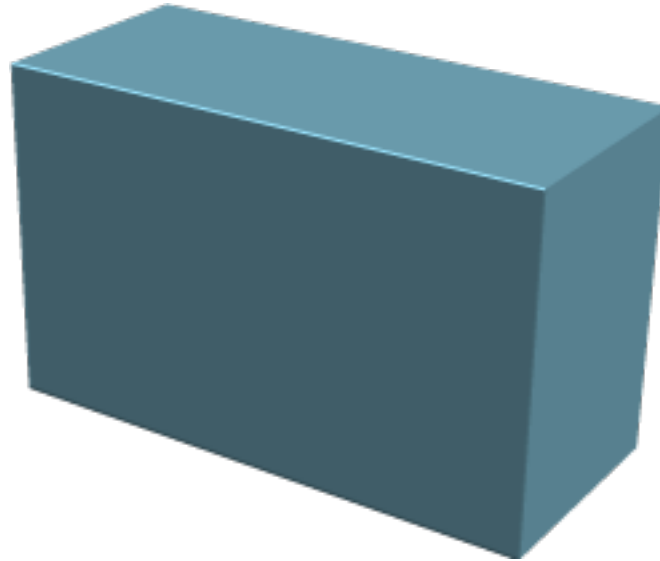
- Laser diffraction measures particles in the size range from nanometres to millimetres

Particles come in many different shapes (as well as sizes)



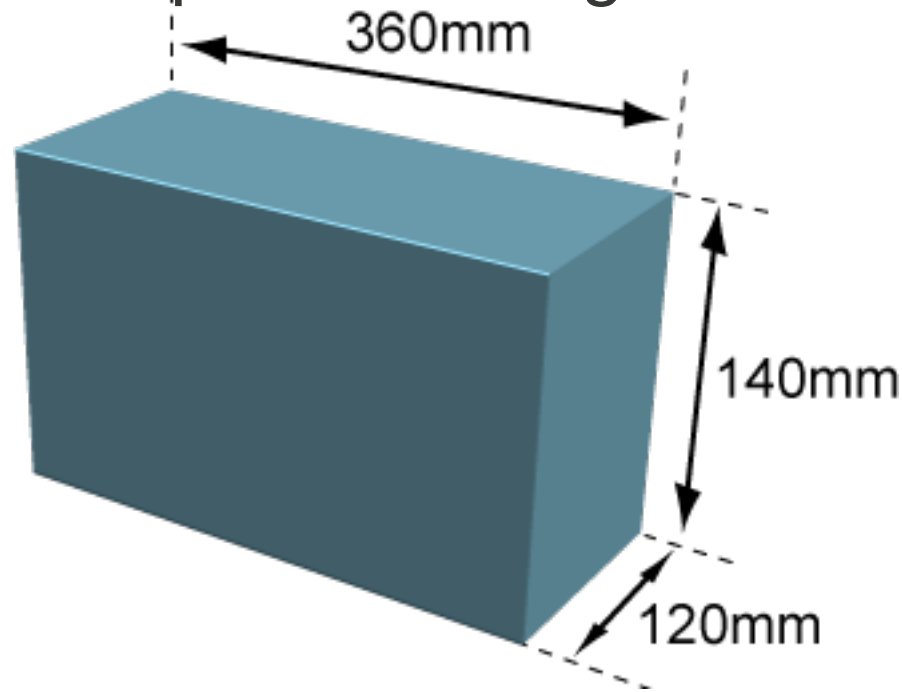
- How do we describe the size of these particles?

Basic concepts of particle sizing



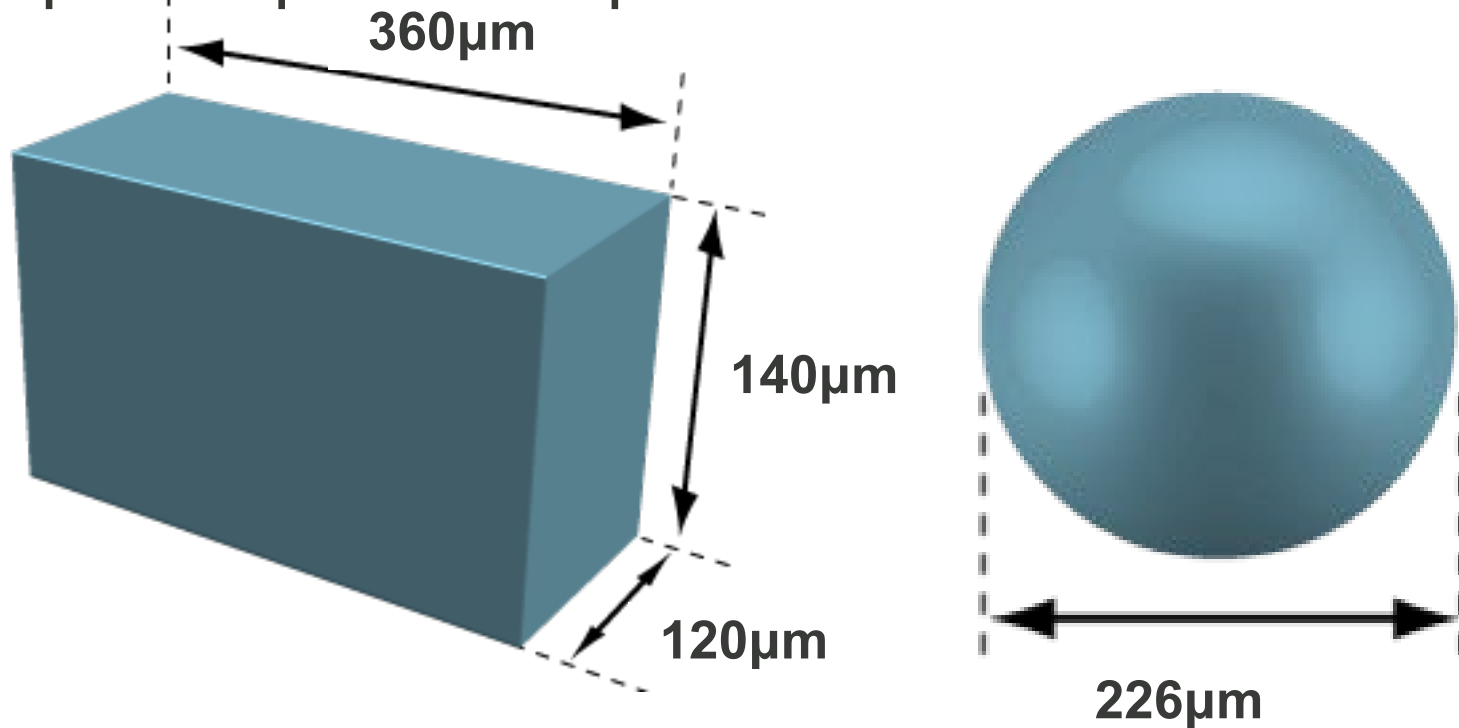
- You are given a regular-shaped object and a ruler and asked to give a one-number indication of its size
- What would your reply be ?

Basic concepts of particle sizing



- You may reply: “360x140x120mm”
 - Which might be correct but it is not **one** number.
 - It is not possible to describe the size of this 3-dimensional object with a single number

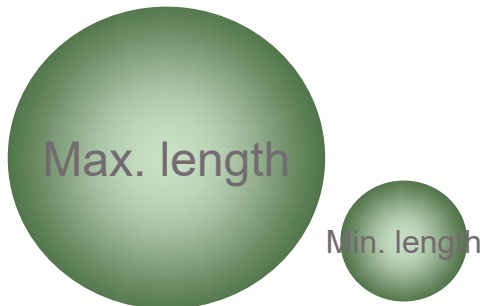
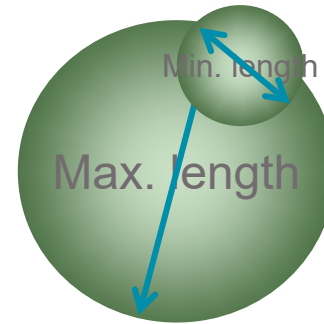
Concept of equivalent spherical diameters



- The rectangular box has the same volume as a sphere of 226 μm diameter.
 - The volume equivalent spherical diameter is 226 μm

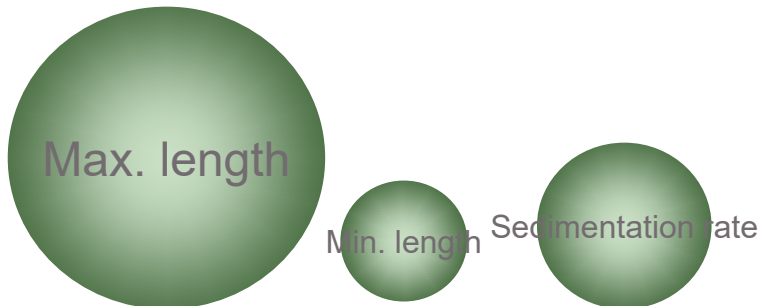
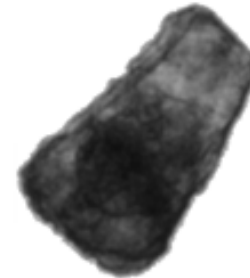
How do we describe the size of these particles

- Equivalent spheres
 - Maximum length
 - Minimum length



How do we describe the size of these particles

- Equivalent spheres
 - Maximum length
 - Minimum length
 - Sedimentation rate



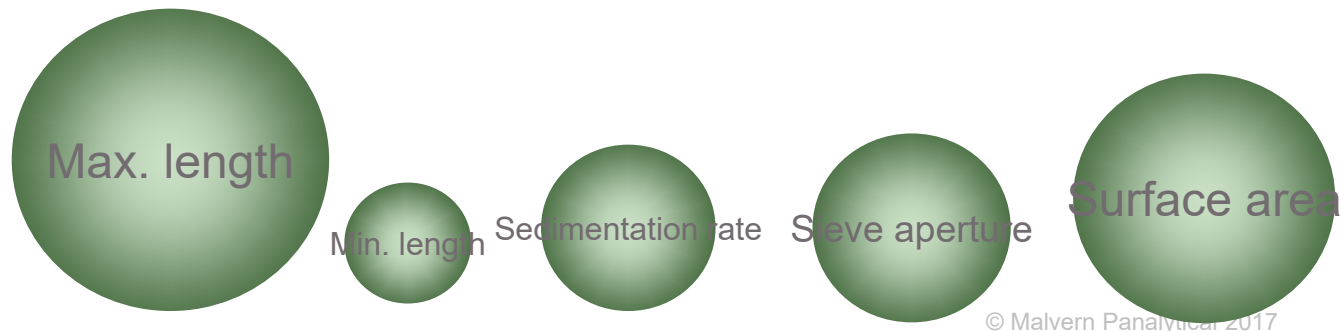
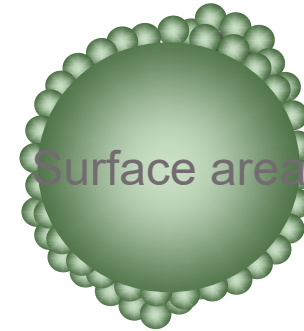
How do we describe the size of these particles

- Equivalent spheres
 - Maximum length
 - Minimum length
 - Sedimentation rate
 - Sieve aperture



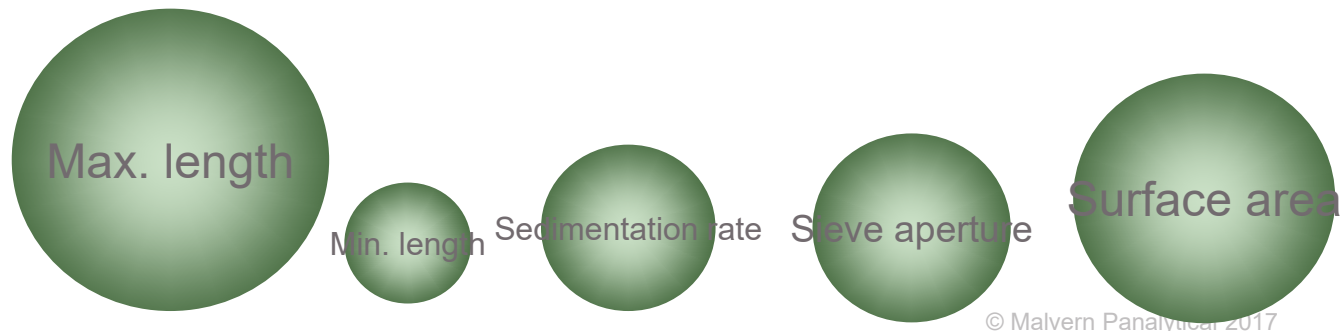
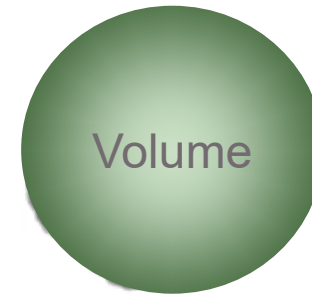
How do we describe the size of these particles

- Equivalent spheres
 - Maximum length
 - Minimum length
 - Sedimentation rate
 - Sieve aperture
 - Surface area



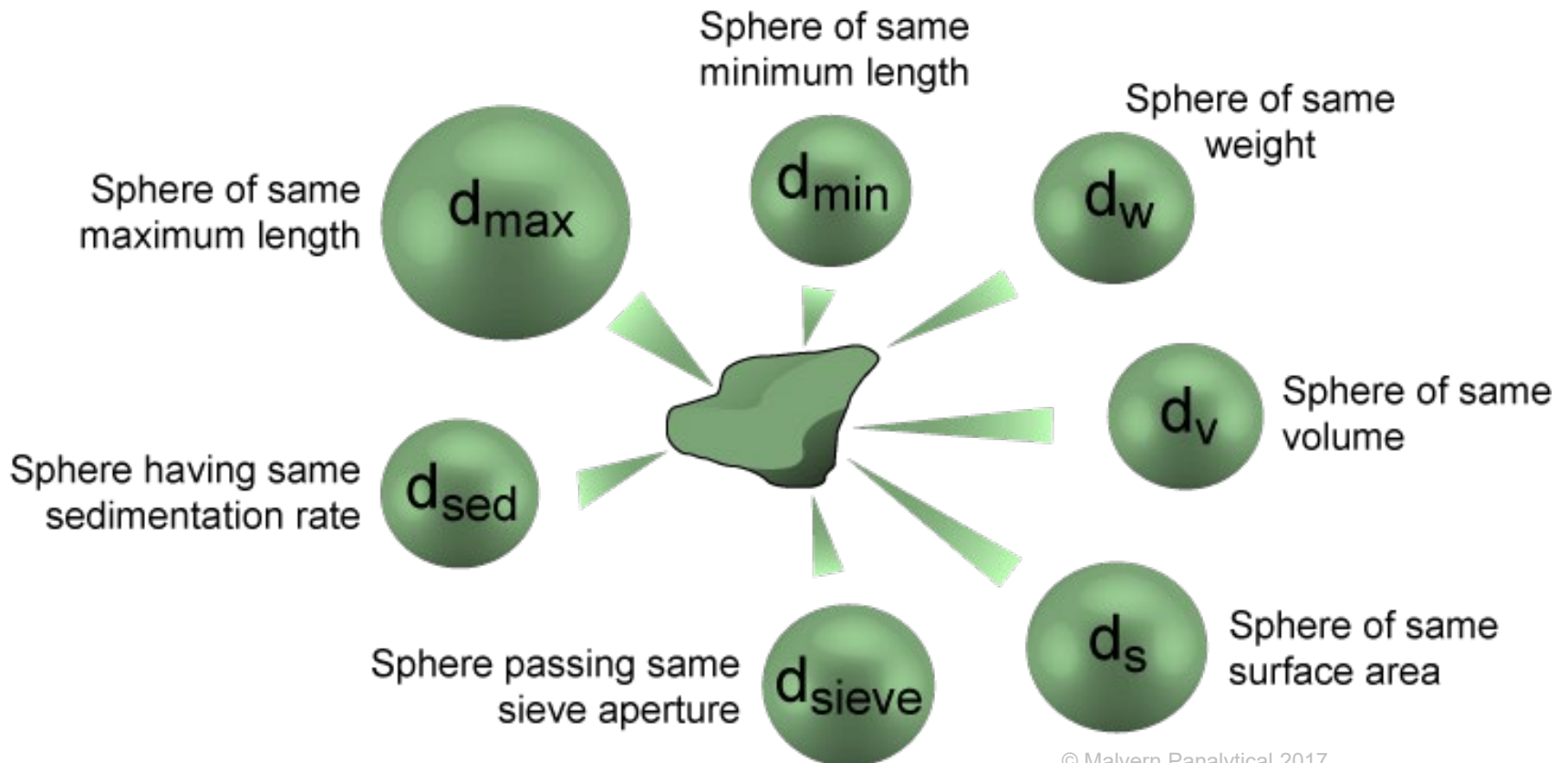
How do we describe the size of these particles


- Equivalent spheres
 - Maximum length
 - Minimum length
 - Sedimentation rate
 - Sieve aperture
 - Surface area
 - Volume



Concept of equivalent spherical diameters

- Different particle sizing techniques report different equivalent spherical diameters
 - Dependent on the physical property that is measured

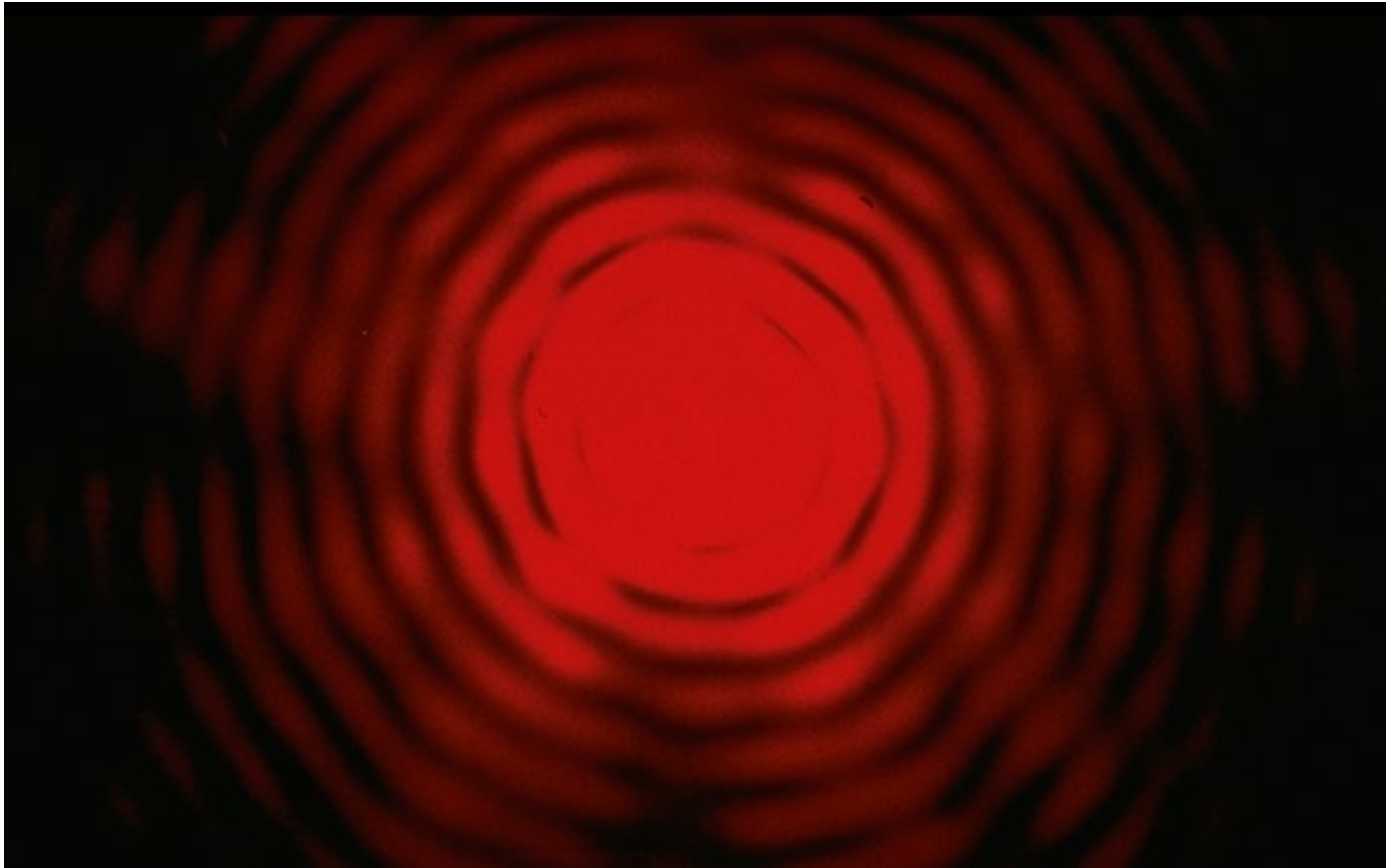


The background of the slide features a series of concentric, slightly blurred red circles that create a ripple effect, centered on the left side of the frame. The circles are more prominent in the center and fade towards the edges.

How does laser
diffraction work?



Laser Diffraction – The light scattering pattern from a group of particles





Malvern
Panalytical

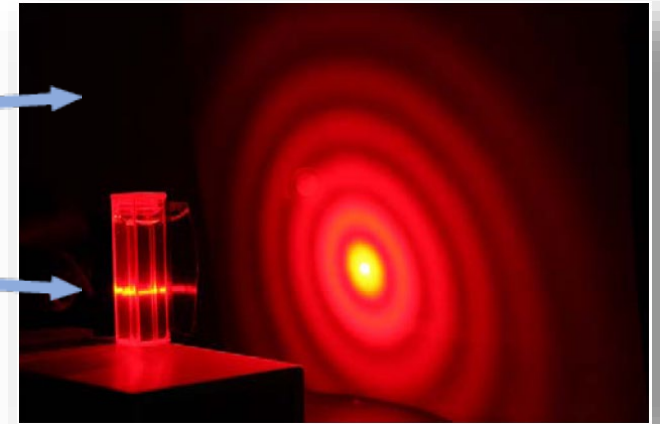
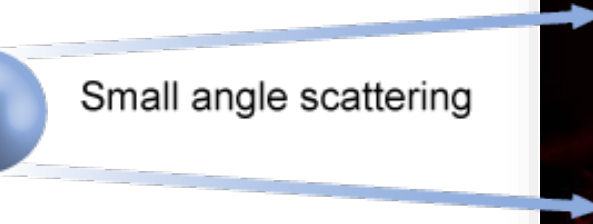
Large particles

Laser diffraction: light scattering

Incident light



Small angle scattering

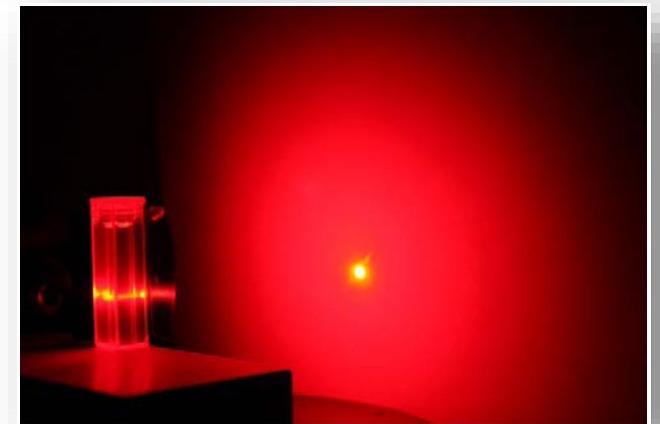
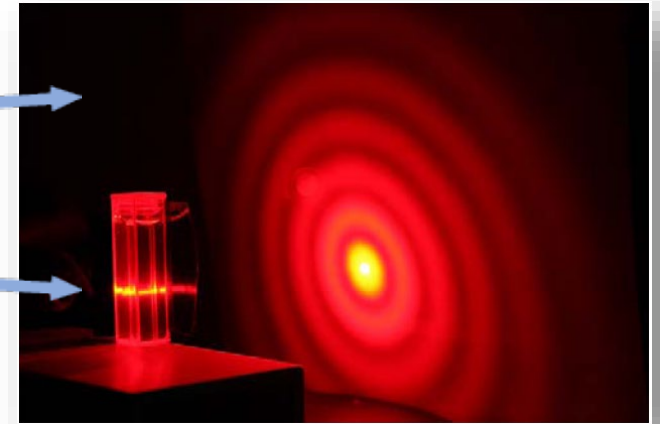




Malvern
Panalytical

Laser diffraction: light scattering

Large particles



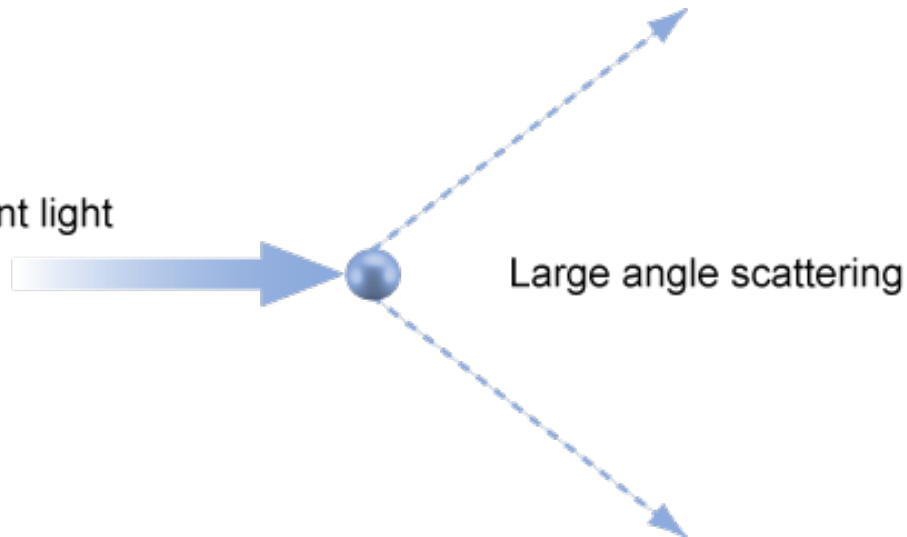
Submicron
particles

Incident light



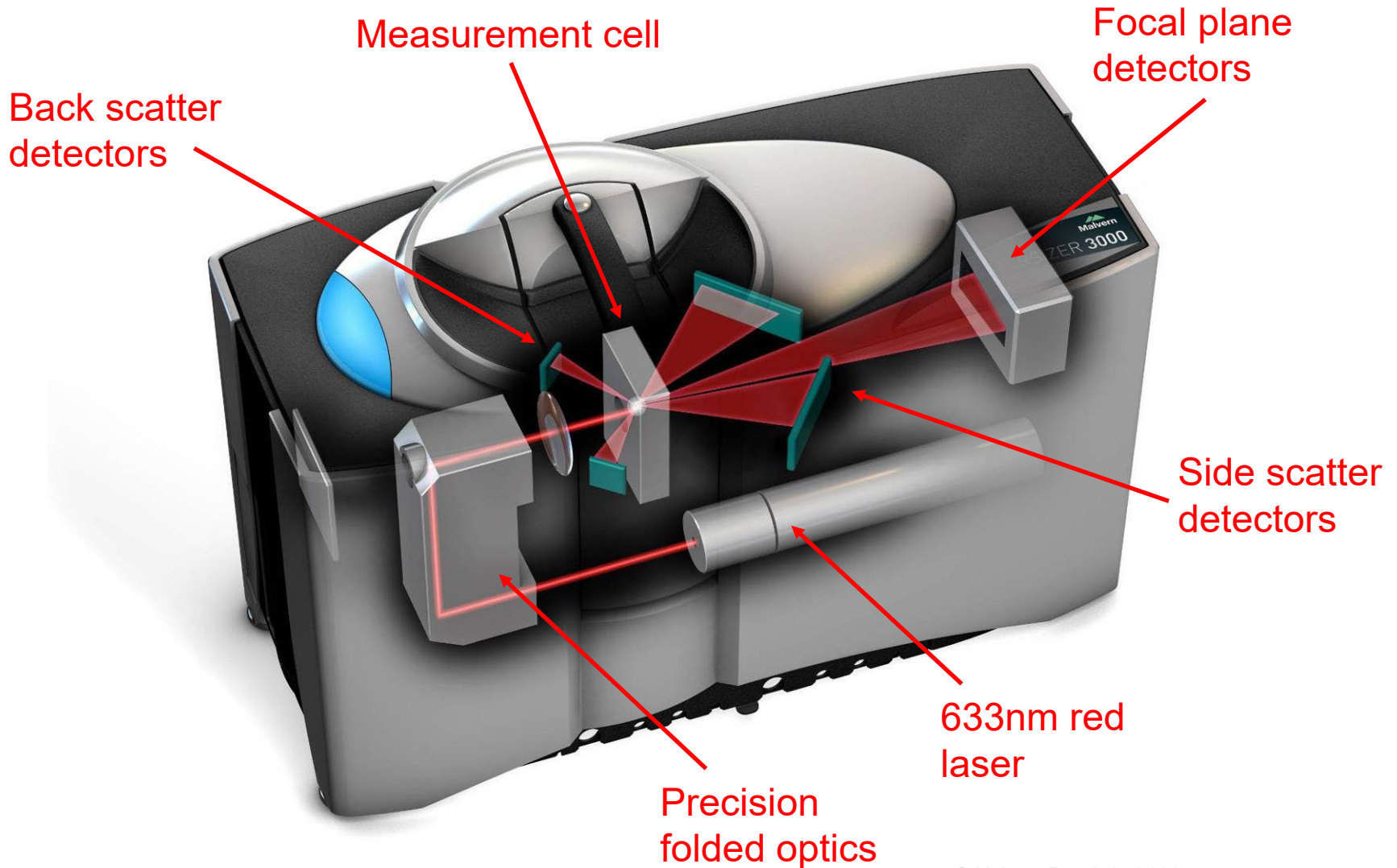
Small angle scattering

Incident light

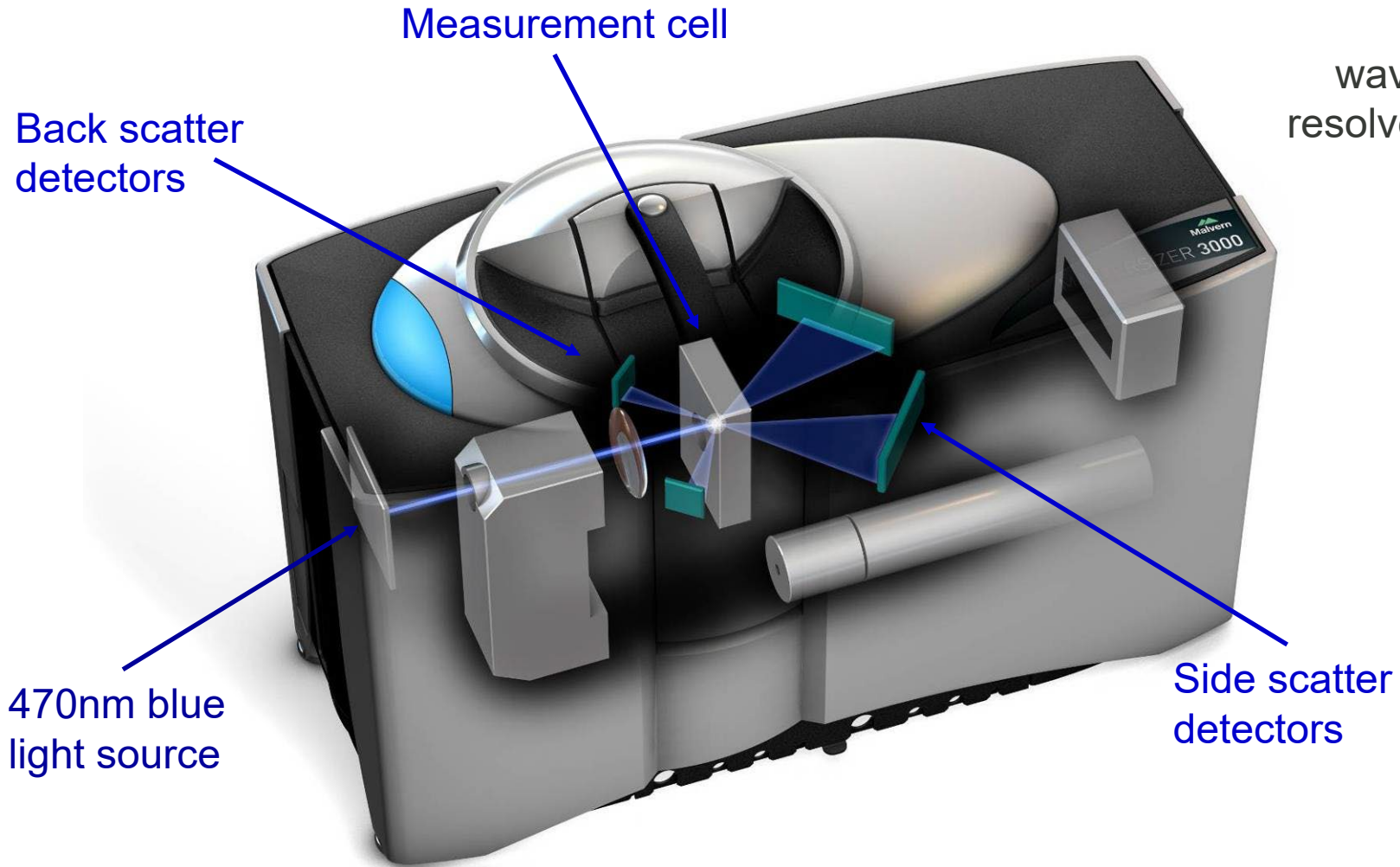


Large angle scattering

Mastersizer 3000 optics: red light measurements

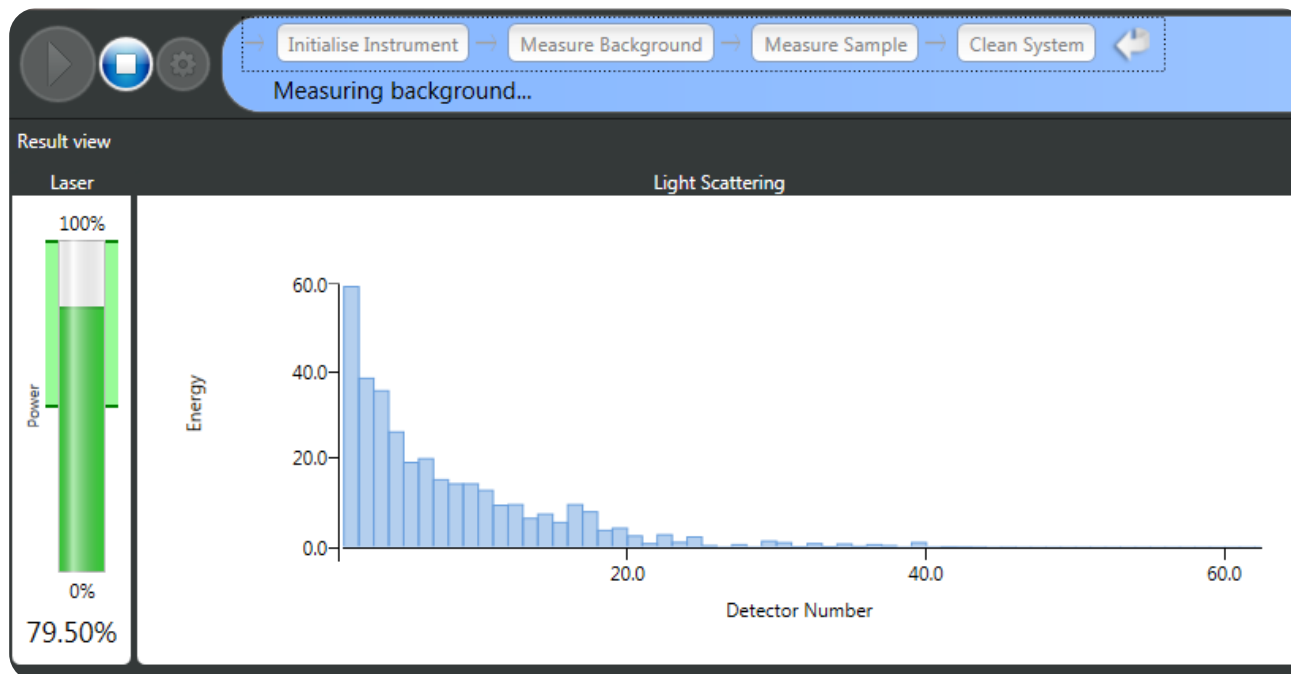


Mastersizer 3000 optics: blue light measurements

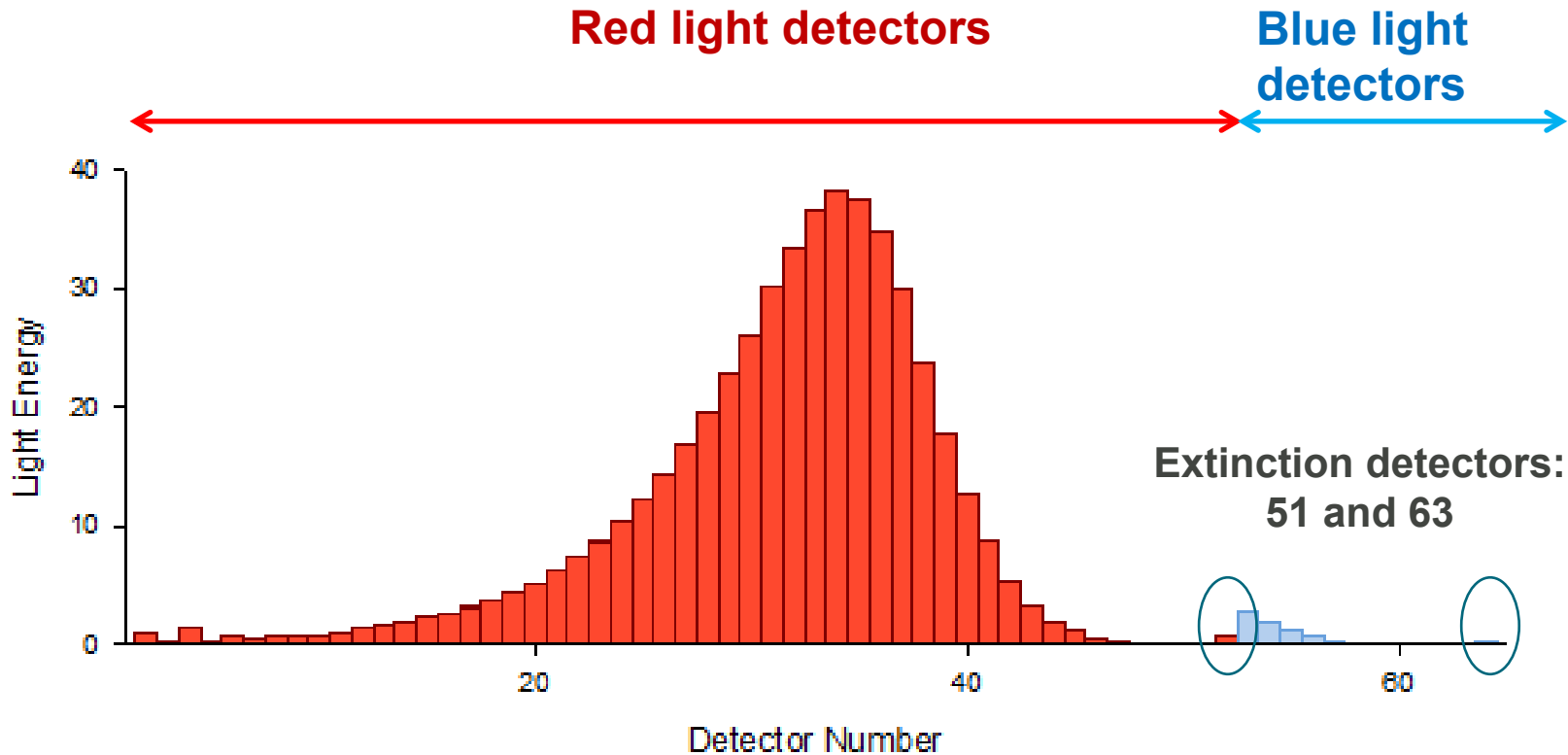


Measuring the scattering data

- Angular scattering data is presented in real-time in the measurement manager
- Increasing detector numbers represent increasing angle

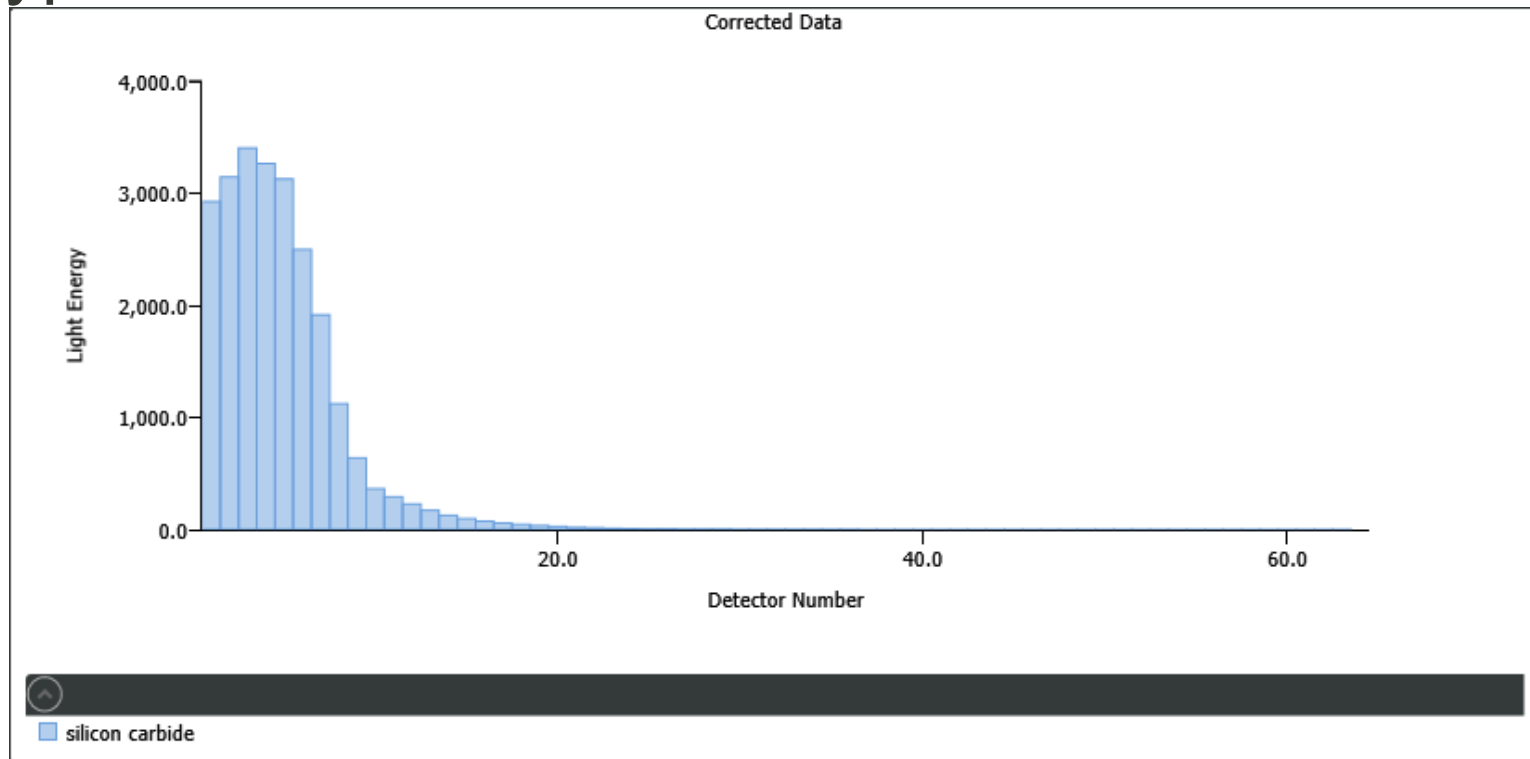


The measured scattering data



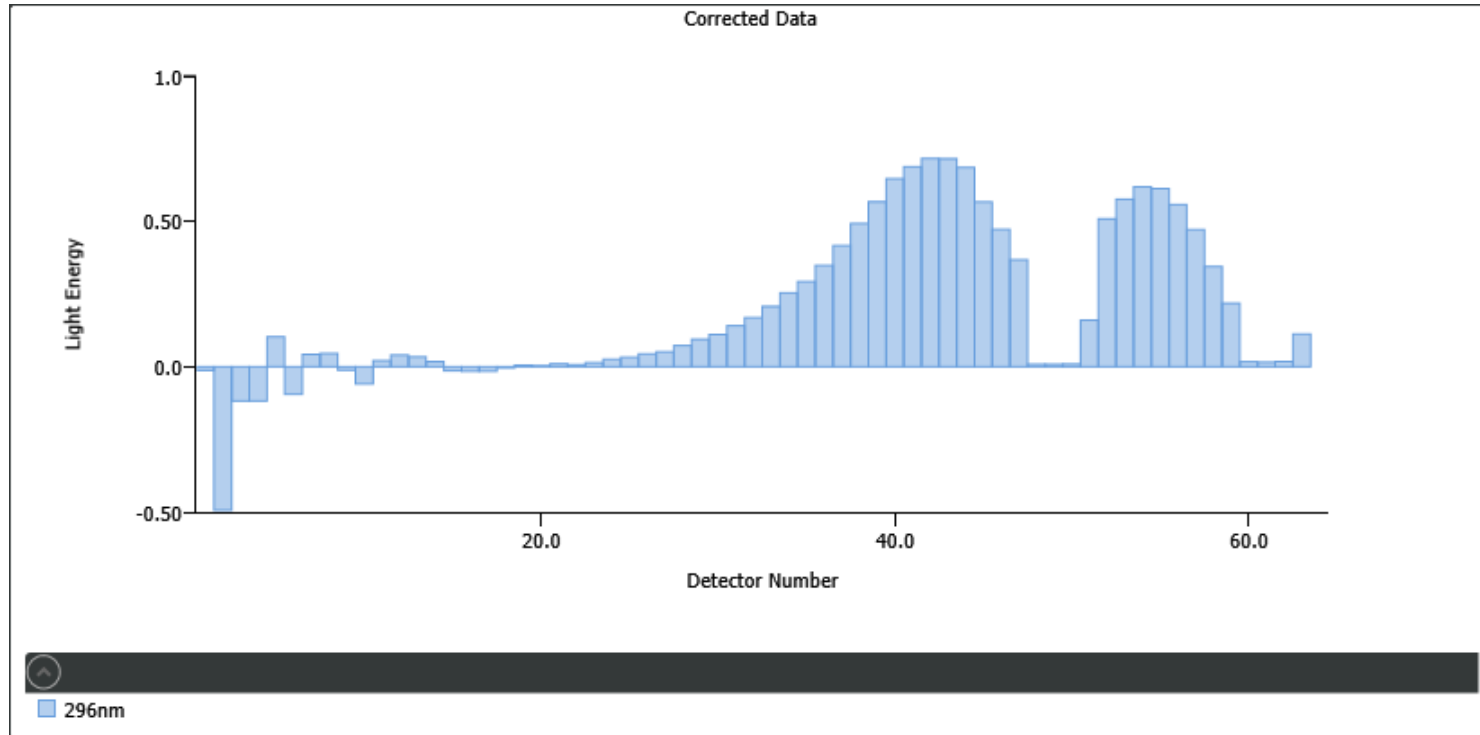
Increasing angle / Decreasing particle size

The Mastersizer 3000 Typical Data Set – Coarse Particles

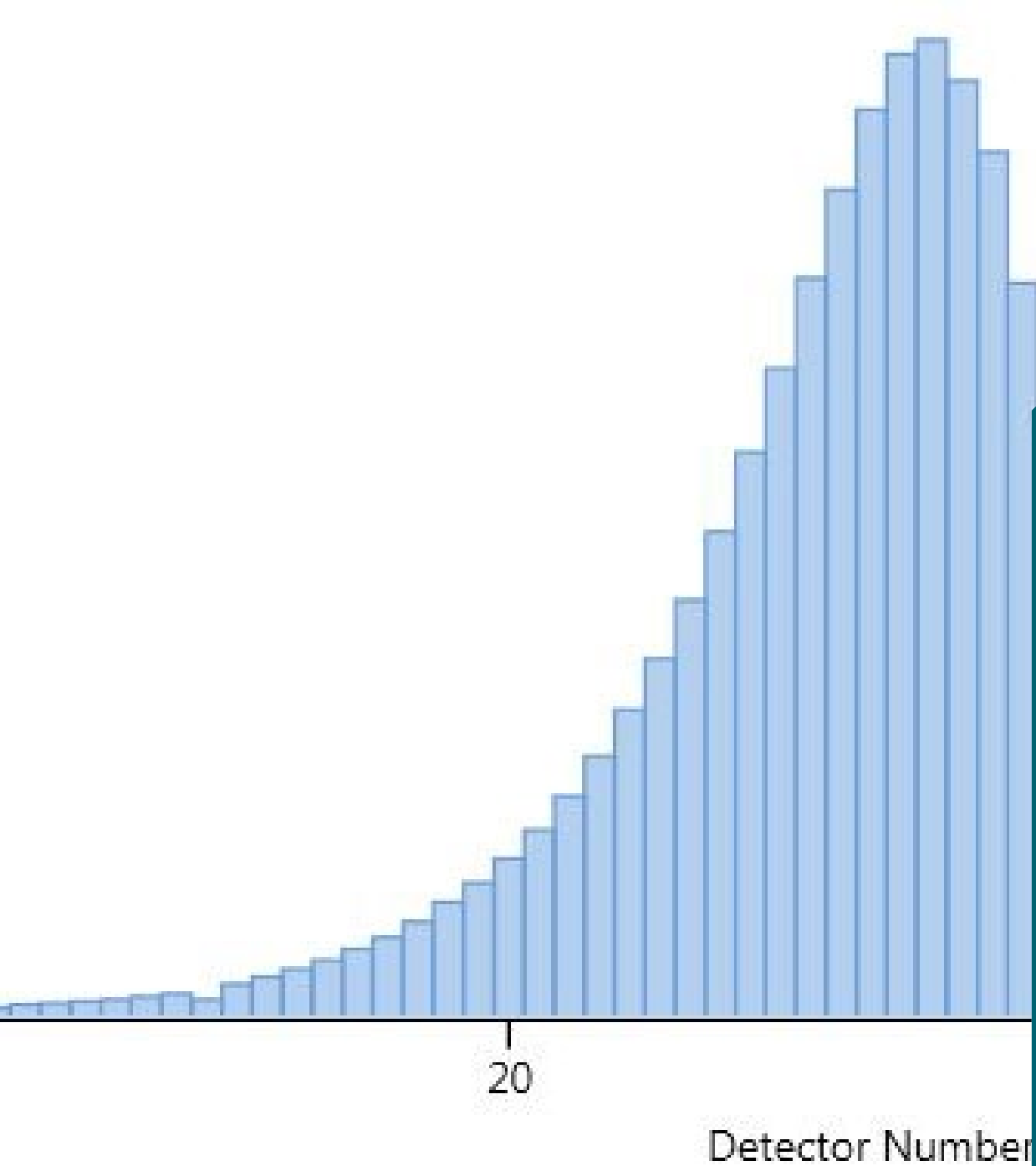


Large particle scattering is concentrated in the low angle region which corresponds to low detector numbers.

The Mastersizer 3000 Typical Data Set – Sub-Micron Particles



Small particles scatter light at high angles which produces data in the high detector number region.

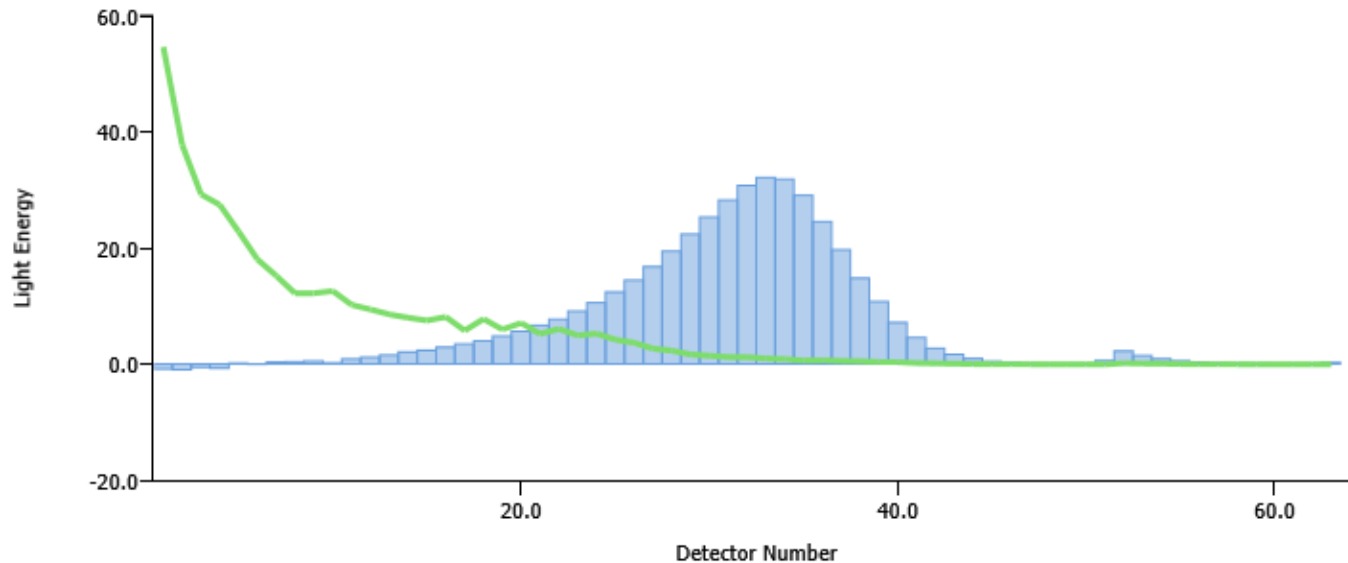


Data quality



Data quality - introduction

- Data is the fundamental light scattering caused by the sample

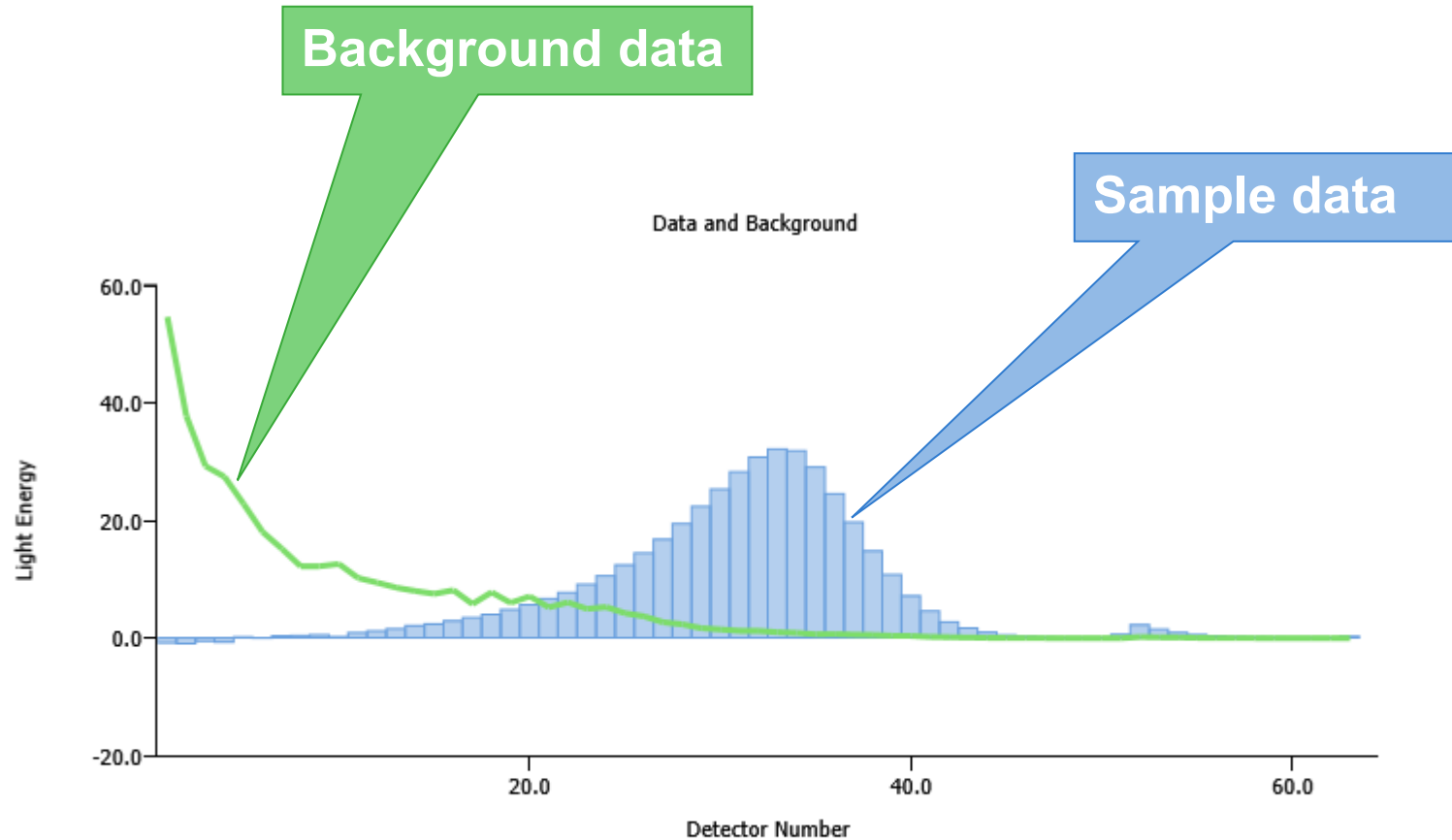


- Data is not the particle size result (.pdf)
- Data is independent of the optical model
- A stable result requires stable data

What is good quality data?

- A good background measurement shows:
 - Clean cell windows and dispersant
 - Good alignment of the system
 - Stability of the dispersant
- Good sample measurement should have:
 - Sufficient signal to noise ratio
 - Limited negative data
 - No multiple scattering
 - No beam steering

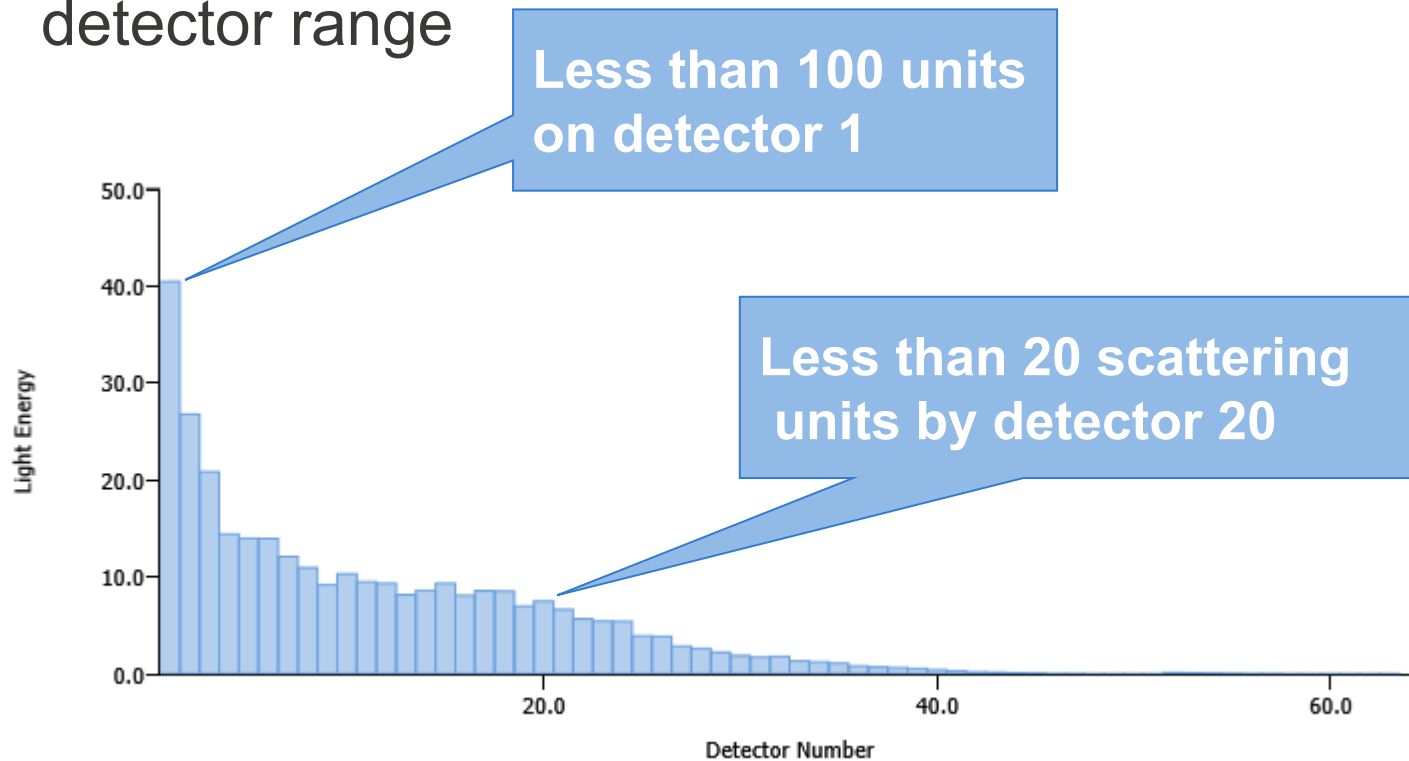
Data components - types of data



Corrected Data-sample 1 post ultras Background data-sample 1 post ultra

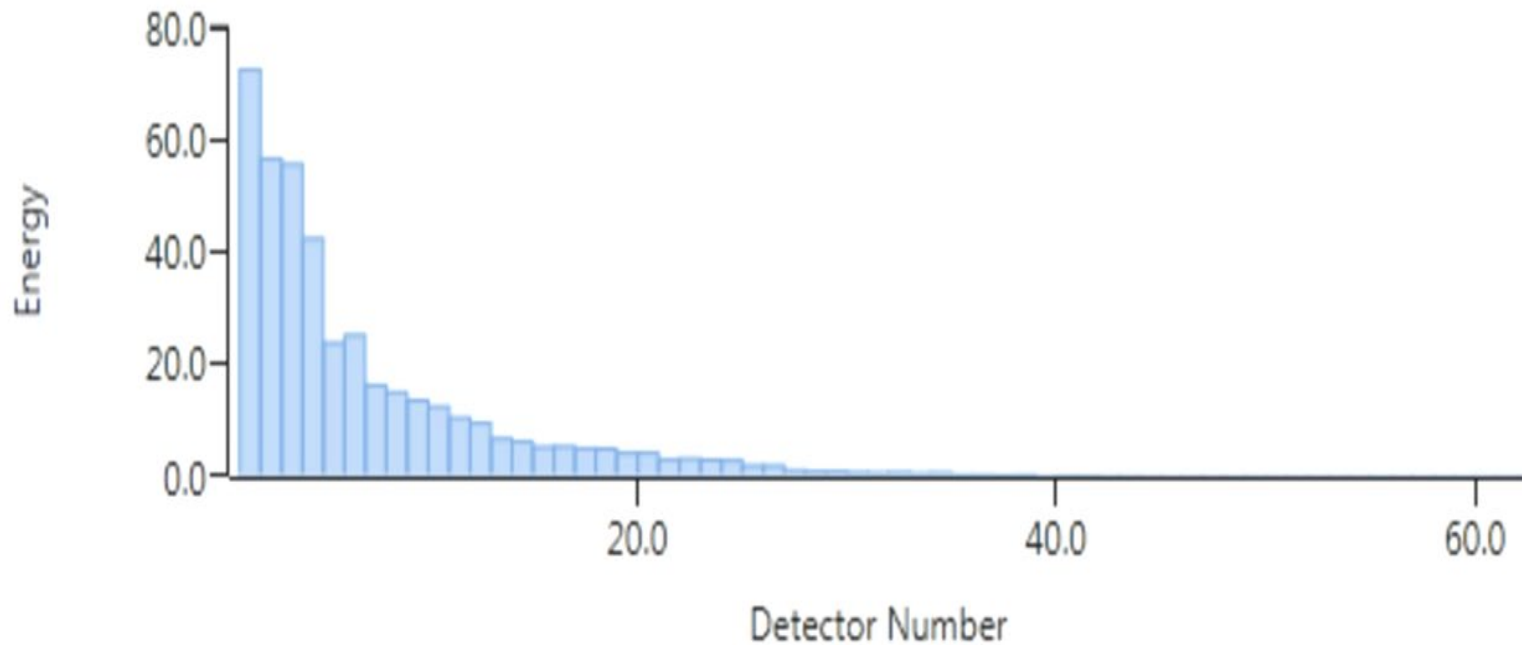
Background data and system cleanliness

- A good measurement requires a clean, stable background
- This should show progressive decrease across the detector range



A clean background – wet system

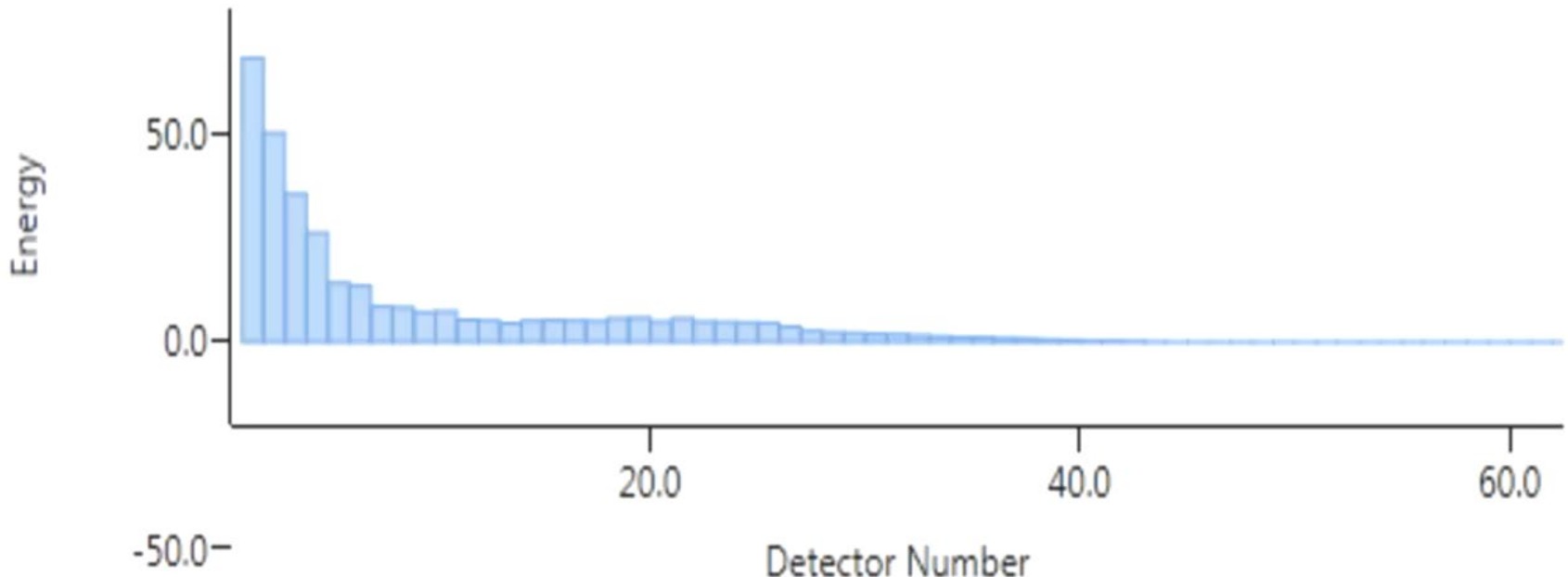
- A good clean background on a wet system should look very similar to this...



- Less than 100 on detector 1
- Less than 20 on detector 20
- Decreasing curve
- Limited fluctuations

A clean background – dry system

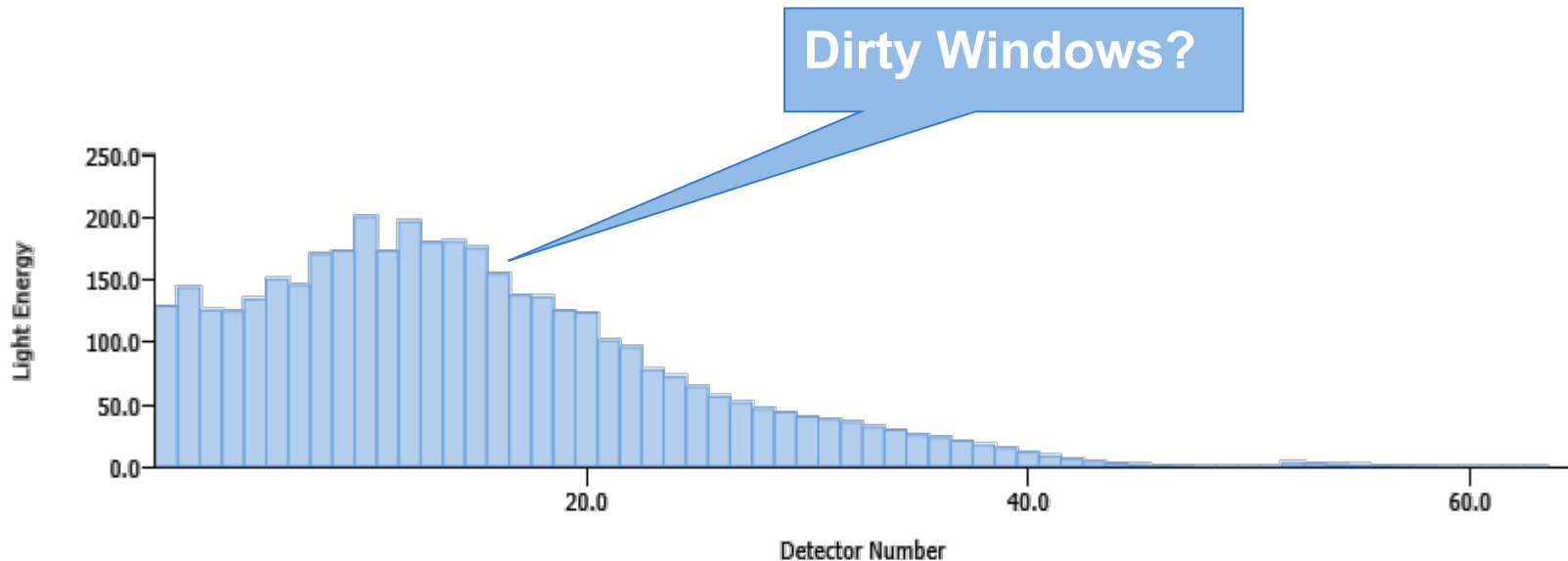
- The air flow causes larger fluctuations in the background than in a wet measurement.
- The same rules apply!



- Less than 100 on detector 1
- Less than 20 on detector 20
- Decreasing curve
- Limited fluctuations

Poor background - material stuck to the windows

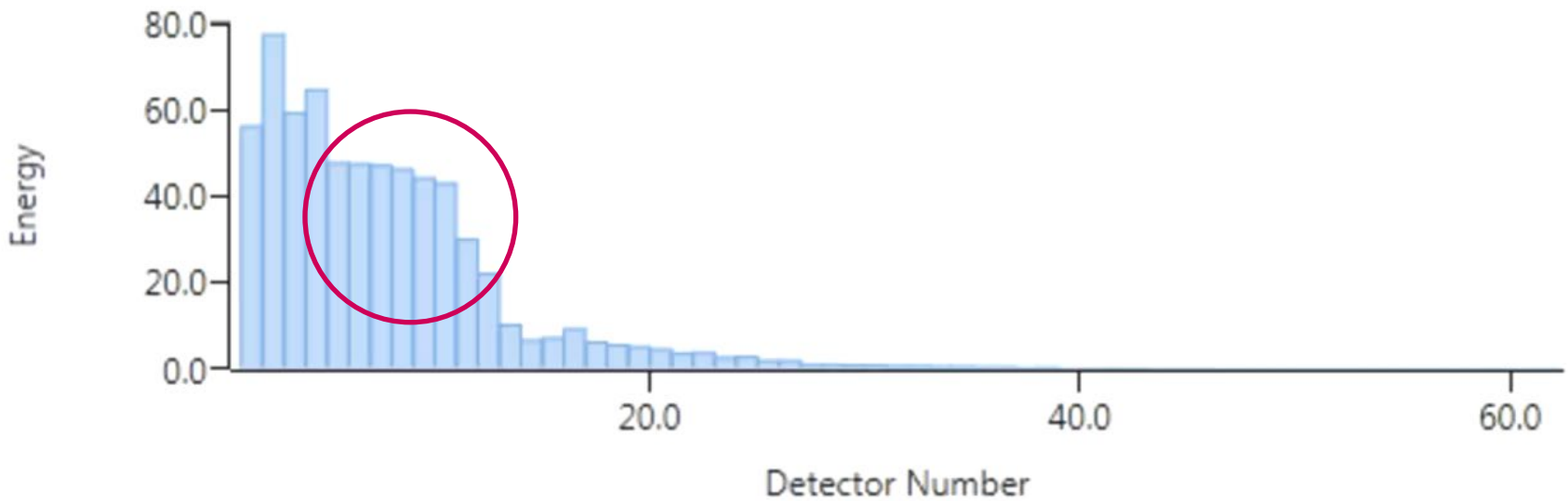
- A 'hump' in the data is often an indication of material stuck to the cell windows



- All scattered light is included in the measurement

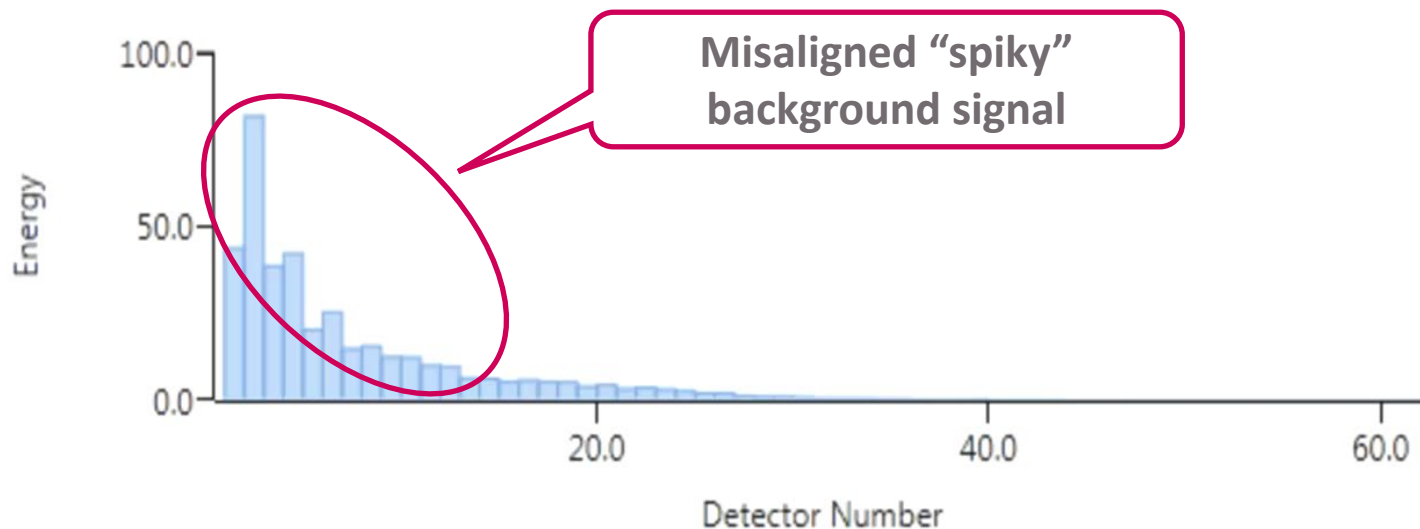
Poor background – contaminated dispersant

- Intermittent peaks in the background may be caused by contaminants in the dispersant
 - Particulates (rinse the dispersion unit)
 - Bubbles (degas dispersant, stop-start pump)



Poor background – misaligned system.

- A spiky background signal indicates misalignment
 - Detectors are arranged on opposite sides of the pinhole
- Misalignment can be caused by
 - Contamination on the cell windows
 - A change in the refractive index of the dispersant



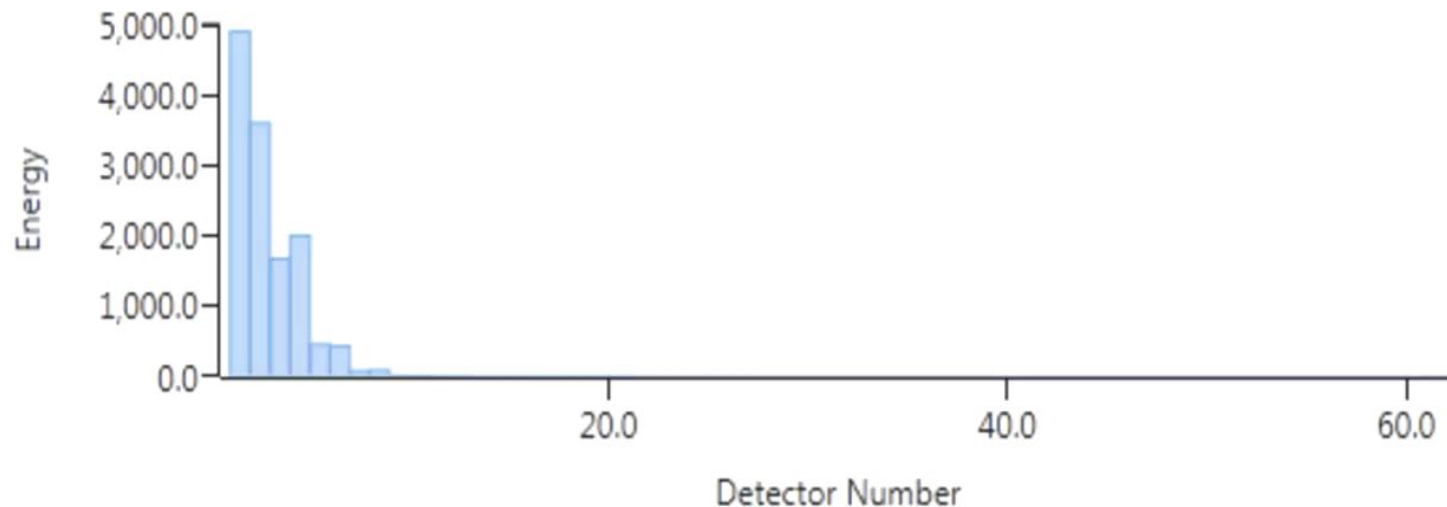
Poor background: instability due to thermal gradients

- The dispersion unit may be warmer than the dispersant
- This temperature difference causes thermal gradients in the dispersant
 - And high backgrounds and possible alignment problems



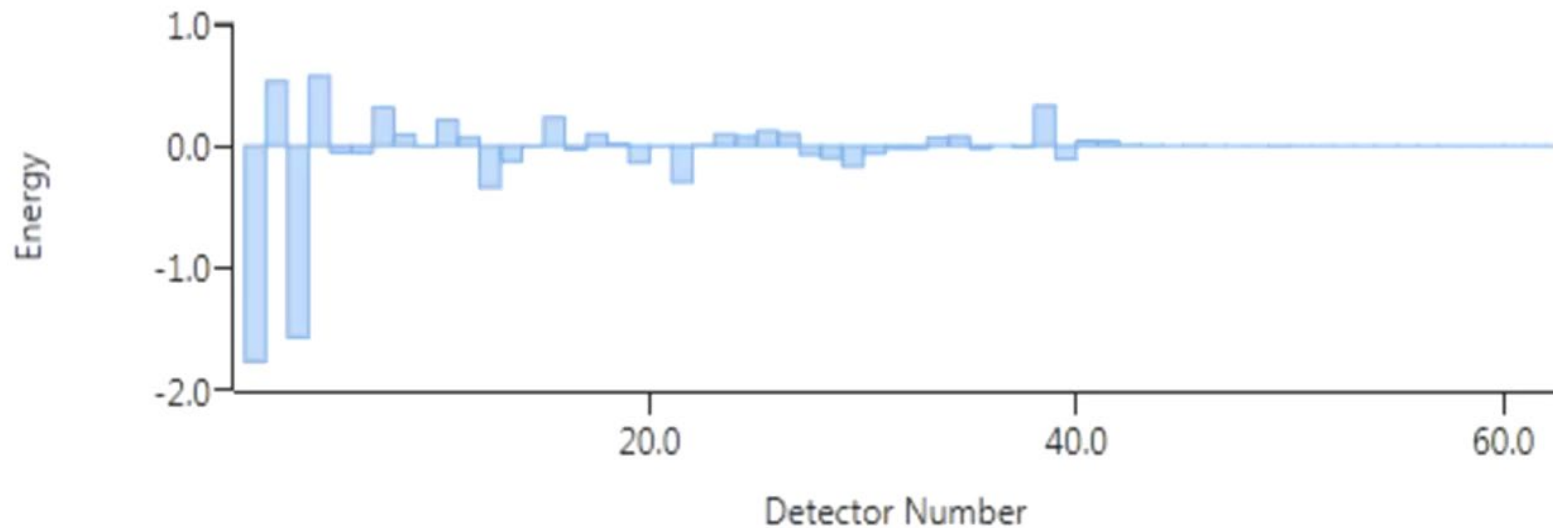
How will I recognise thermal gradients?

- Thermal gradients cause high background signals, and large fluctuations
 - The background signal decreases as the temperature stabilises and refractive index gradients disappear
 - This will take longer for more volatile dispersants



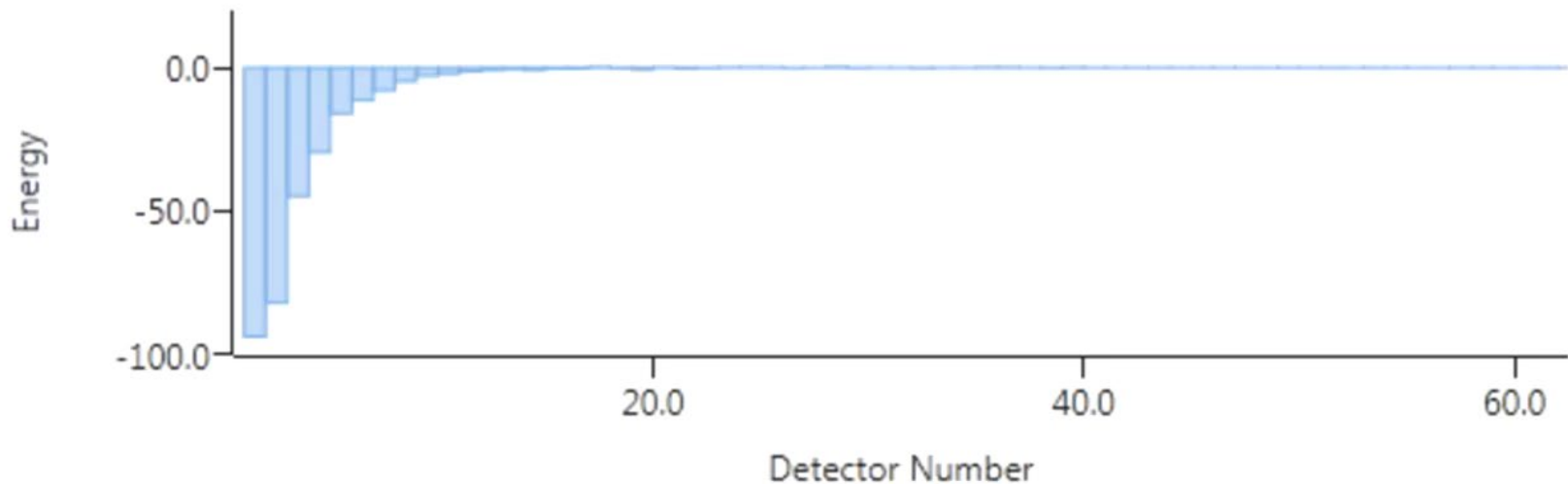
The Add Sample stage

- The background signal as been subtracted
- The live data should then show random fluctuations around zero
- Any 'blocks' of scattering will indicate dispersant contamination



The Add Sample stage – negative data

- The background signal has been subtracted
- The live data should then show random fluctuations around zero
 - Significant negative data suggest that the background had not stabilised before it was measured
 - If you see this signal, re-measure the background



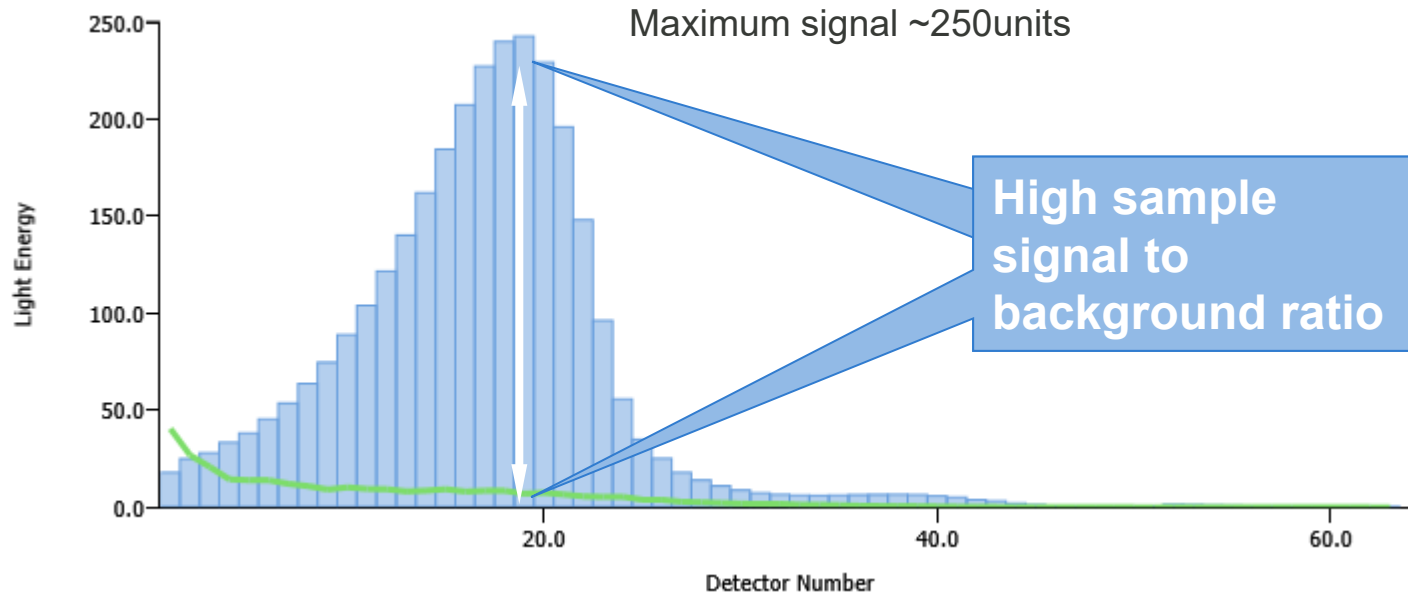
Sample addition

- How much sample should be added to the dispersion unit?
 - **Too little:**
 - Signal to noise ratio may be poor, or
 - Not enough sample may have been added to be representative of the bulk – particularly if the sample is very polydisperse
 - **Too much:**
 - Multiple scattering may affect the reported particle size distribution – particularly if the material is small (typically < 10 microns)
- What is the *correct* obscuration range?

*Obscuration = amount of laser light blocked and/or scattered by the sample,
a guide to concentration*

Low obscuration limit: signal to noise ratio

(42.58 micron glass beads measured at an obscuration of **only 7%**)

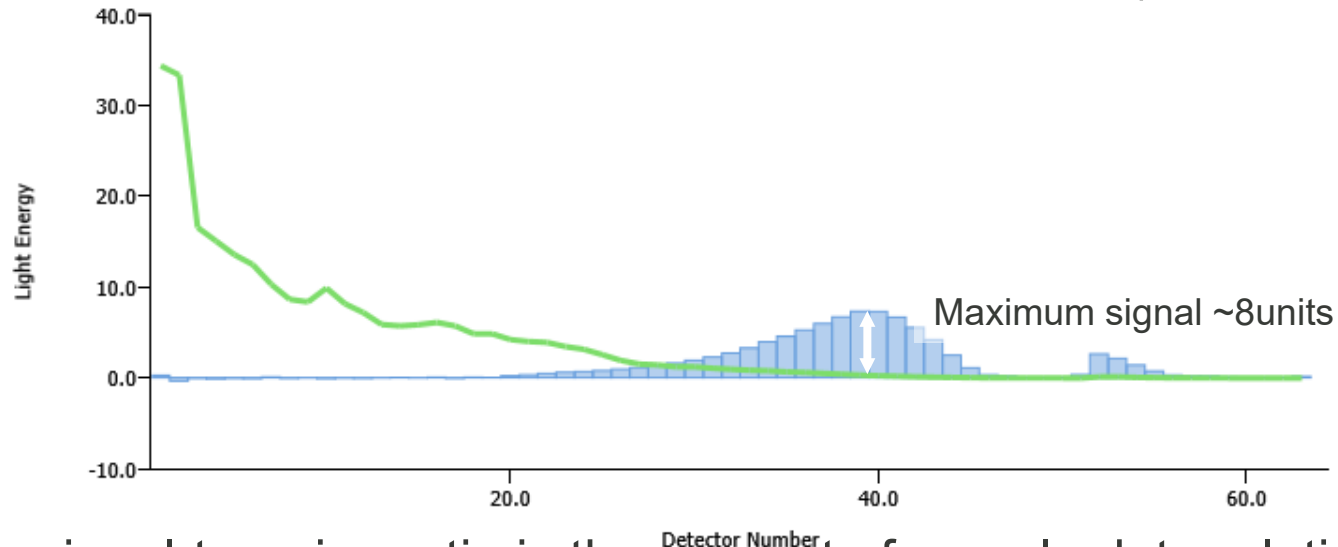


Note: the signal to noise ratio is usually high for large particles because these scatter light more strongly.

Consequently, signal-to-noise ratio is less of an issue for large particles.

Low obscuration limit: signal to noise ratio

(1 micron Latex measured at an obscuration of 5%)



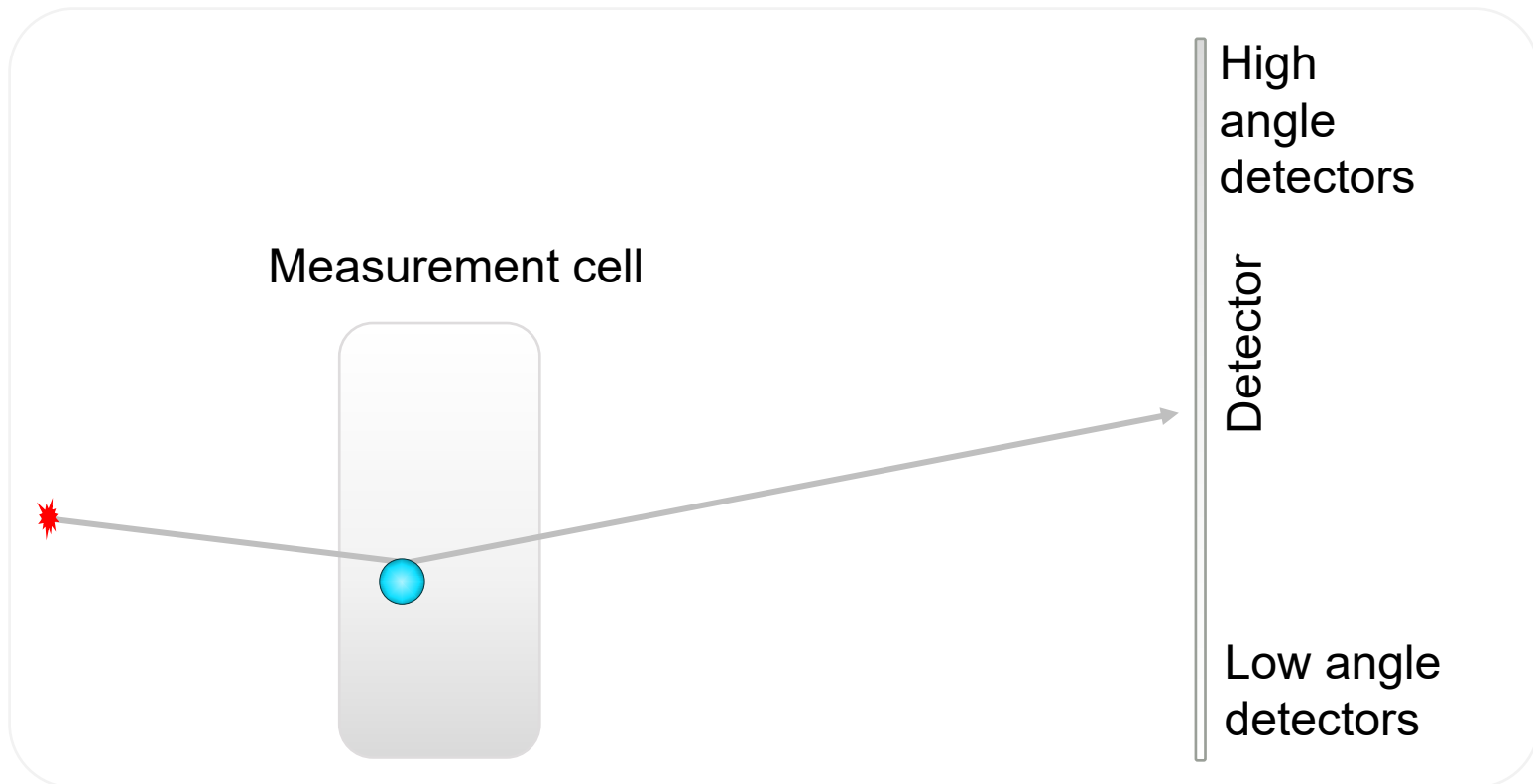
The signal-to-noise ratio is the amount of sample data relative to the background data.

Because small particles scatter light weakly, it is important that the background does not swamp the data signal.

However, in this graph, the data is good since it falls where there is little or no overlap between the sample data and the background data.

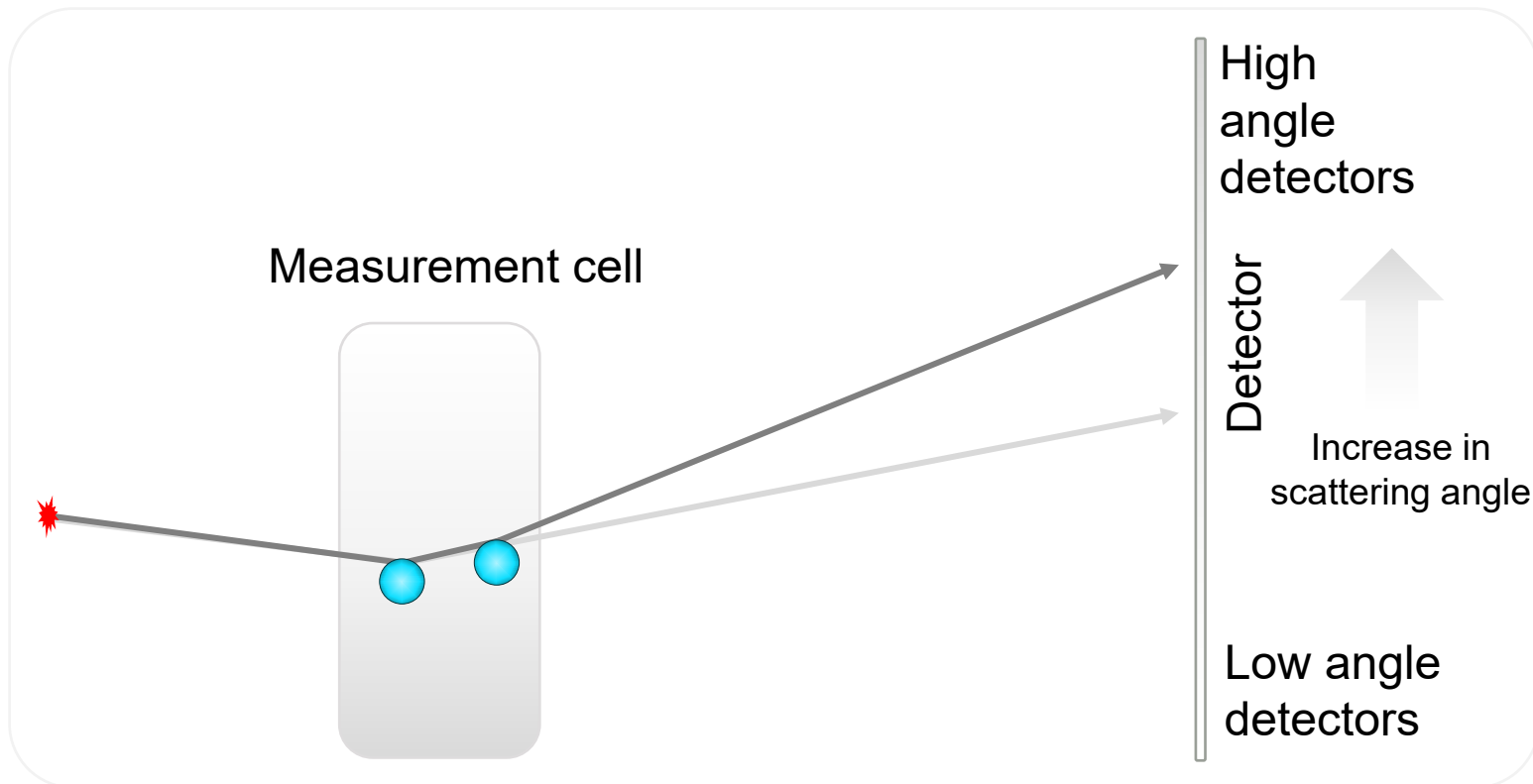
What defines the upper obscuration limit?

- If we add too much sample the results may be affected by multiple scattering
 - This generally affects samples smaller than $10\mu\text{m}$



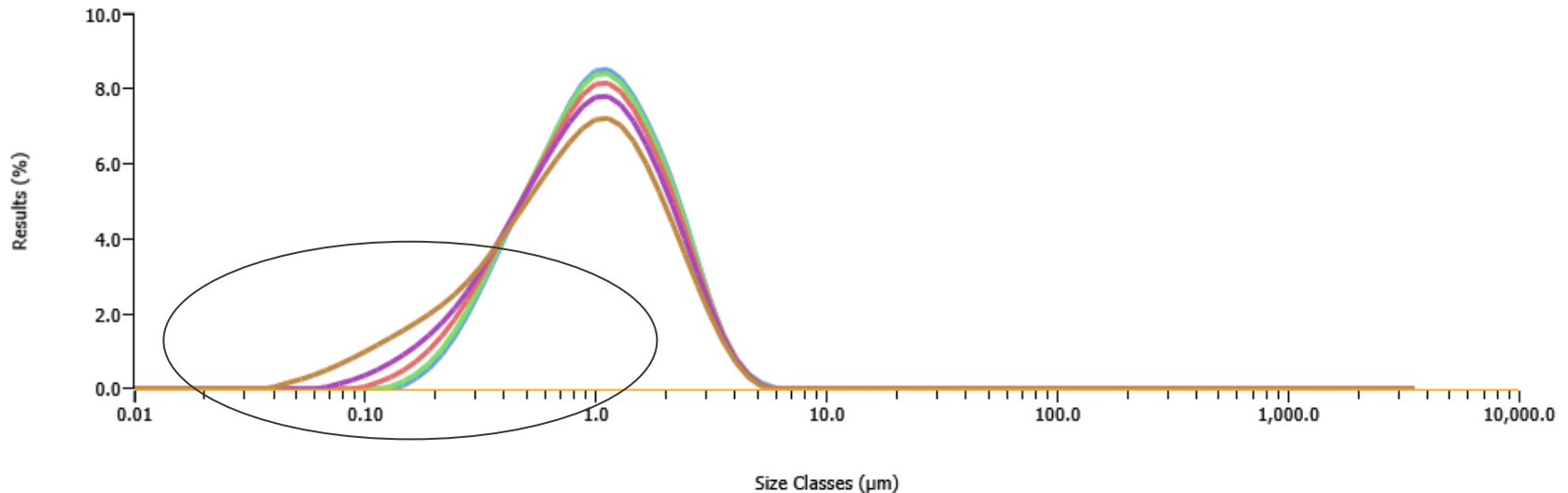
What defines the upper obscuration limit?

- If we add too much sample the results will be affected by multiple scattering
 - This generally affects samples smaller than $10\mu\text{m}$



Wet analysis - multiple scattering

....leading to exaggerated fines being interpreted



3.32 %

5.30 %

9.16 %

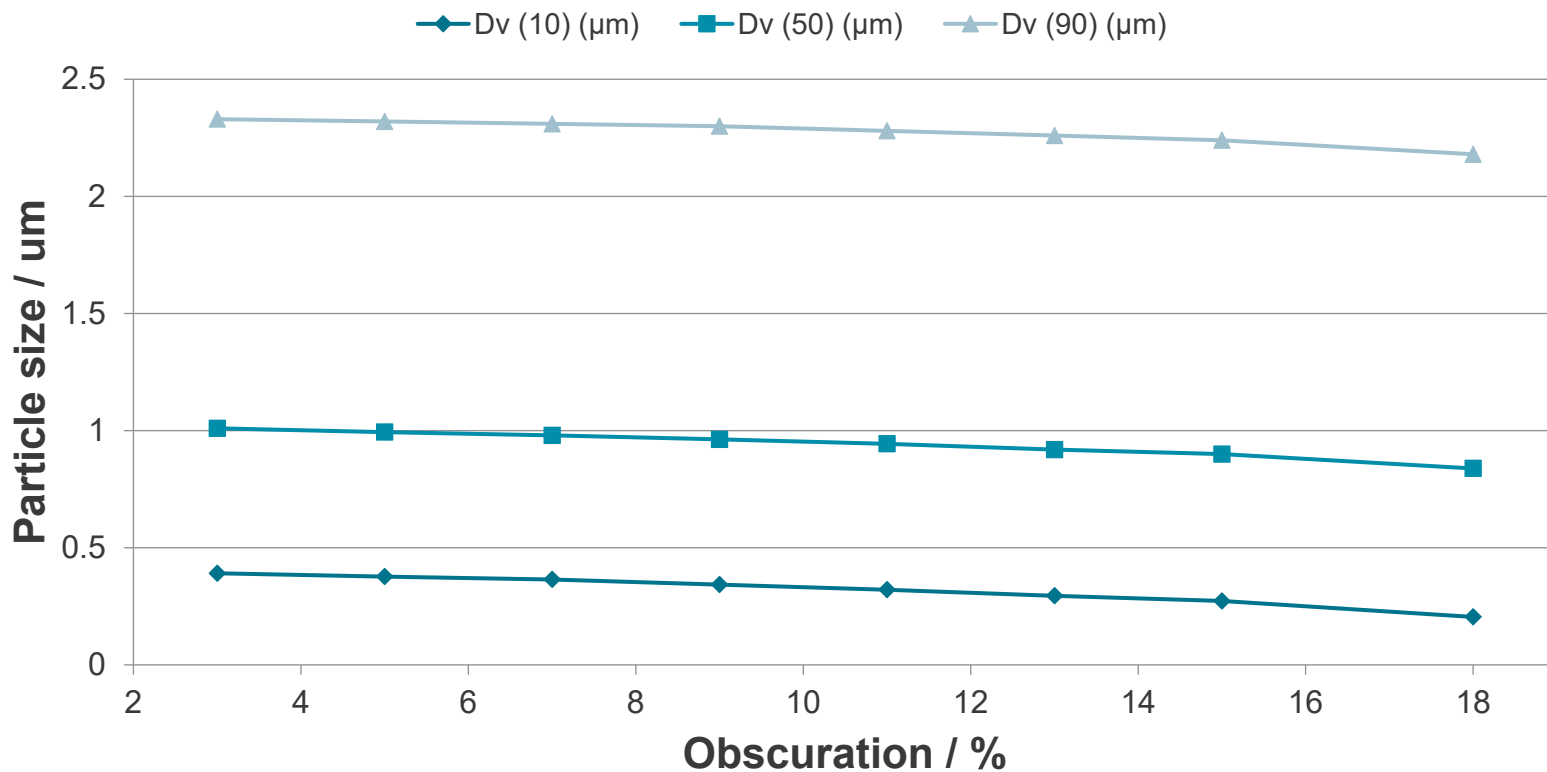
13.28 %

18.81 %

- If in doubt, carry out an obscuration titration to determine the effect of measuring at increasing obscurations on the particle size distribution.

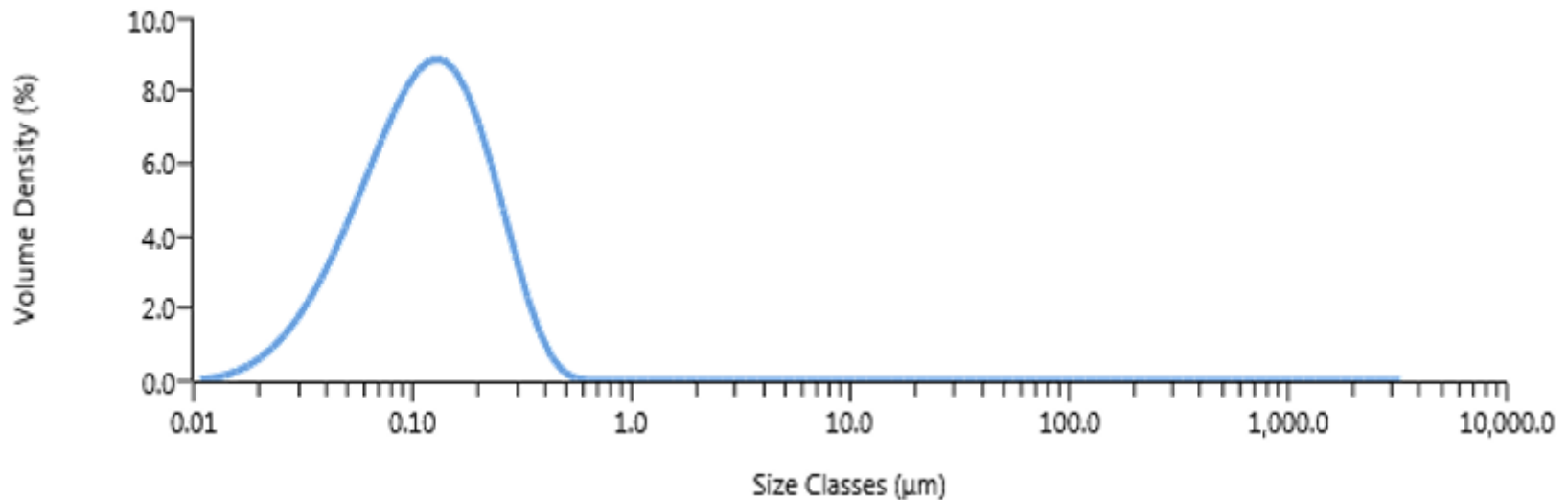
High obscuration limit: multiple scattering

- The upper limit of the obscuration range depends on multiple scattering:
 - Sample should be measured in the range where size is stable with obscuration.



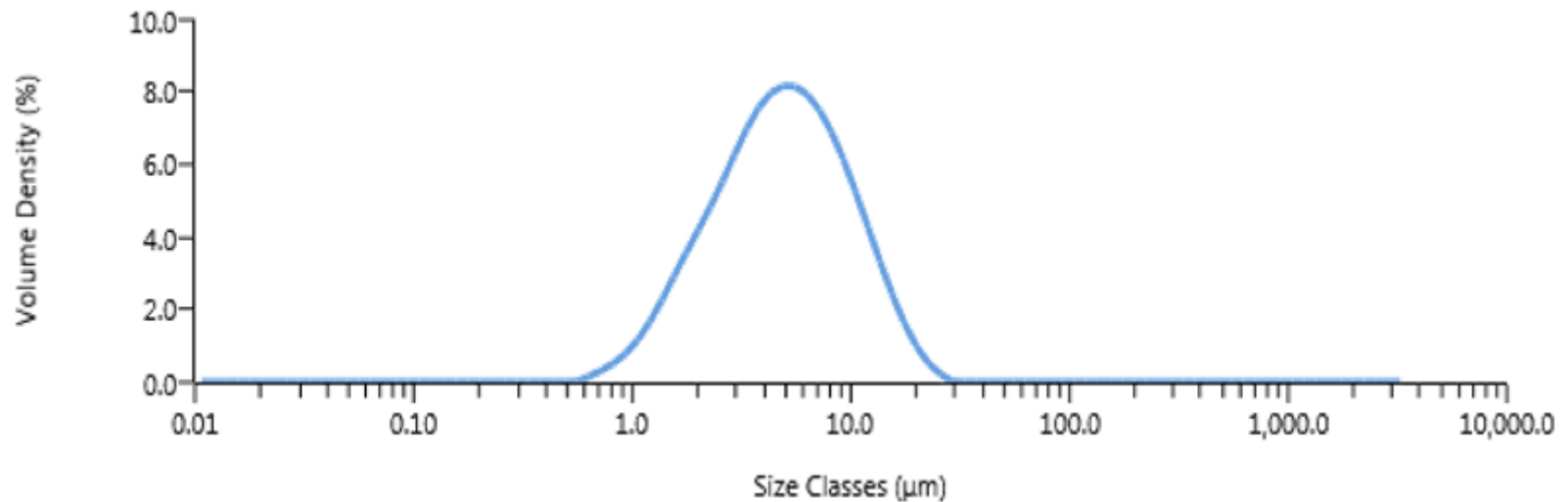
Target obscuration ranges: Wet measurements

- Very fine particles
- $<1\mu\text{m}$
- $<5\%$ obscuration



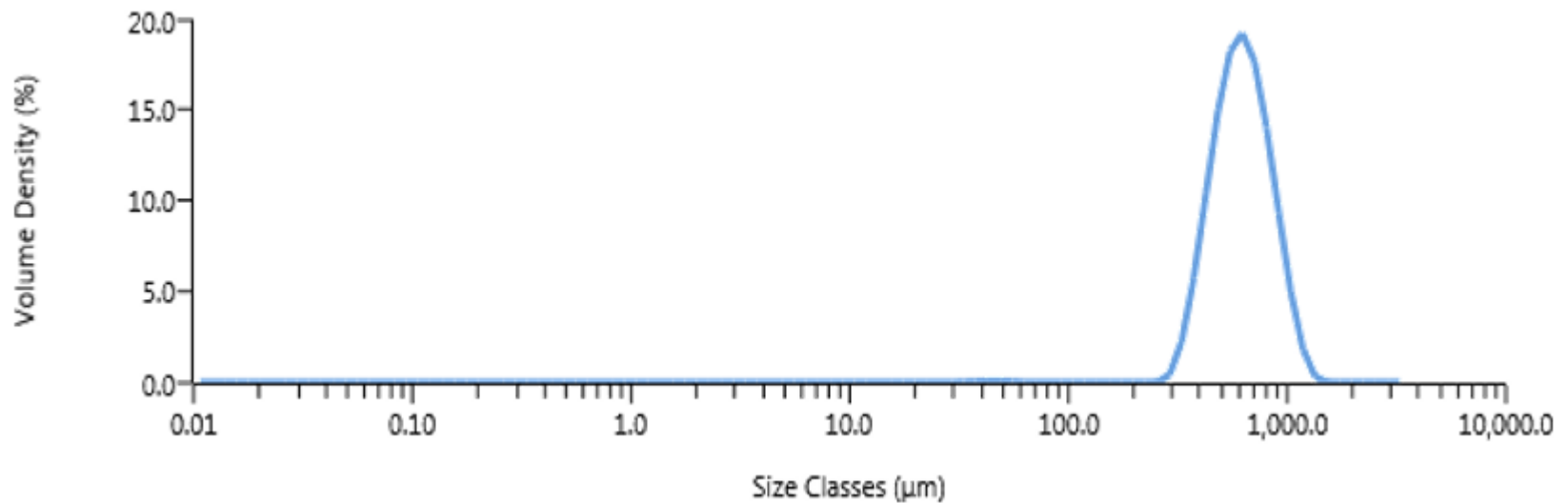
Target obscuration ranges: Wet measurements

- Fine particles
- 1-100 μm
- 5-10% obscuration



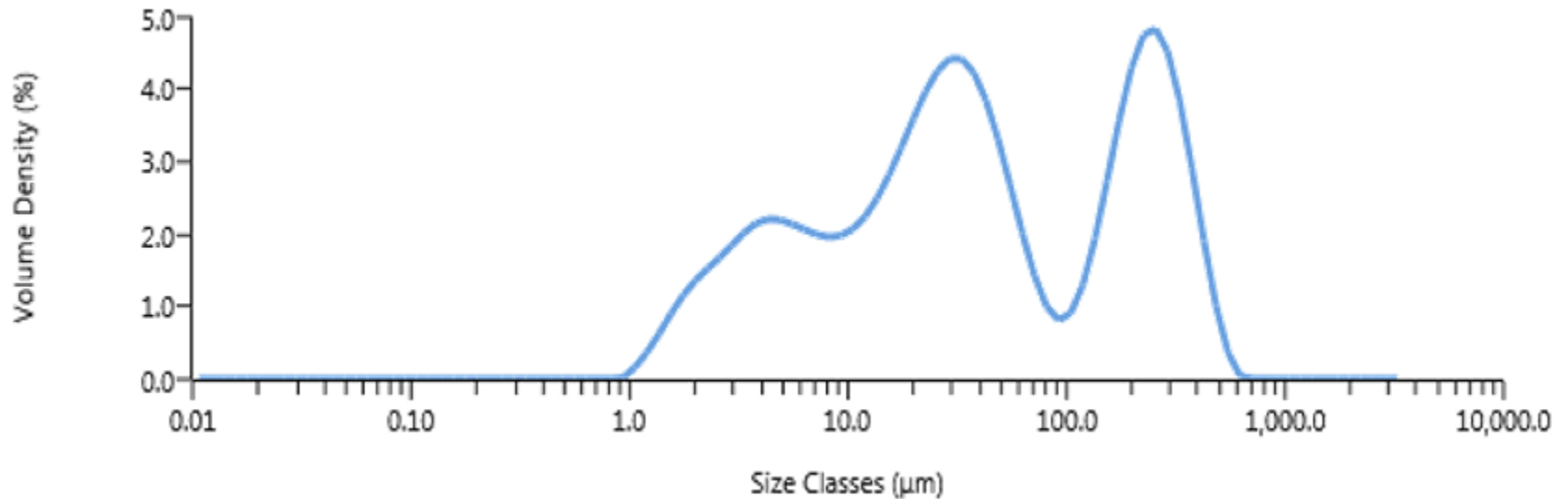
Target obscuration ranges: Wet measurements

- Coarse particles
- >100 μm
- 10-20% obscuration



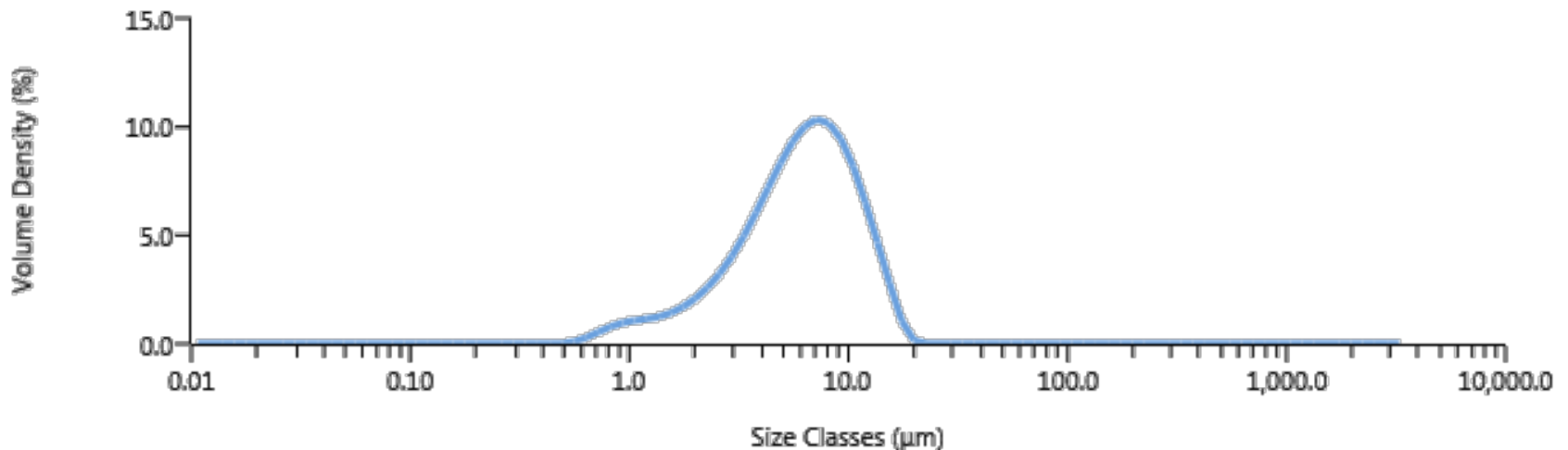
Target obscuration ranges: Wet measurements

- Polydisperse samples
- eg 1-500 μ m in one sample
- 10-20% obscuration



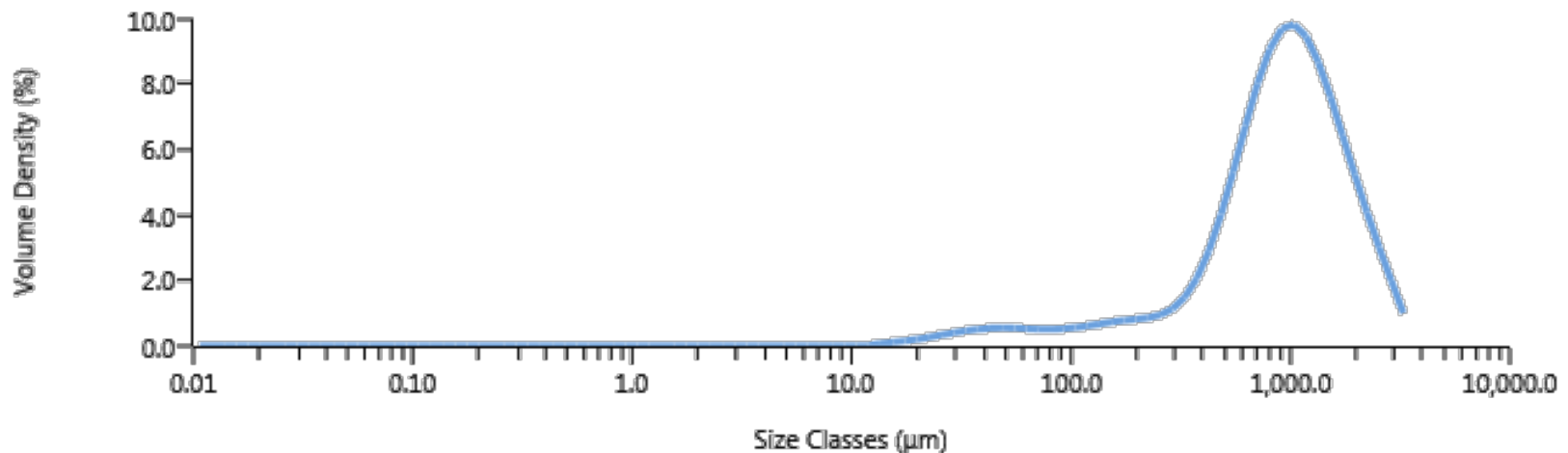
Target obscuration ranges: Dry measurements

- Fine and cohesive powders
- 0.5 to 3-5% obscuration
- Obscuration filtering ensures that only detector scans within the set obscuration range are included in the results



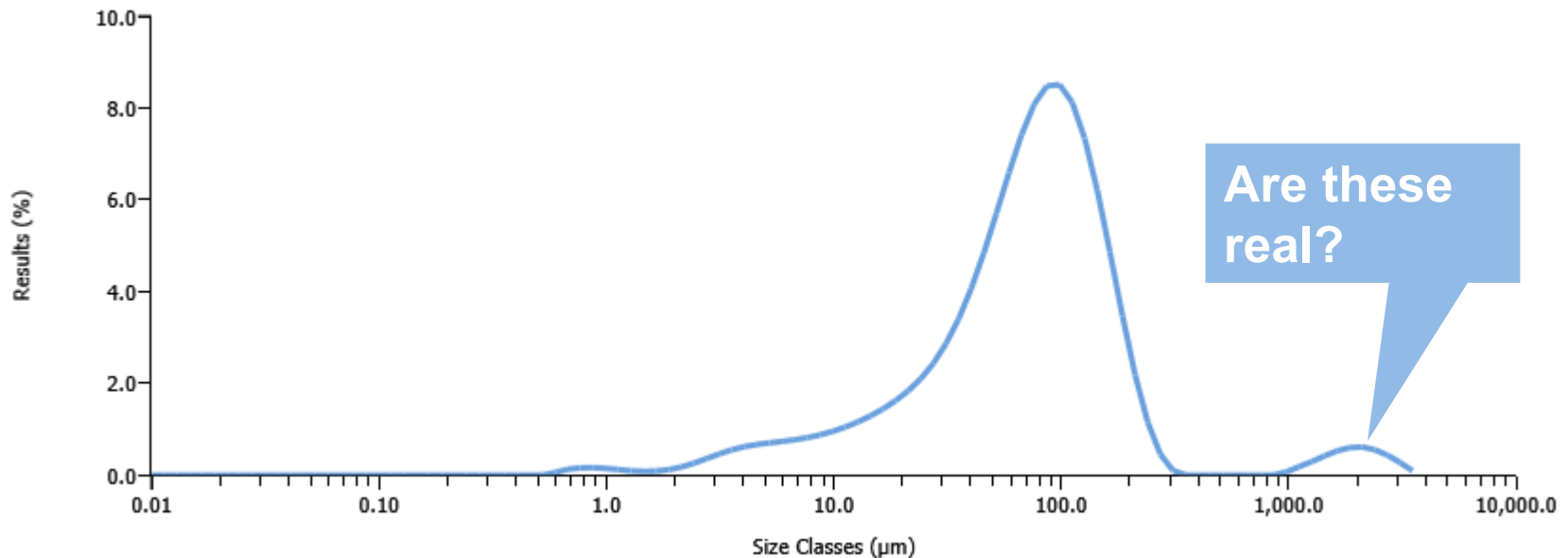
Target obscuration ranges: Dry measurements

- Coarse and free-flowing powders
- 0.1 to 6-8% obscuration
- Obscuration filtering ensures that only detector scans within the set obscuration range are included in the results



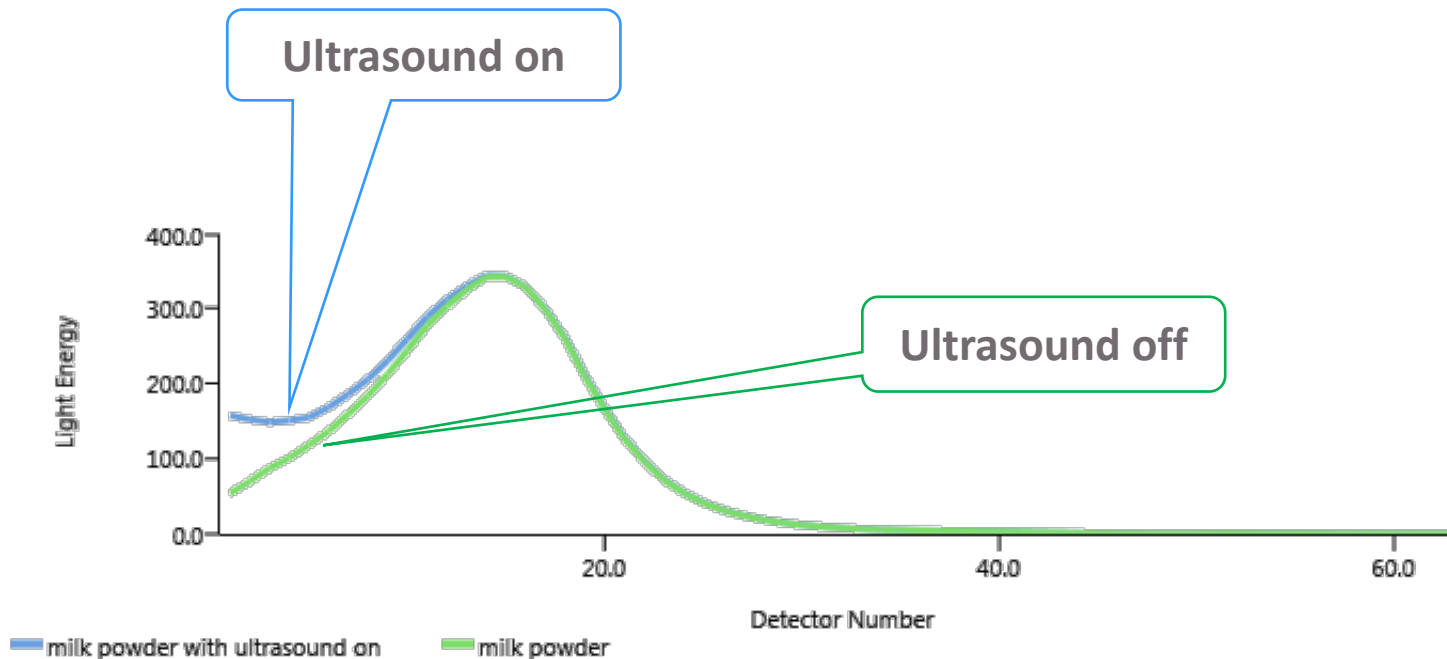
Beam Steering

- The peak on the right of this graph suggests the presence of large particles.
- However this could be caused by Beam Steering, resulting from:
 - Thermal instability following the use of ultrasound in solvents
 - Partial dissolution of the sample in the dispersant - changing its refractive index
- Or real large particles?



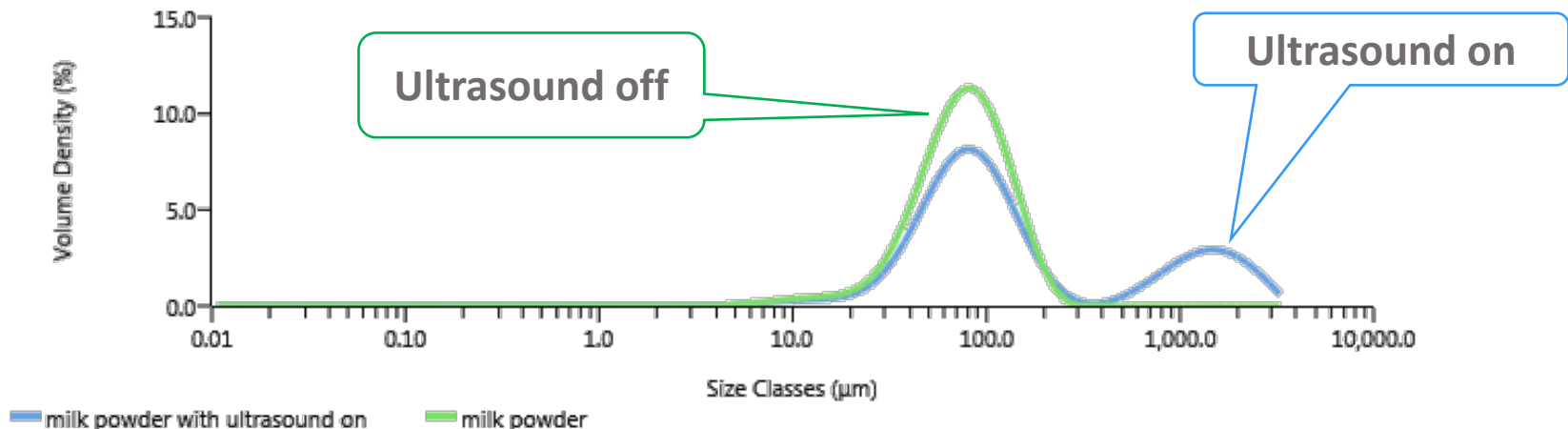
Unexpected large particles: Beam steering

- Ultrasound generates heat in the dispersant
 - Causing scattering signal on low angle detectors
 - Particularly in more volatile dispersants



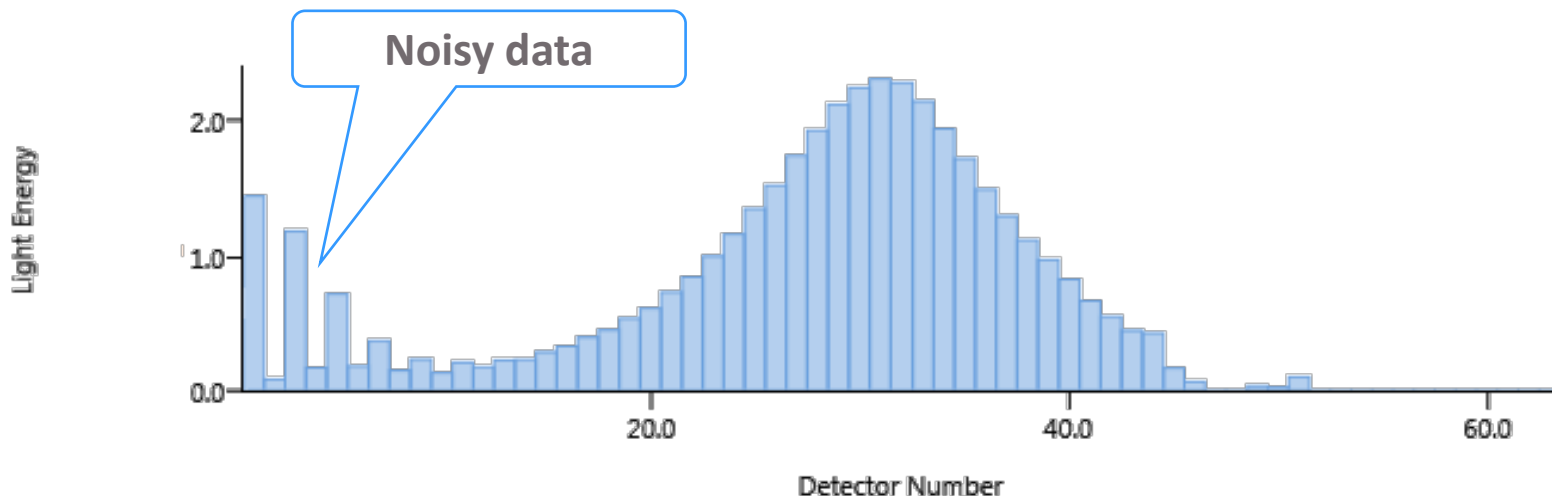
Unexpected large particles: Beam steering

- Ultrasound generates heat in the dispersant
 - Causing scattering signal on low angle detectors
 - Particularly in more volatile dispersants
 - This low angle scattering is interpreted as large particles
- Use a pre-measurement delay after ultrasound
 - Allows thermal gradients to dissipate



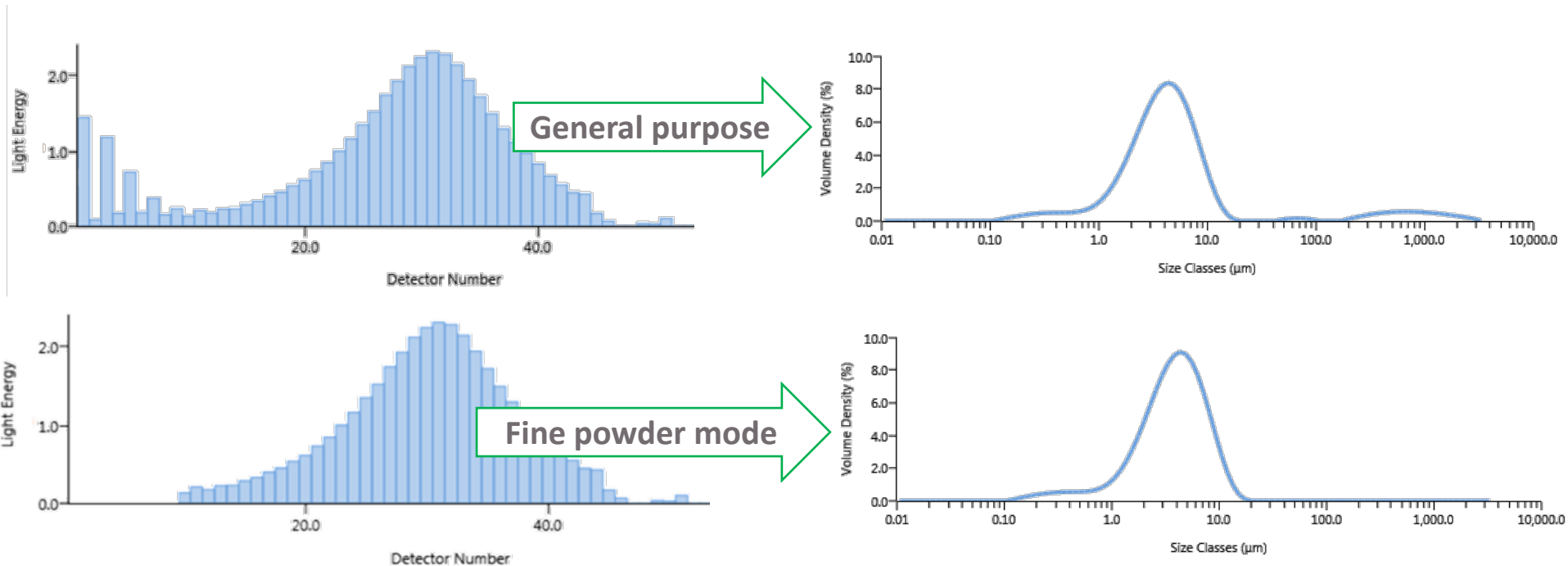
Unexpected large particles: Dry measurements

- Noise on the low angle detectors (1 to 10) can be significant when measuring fine particles dry
 - The noise is caused by thermal fluctuations in the air
 - This noise can be interpreted as large particles
- Measuring a longer background may help
 - Otherwise use the fine powder analysis mode.



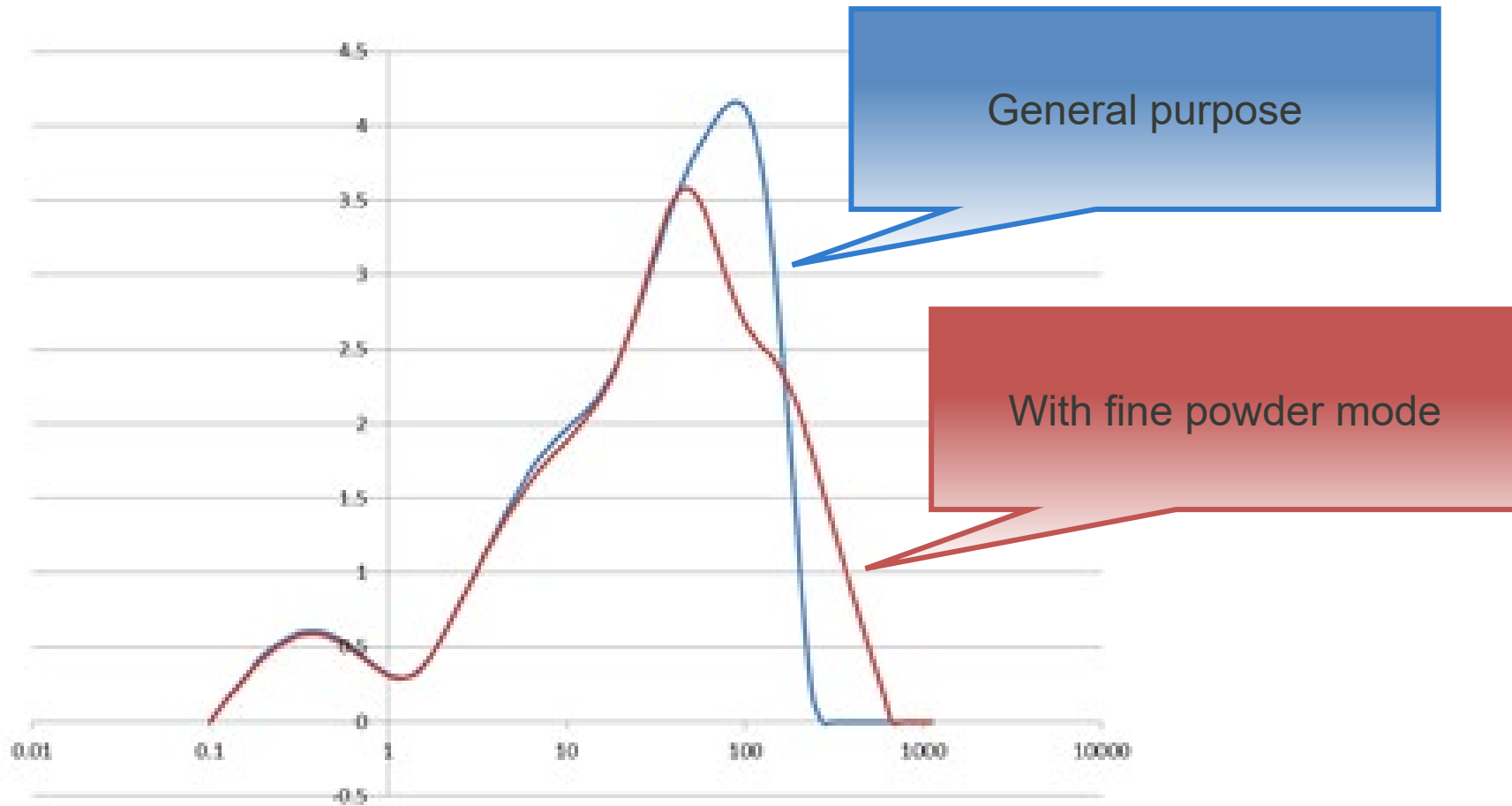
Dry measurements: Fine powder mode

- Fine powder mode removes the first 9 detectors
 - Eliminating the noise which may affect measurement of samples smaller than 10 micron
 - And removing the large modes that it produces
 - Limits the top end of the dynamic range to 600 micron



And when fine powder mode is not appropriate...

- ... strange results can be generated.





Further considerations for dry measurements:

- Check:
 - Are the cell windows clean?
 - Is the air pressure correct?
 - Does the air filter need changing?
 - Is there oil droplet contamination or moisture in the air supply?
 - Is the feeder earthed against static electricity ?
 - Is the vacuum bag full?
- **Is the sample flow even?**
 - if the sample obscuration is high, try lowering the feed rate or adjusting the hopper height
 - try changing the height of the hopper, different basket, ball bearing
 - try a different feed tray: often one tray design will deliver a more even sample flow for a particular material
- Use Fine Powder Mode when material less than 10 microns is present

Summary - data quality

- **Background data**

- Make sure that:
 - Material is not stuck to the cell windows
 - There is no dispersant contamination
 - There are no thermal gradients
 - That the system has been properly aligned

- **Sample data**

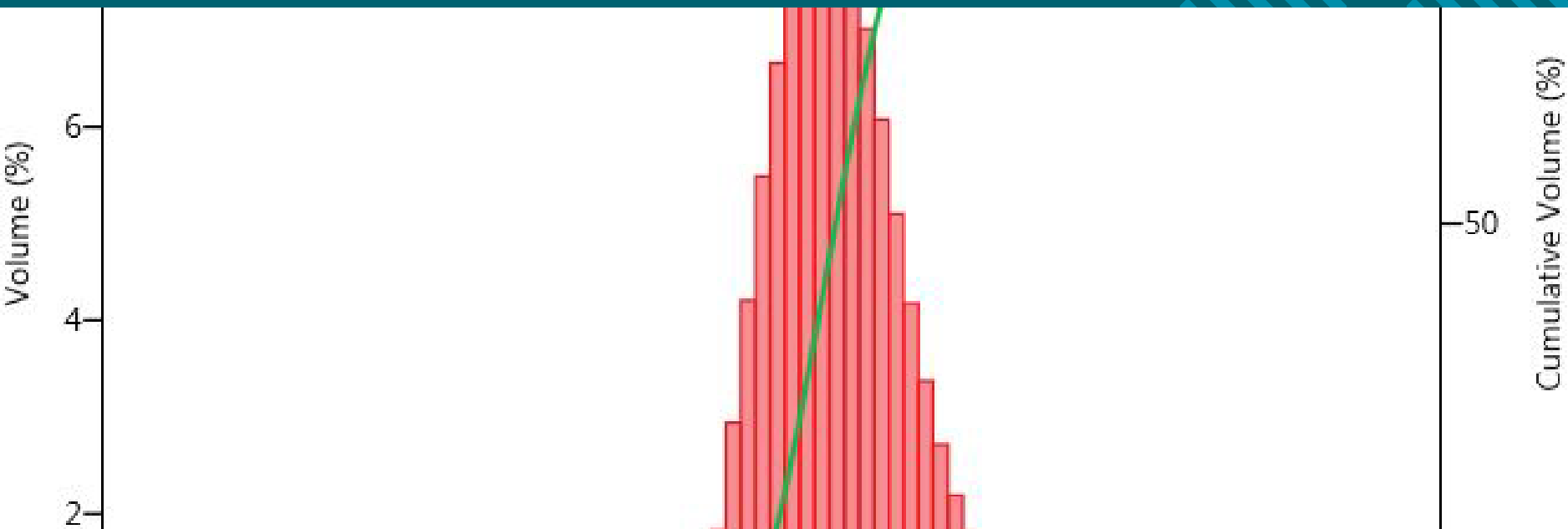
- Check that
 - There are reasonable signal to noise levels
 - There is no multiple scattering
 - There is no negative data
 - There is no noisy data
 - The inner detector data is free from castellation
 - There is no beam steering



**Malvern
Panalytical**

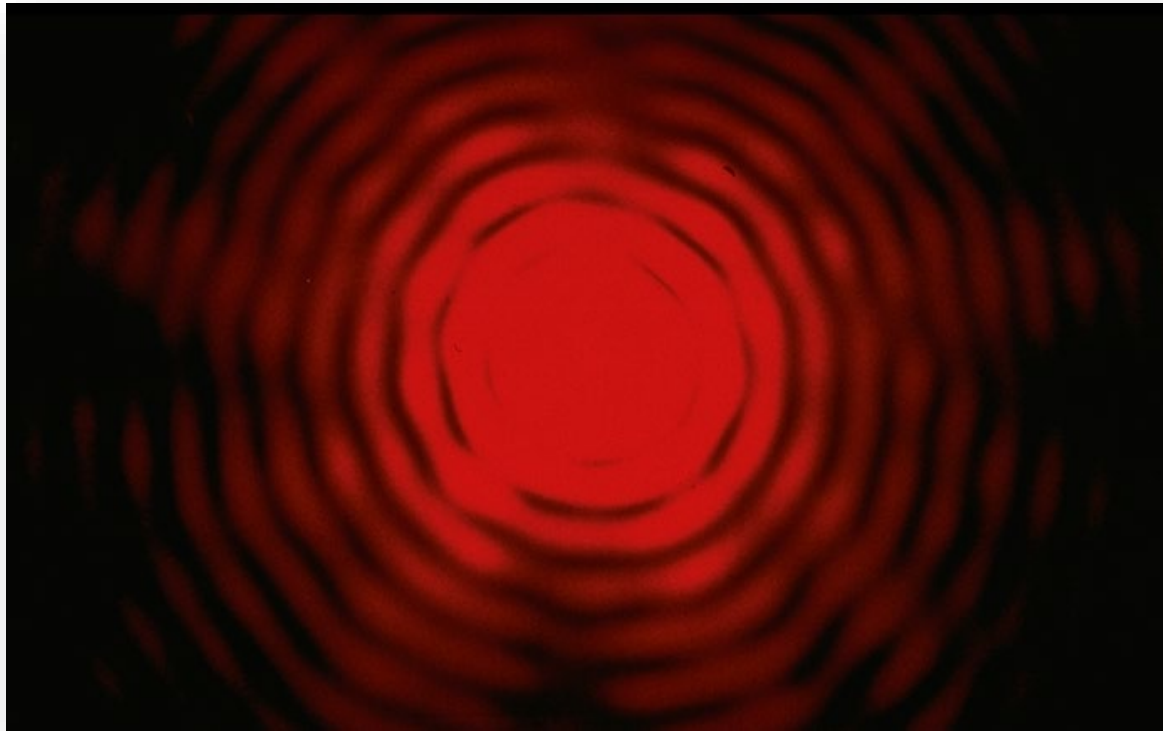
Mastersizer 3000 Part 2:

Obtaining and Understanding the Size Distribution



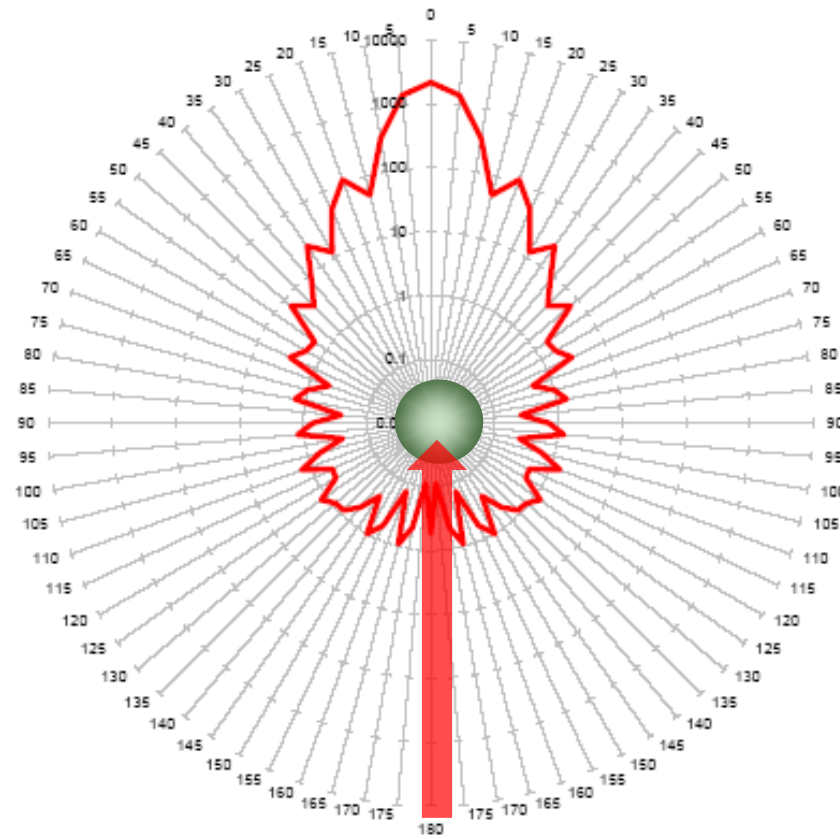
What does laser diffraction measure?

- Laser diffraction systems measure the scattering pattern produced by an ensemble of particles suspended in a laser beam



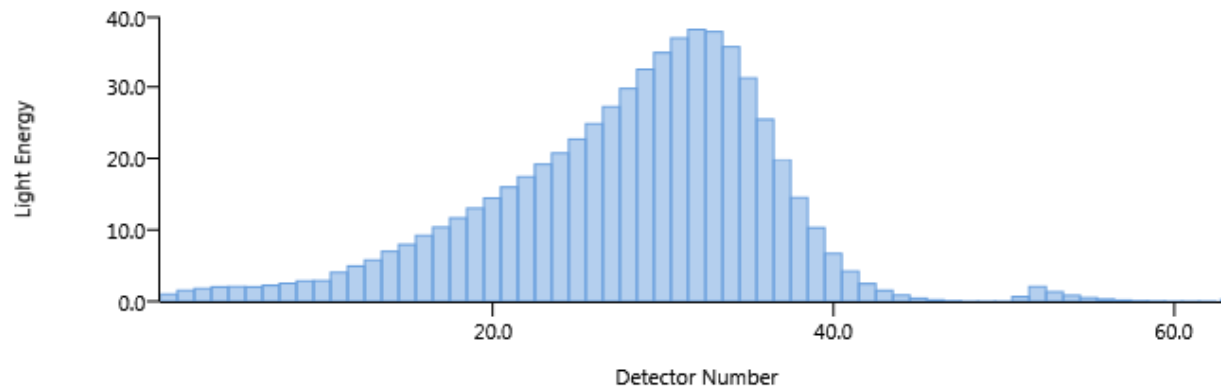
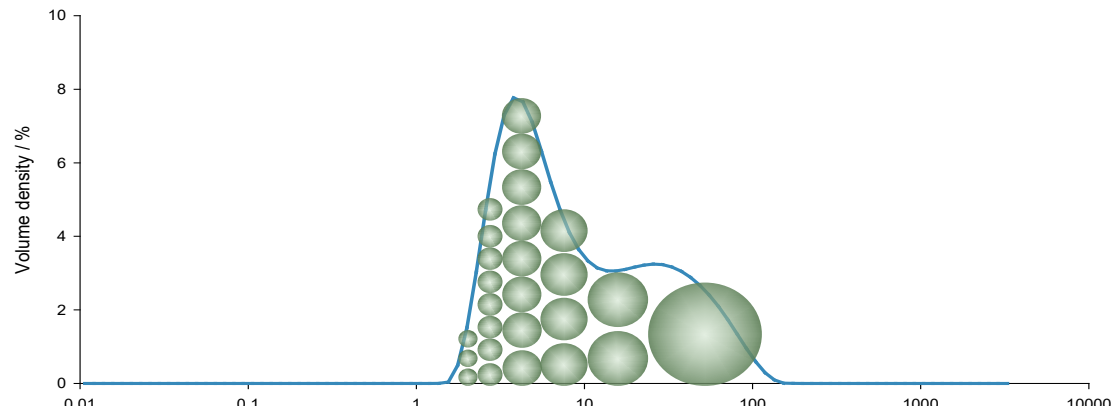
What does an optical model do?

- An optical model predicts the scattering pattern produced by a particle

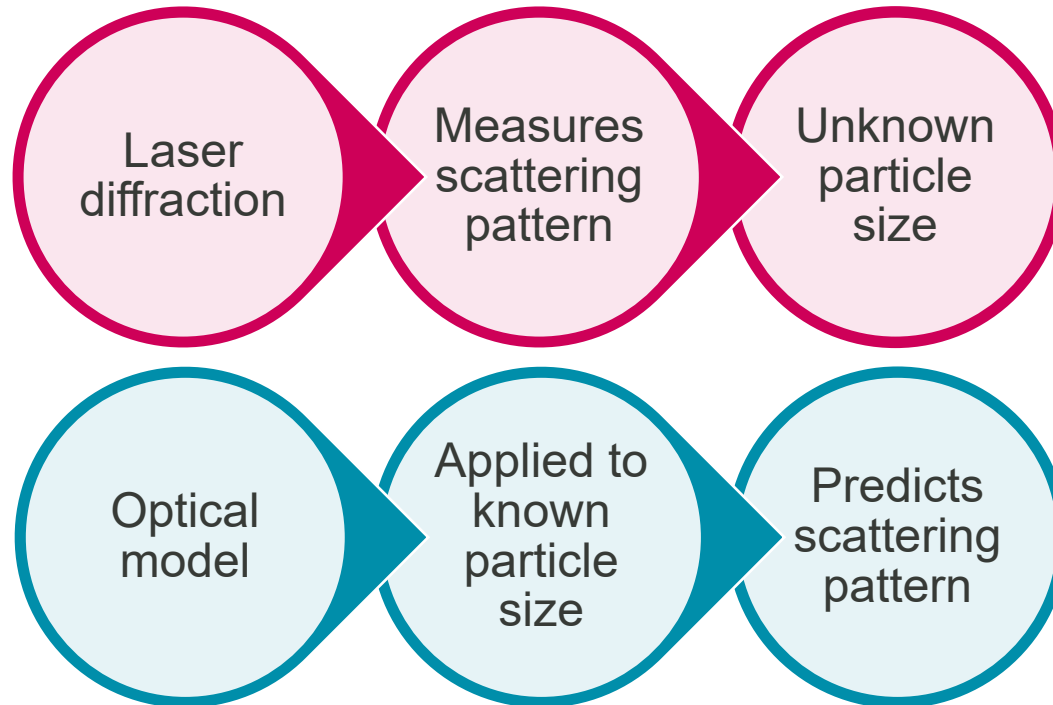


What does an optical model do?

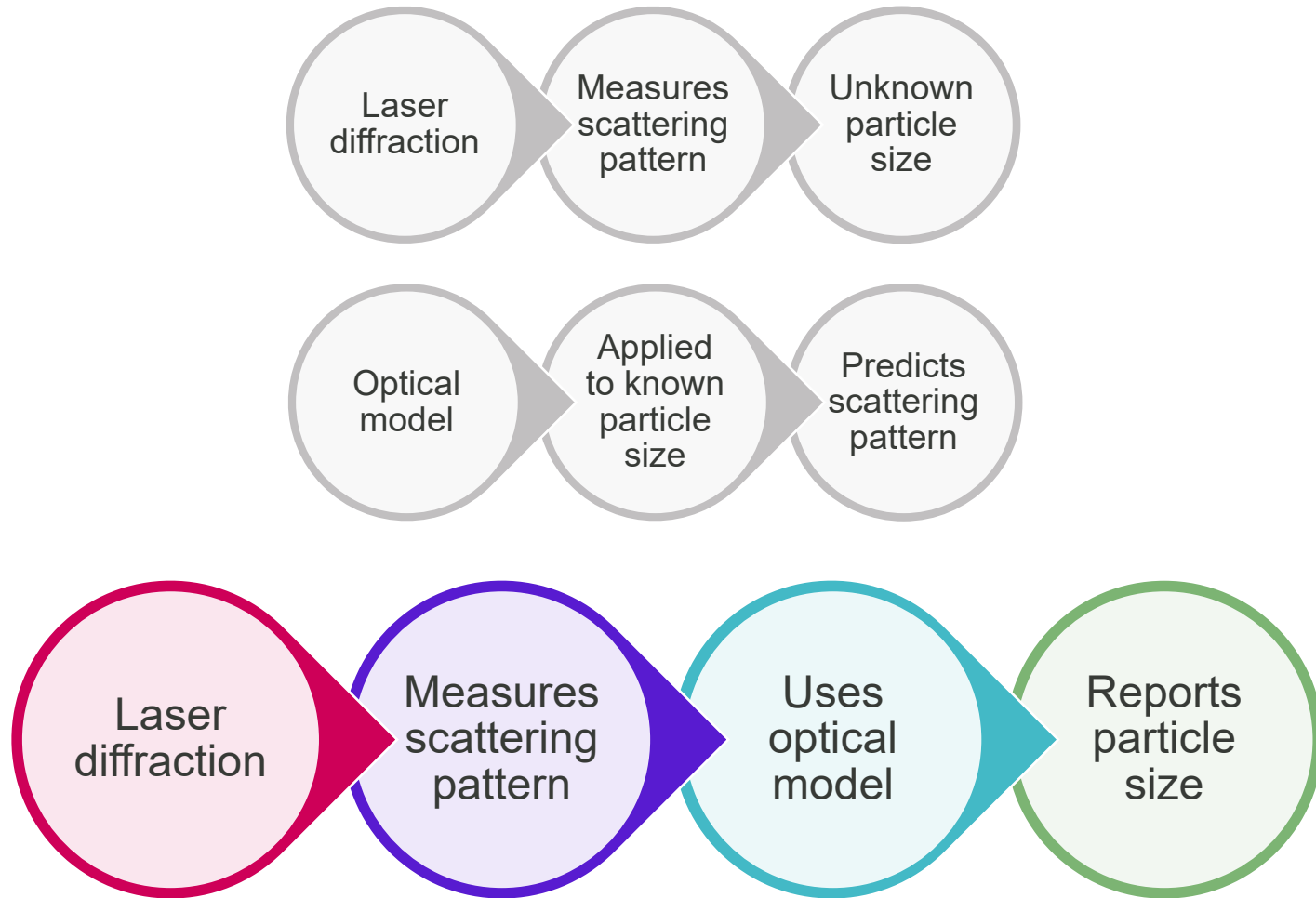
- And can therefore predict the scattering pattern produced by many particles



How do we use the optical model?



How do we use the optical model?

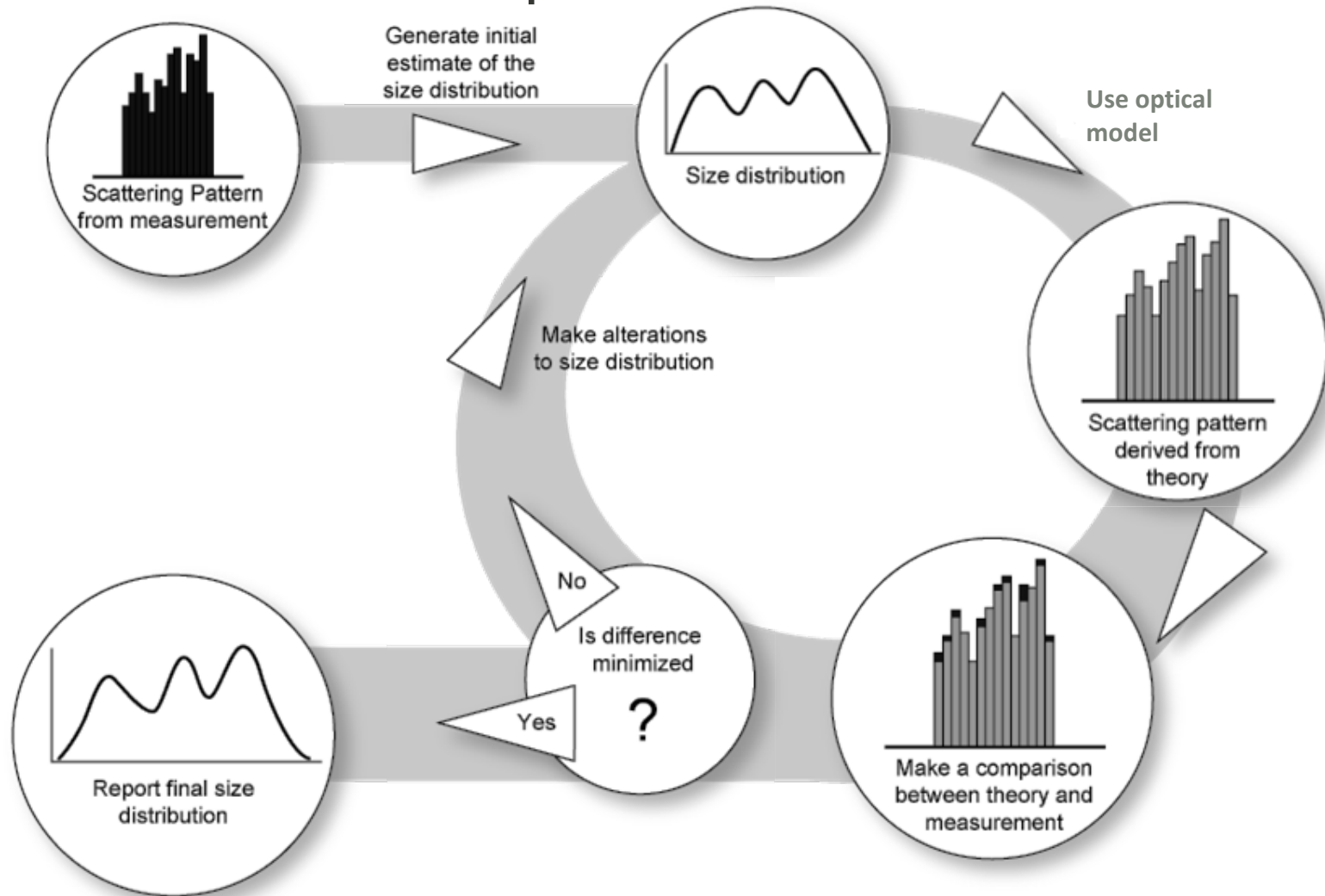


Understanding the data inversion process

Consider the following example:

- What is the answer to 6×7 ?
 - The answer to this is easy, it's 42!
 - Working this way through the problem gives one answer
- But what is 42?
 - 21×2
 - $40 + 2$
 - $126 \div 3$
 - $\sqrt{1764}$
 - We see that working backwards yields many possible solutions

How do we use the optical model?



Scattering models: Mie Theory

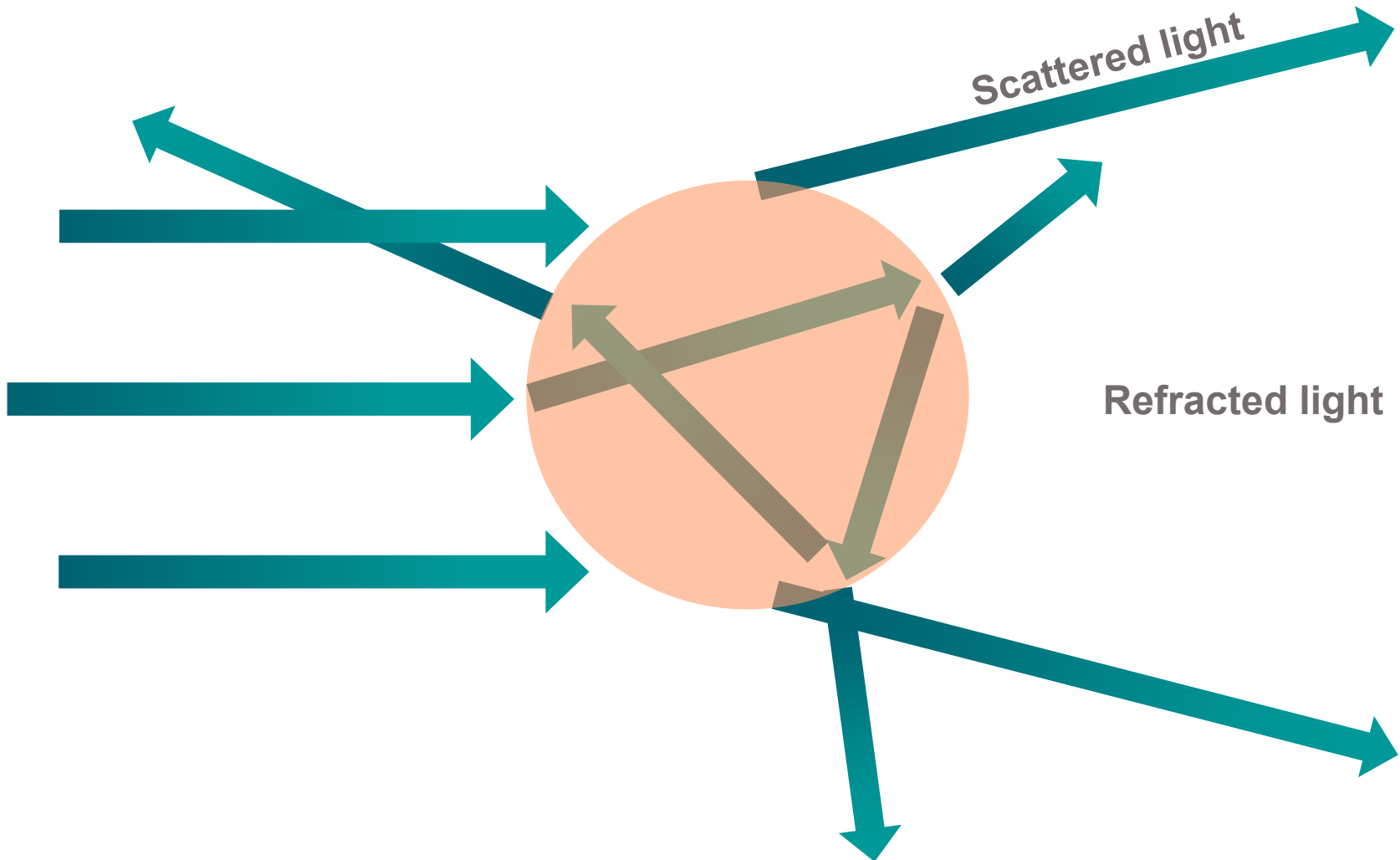
- Models the interaction of light with matter
 - Assuming that the particles are spherical
 - Assuming that it is a two phase system
- Valid for all wavelengths of light and all particle sizes
- Predicts the dependence of scattering intensity on particle size
- Predicts that secondary scattering is observed for small particles



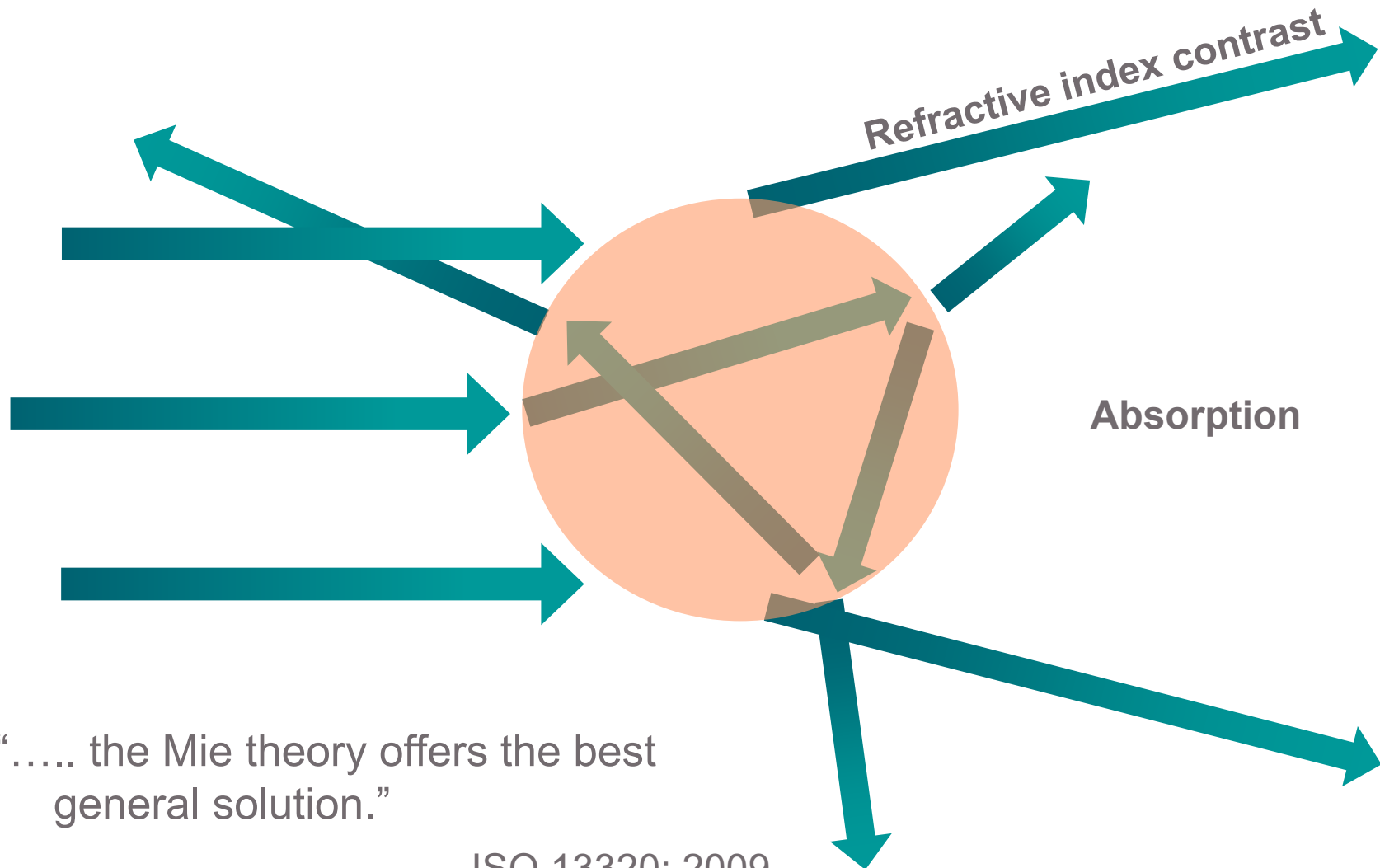
‘For particles smaller than about $50\mu\text{m}$ Mie theory offers the best general solution’

ISO13320

Mie Theory: Predicted scattering



Mie Theory: Optical properties



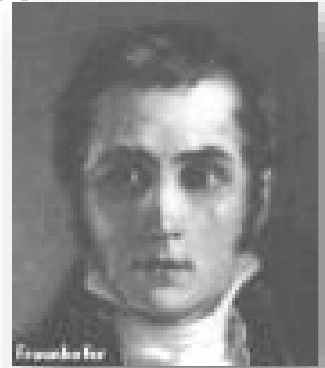
“..... the Mie theory offers the best general solution.”

ISO 13320: 2009



Scattering models: Fraunhofer Approximation

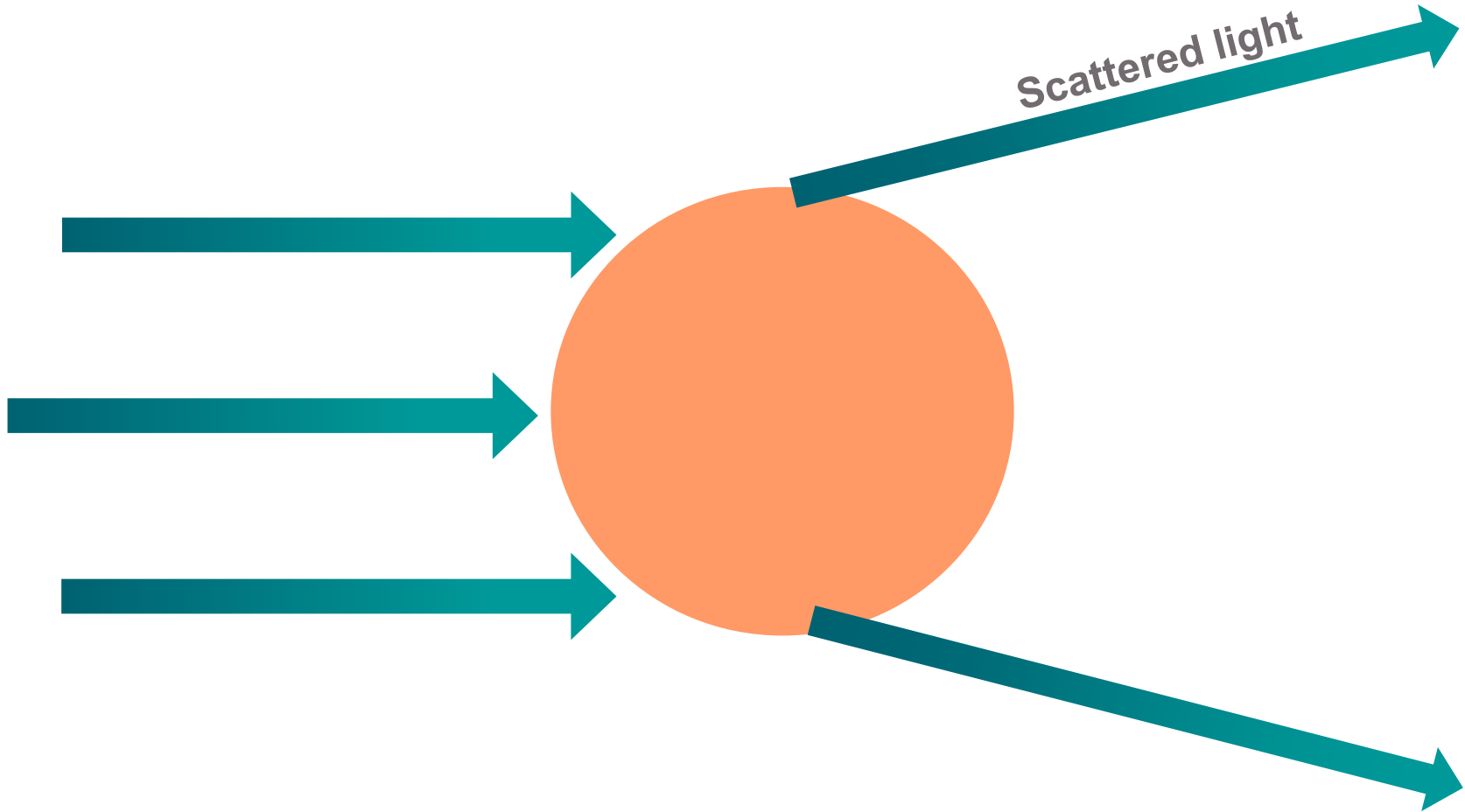
- Same basic assumptions as Mie Theory
 - Assuming that the particles are discs
 - Assuming that it is a two phase system
- Plus the additional assumptions that
 - The refractive index contrast is high ($RRI > 1.73$)
 - The particles are opaque
 - The wavelength of the light is much shorter than the particle size
 - The angle of refracted light is small
- In the MS3000 software the Fraunhofer approximation is available as a particle type.



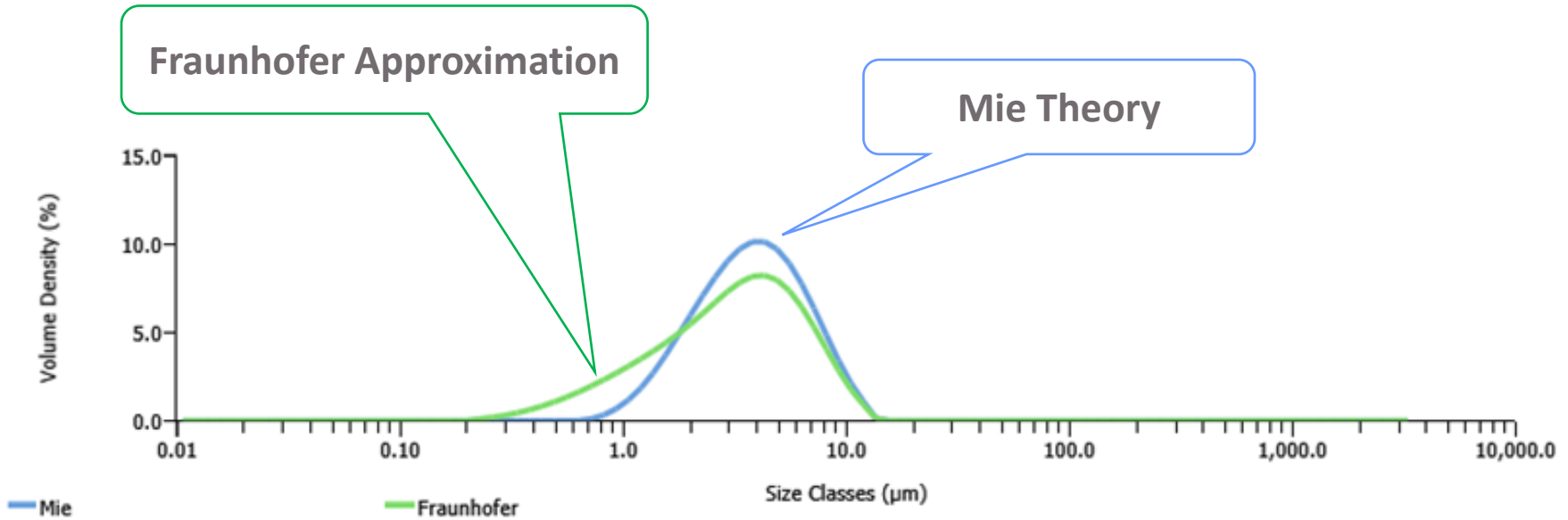
‘The advantage of this equation is that it is relatively simple and quick to calculate’. ‘The Fraunhofer approximation does not make use of any knowledge of the optical properties of the material’.

ISO13320

Fraunhofer Approximation: Predicted scattering



Comparing the results of the scattering models



'If the Fraunhofer approximation is applied for samples containing an appreciable amount of small, transparent particles, a significantly larger amount of small particles may be calculated.'

ISO13320

Scattering models - Fraunhofer Approximation

- **Disadvantages**

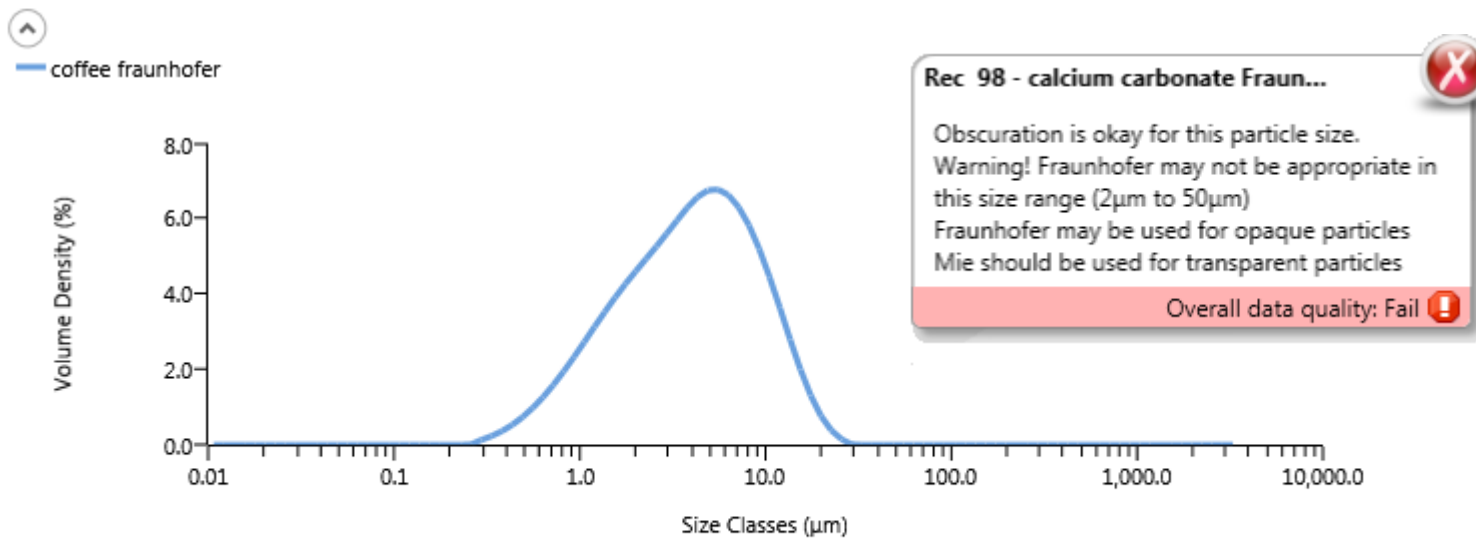
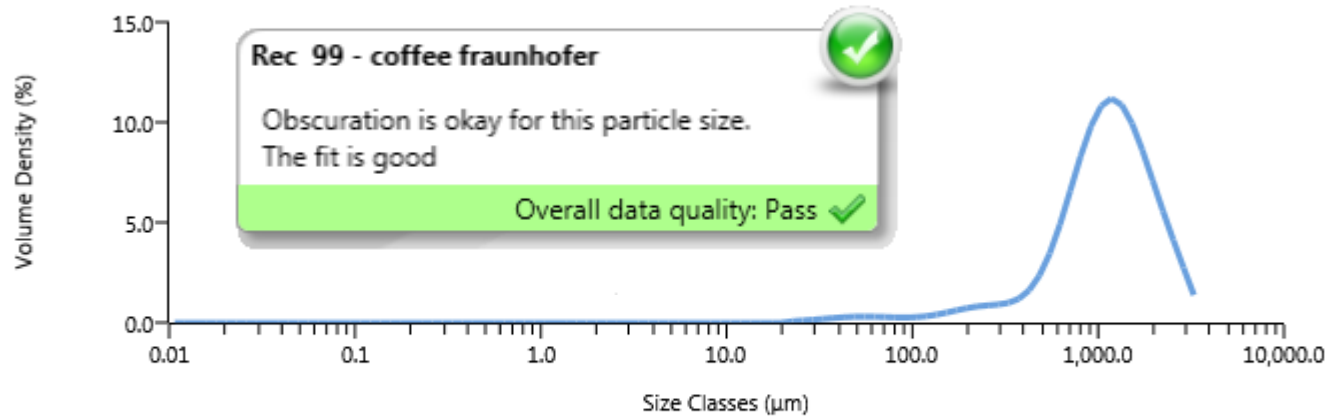
Will produce incorrect answers when...

- Particles are $<50\mu\text{m}$
- The angle of scatter becomes large and secondary scattering occurs.
- The relative refractive index is small (<1.3) - this equates to a particle refractive index of 1.73 in water.

- **Claimed Advantage**

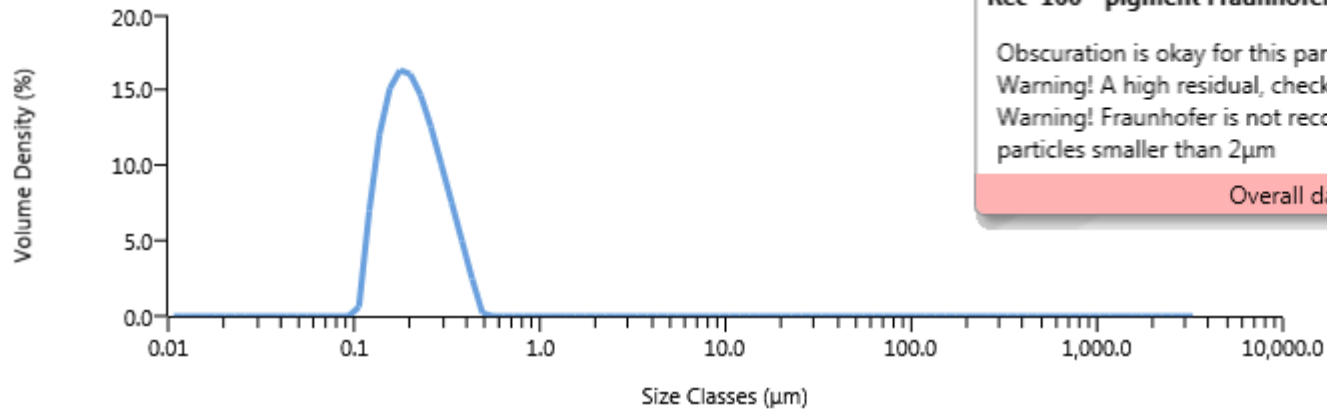
- “No need to know the *optical properties* of your material.”

Mie vs Fraunhofer: Data quality advice



calcium carbonate Fraunhofer

Mie vs Fraunhofer: Data Quality advice



Rec 100 - pigment Fraunhofer

Obscuration is okay for this particle size.
Warning! A high residual, check your fit.
Warning! Fraunhofer is not recommended for
particles smaller than 2µm

Overall data quality: Fail



— pigment Fraunhofer

Which optical properties do we need?

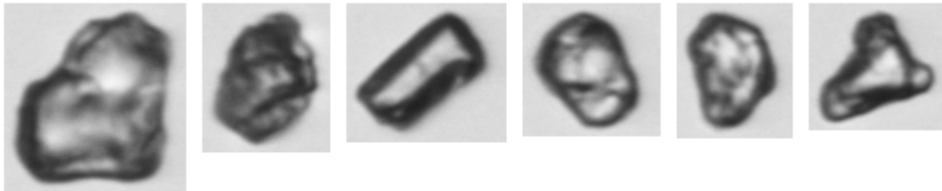
- To use Mie theory correctly we need to know three optical properties:
 - The refractive index of the dispersant
 - The refractive index of the sample material
 - The imaginary part of the refractive index of the sample material
 - Often referred to as the absorption

‘Good understanding of the influence of the complex refractive index in the light scattering from particles is strongly advised in order to apply the Mie theory or the Fraunhofer approximation appropriately.’

ISO13320


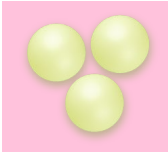



The imaginary refractive index (or absorption)

- The absorption can be determined by looking at the dispersed sample under a microscope and observing its
 - Shape
 - Transparency
 - Internal structure
- Absorption is generally required to a factor of 10
 - E.g. 0.1 or 0.01 (not 0.023)



Images of some calcium carbonate particles, an absorption of 0.01 would be used for these particles.

Estimating absorption from particle appearance

| Appearance | Absorption | Example |
|---|------------|--|
|  | 0 | Latices |
|  | 0.001 | Emulsions |
|  | 0.01 | Crystalline milled powders |
|  | 0.1 | Slightly colored powders |
|  | 1.0+ | Highly colored (complementary) and metal powders |

Methods for determining the refractive index

- Four main routes to refractive index information
 - Reference books and the internet
 - Appendix of ISO 13320
 - Malvern materials database
 - CRC handbook
 - Manufacturers label (for dispersants)
 - Online info
 - Refractometer measurements
 - Microscope observations
 - Empirical/semi-empirical models



Online sources for refractive index information

- www.Luxpop.com

Index of Refraction: fixed and variable ratio at given wavelength

Different presentation approaches: (A) lists all refractive index values (n or k) from 11 to 12 and (B) gives individual refractive index values, along with more information ranging from bandgaps, to temperature dependencies, to fabrication information, depending on the material. (C) gives refractive index for variable atomic content compound materials. [Explanation of materials and references](#)

(A) GaAs I1: 400 nm I2: 700 nm step: 50 nm n · go

(B) Ag (Silver) I: 1047 nm temp: 25 deg C go

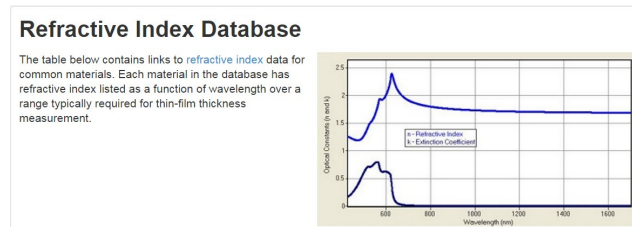
(C): Al(x)Ga(1-x)As I: 1047 nm temp: 25 deg C x: 0.07 go

* See also our [Long List of Index of Refraction Values \(A-Z\)](#), for other materials.
* [Click here for more index of refraction terminology](#) and [variable compound discussion points](#).

- RefractiveIndex.info



- www.Filmetrics.com/refractive-index-database



- Google Scholar



- Remember to look for a wavelength of 633nm for the red light and 470nm for the blue light.

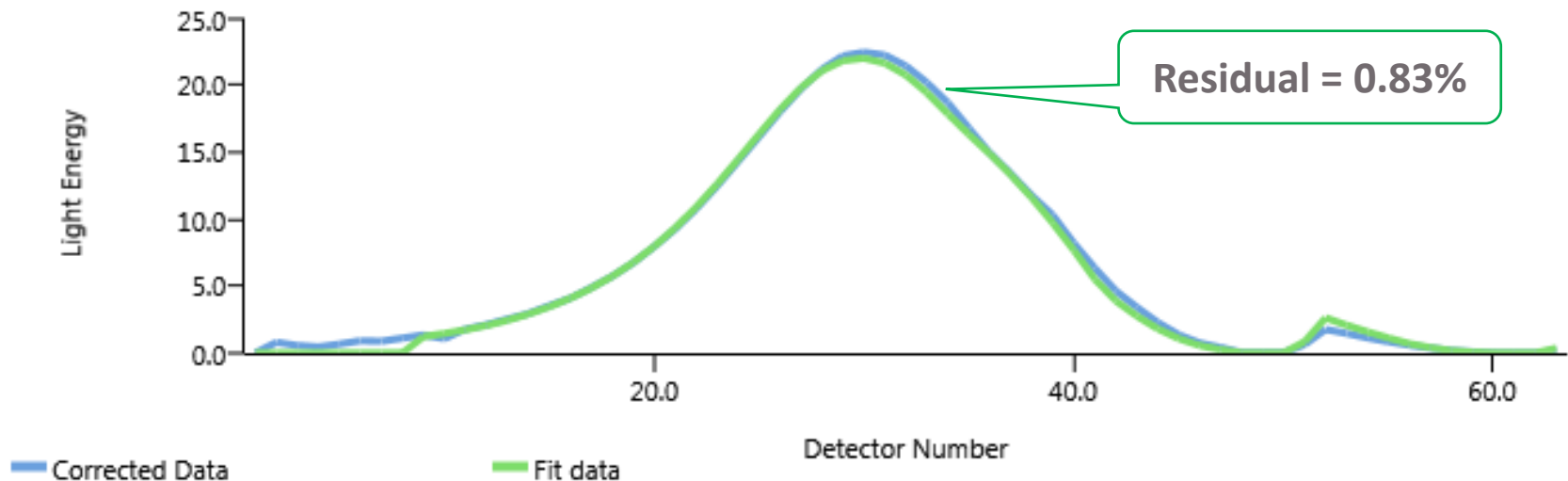
Choosing the refractive index



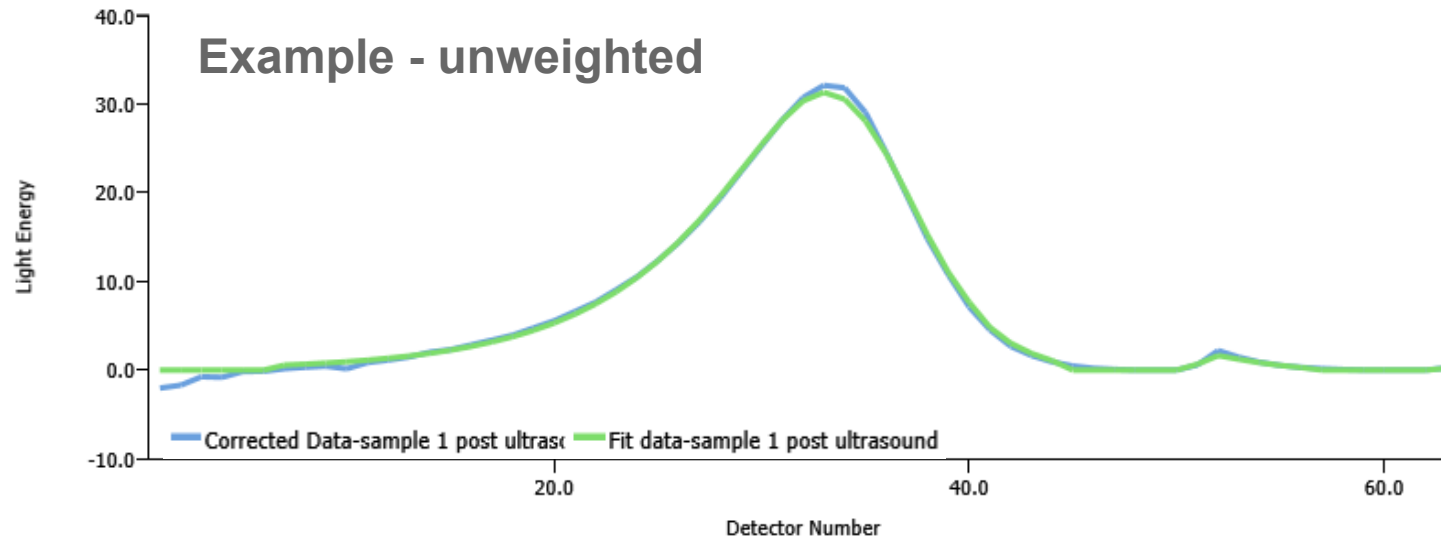
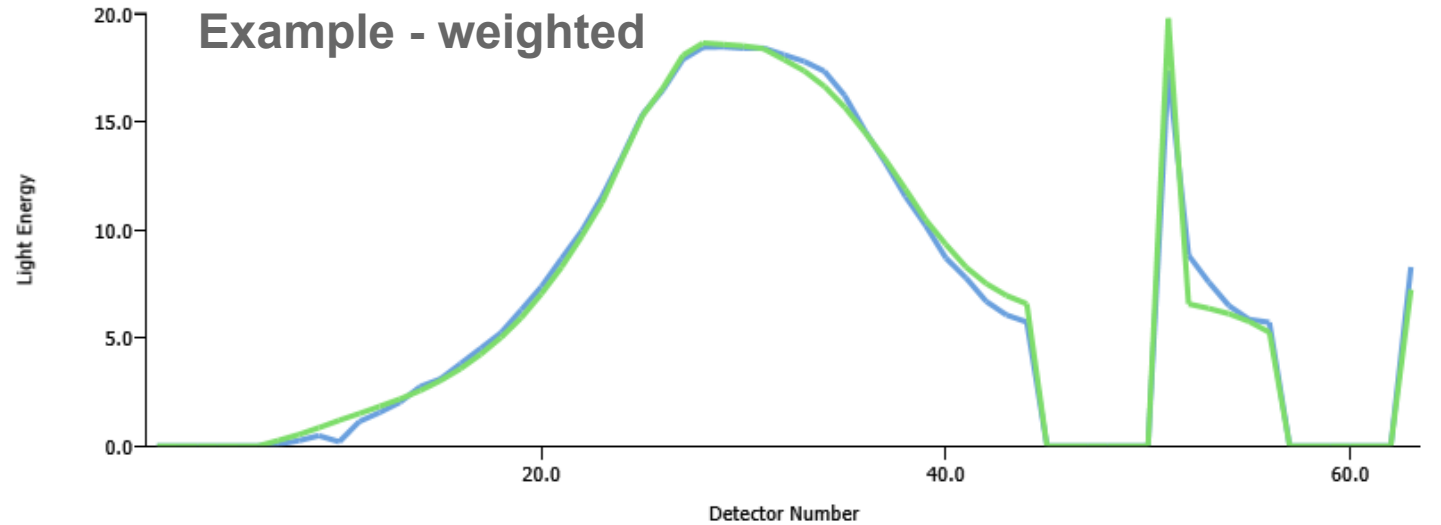
- You can estimate the optical properties based on typical values of similar materials.
 - A Refractive Index is generally only required to 2 decimal places e.g. 1.42 not 1.4234
- Some families of similar materials are:
 - Plastics and elastomers = 1.38 - 1.57
 - Organic compounds = 1.4 - 1.7
 - Inorganic salts = 1.52 - 1.8
 - Metal Oxides = 1.6 - 2.5
- Use the estimated refractive index as a starting point and examine the fit to confirm the suitability of the value chosen.

Assessing the data fit

- The fit report shows the measured and calculated scattering data
- How well these overlay is known as the data fit
- The residual quantifies how good the fit is
 - Residual = area between the two curves

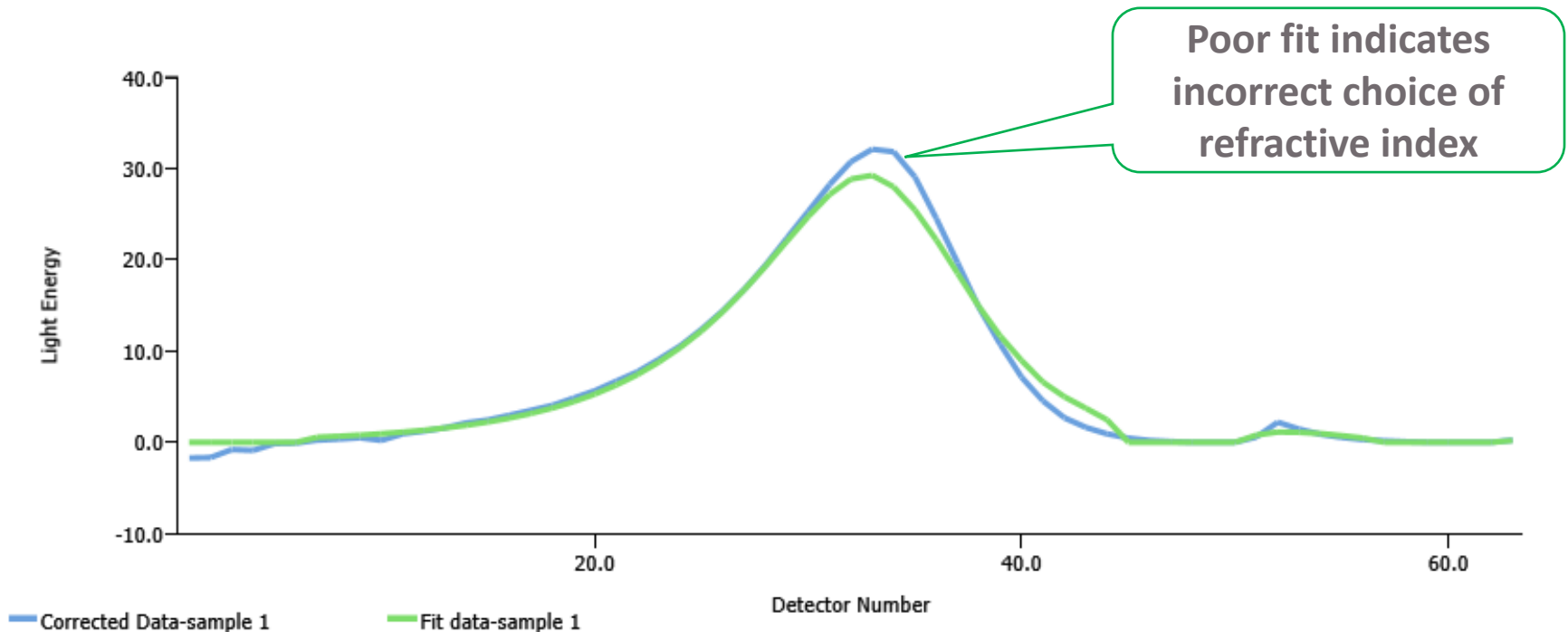


Weighted and un-weighted data fits



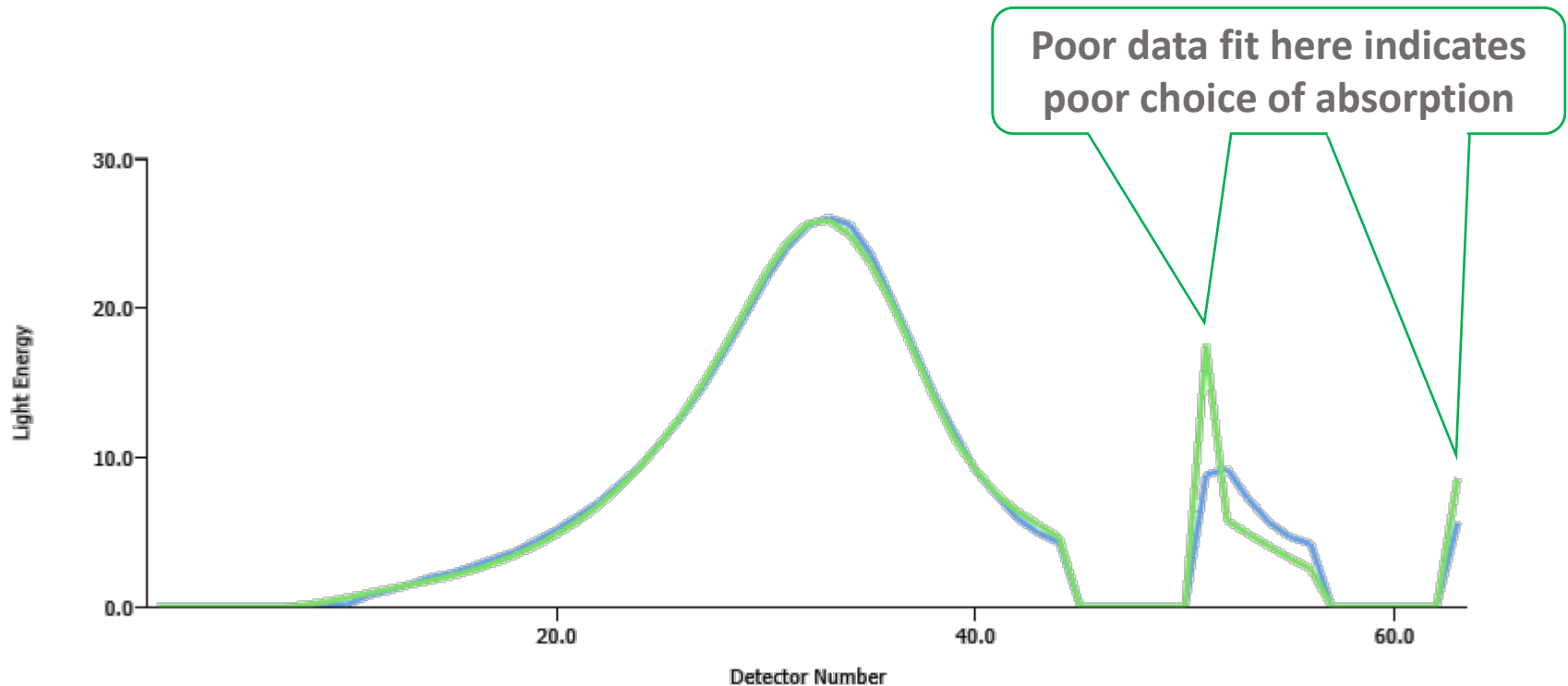
Inspecting the data fit: refractive index

- A poor fit to the focal plane detectors (< 40) suggests an incorrect choice of refractive index



Inspecting the data fit: absorption index

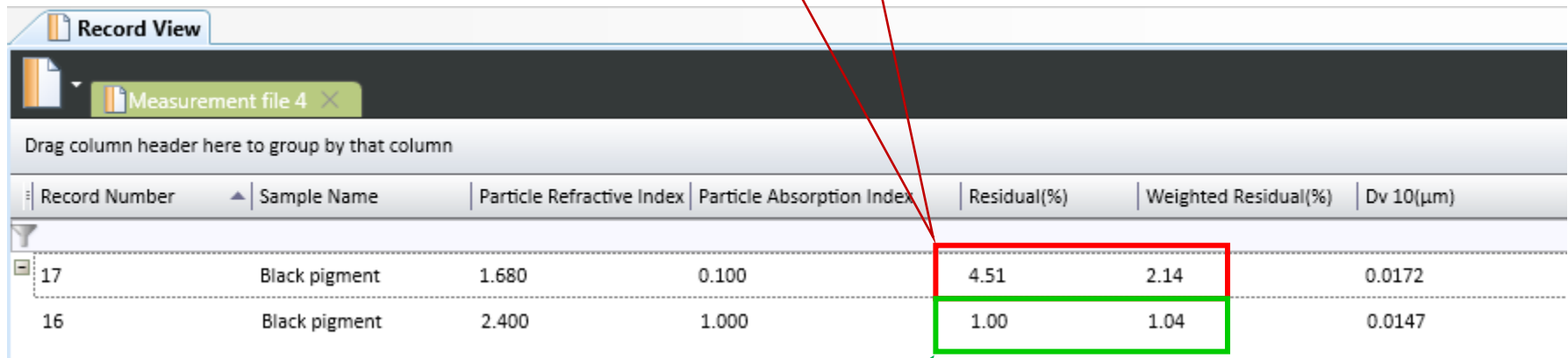
- Misfits to the extinction detectors indicate an incorrect absorption value
 - 51 in the red light
 - 63 in the blue light



Residuals - good “rule of thumb”

- For most size distributions, the residual should typically be less than 1%
- For a **good** fit, the weighted and unweighted residuals should be of similar orders of magnitude
- A low residual is good, but the distribution should be believable

Poor agreement

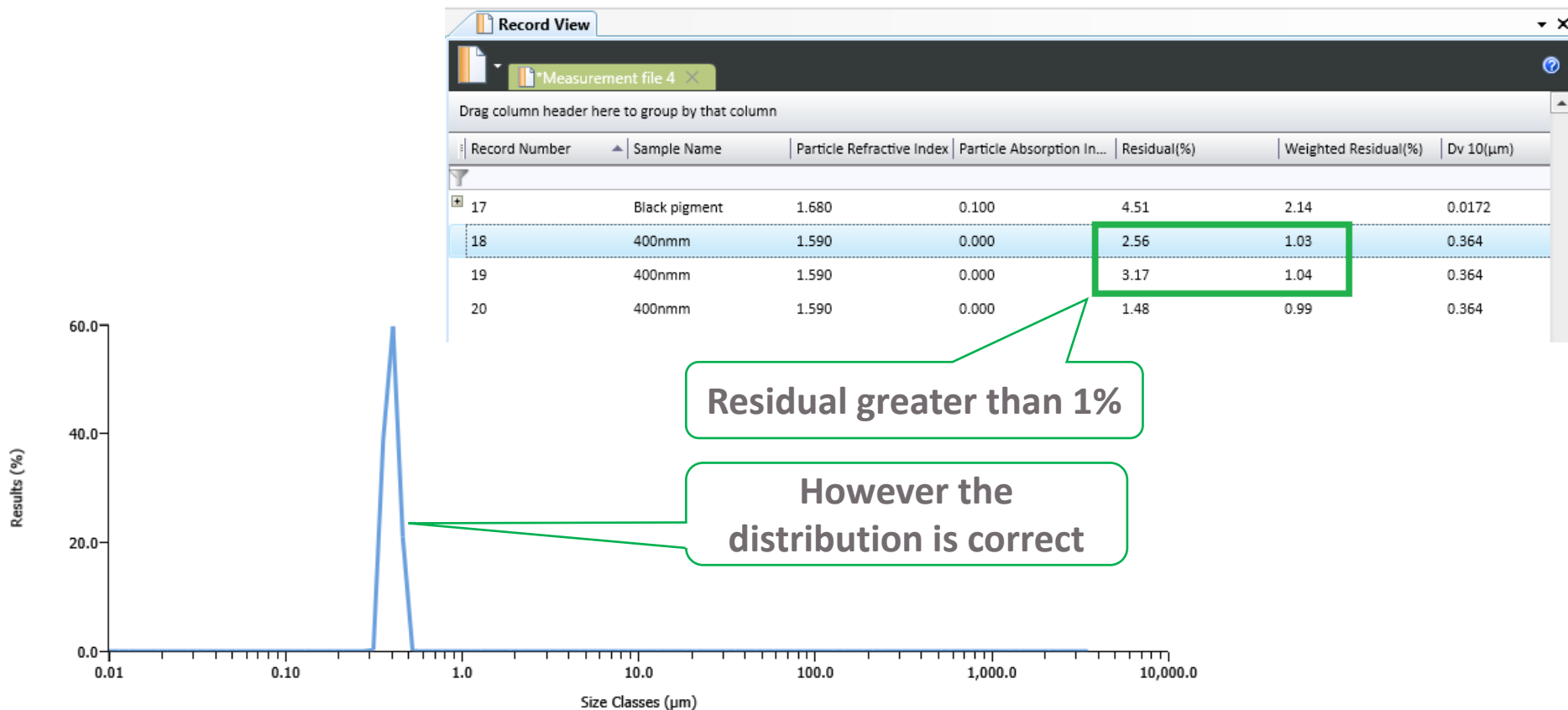


| Record Number | Sample Name | Particle Refractive Index | Particle Absorption Index | Residual(%) | Weighted Residual(%) | Dv 10(μm) |
|---------------|---------------|---------------------------|---------------------------|-------------|----------------------|-----------|
| 17 | Black pigment | 1.680 | 0.100 | 4.51 | 2.14 | 0.0172 |
| 16 | Black pigment | 2.400 | 1.000 | 1.00 | 1.04 | 0.0147 |

Good agreement

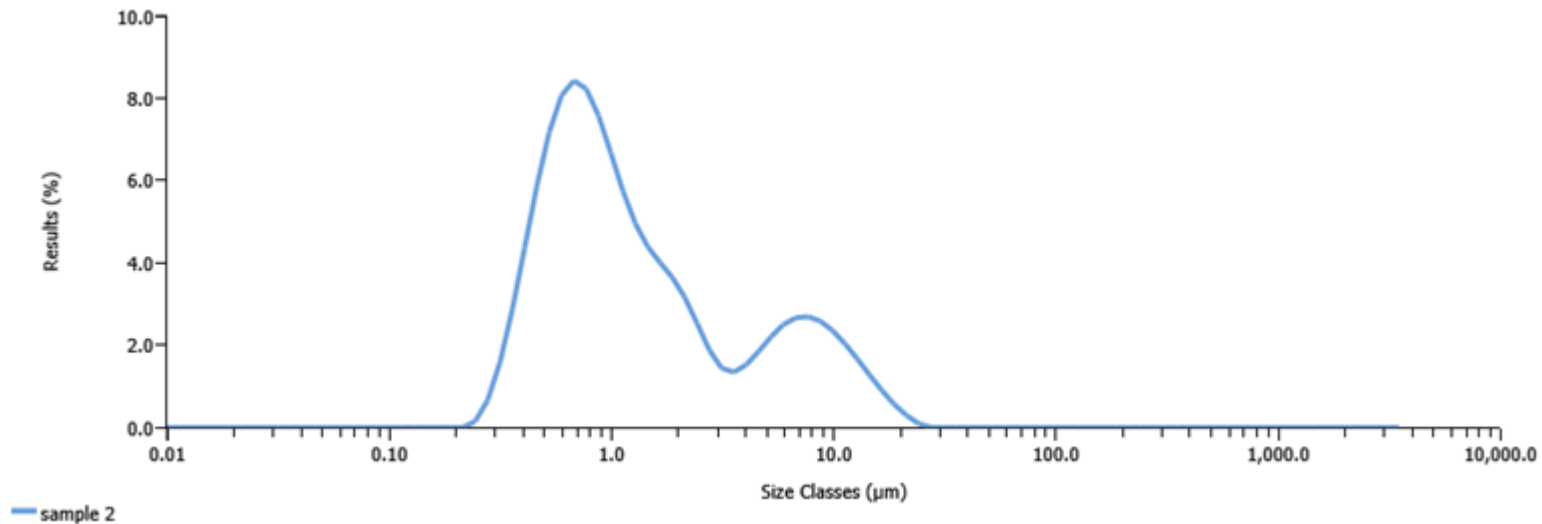
Residuals for very narrow distributions

- Residuals less than 1.0 are not always achievable.
- Narrow distributions can give higher residuals than expected



Assessing the data fit: Example

- The user is seeing an “unexpected” mode of small material.
- The optical properties used were:
 - RI:1.4, Absorption: 0.01





Assessing the fit using 1.4/0.01

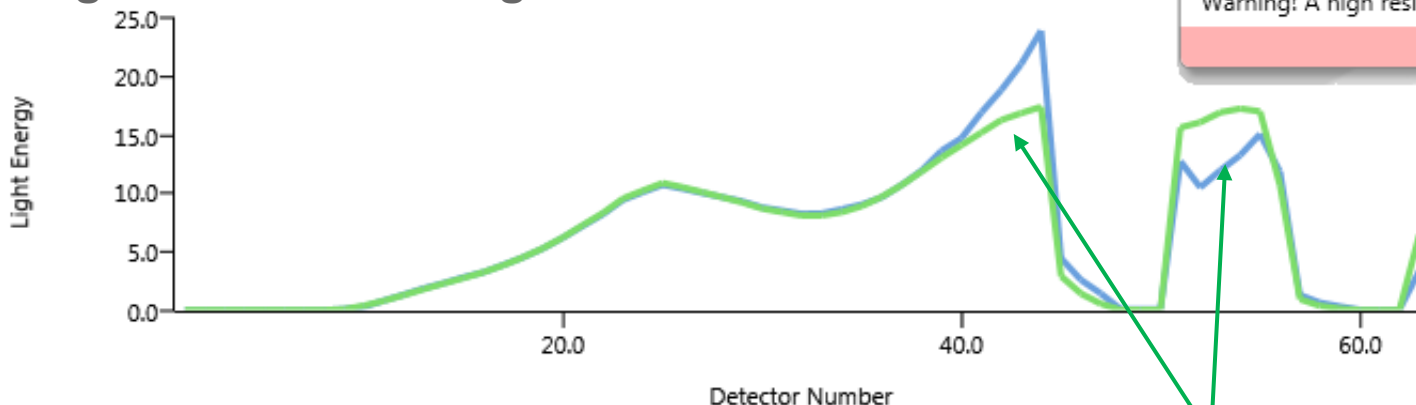
Rec 2 - calcium carbonate

Obscuration is okay for this particle size.
Warning! A high residual, check your fit.

Overall data quality: Fail

Weighted fit

Weighted Residual = 3.26

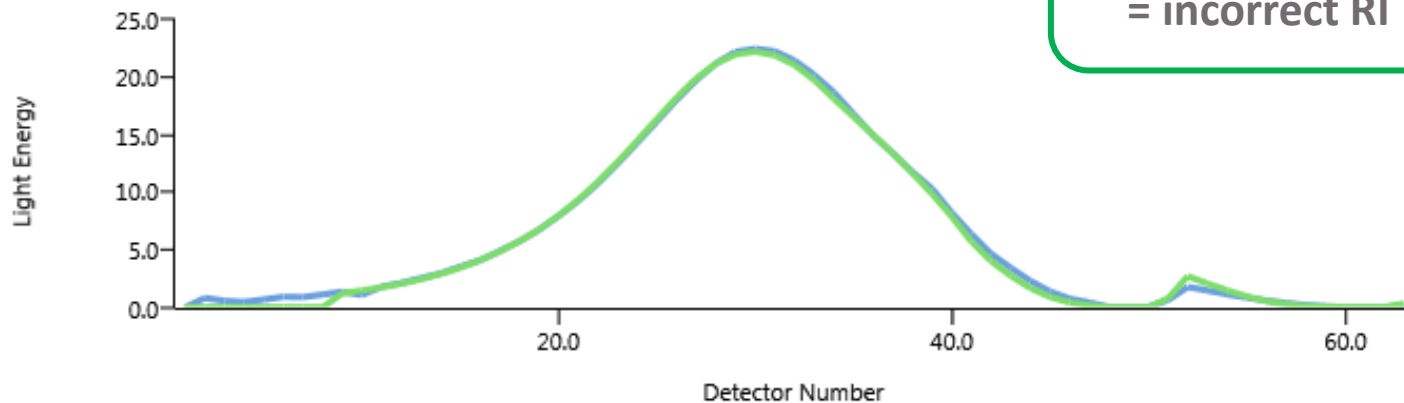


Weighted Data-calcium carbonate Fit data(weighted)-calcium carbonate

Poor fit
= incorrect RI

Un-Weighted fit

Residual = 0.82




Corrected Data-calcium carbonate Fit data-calcium carbonate




Assessing the fit using 1.54/0.01

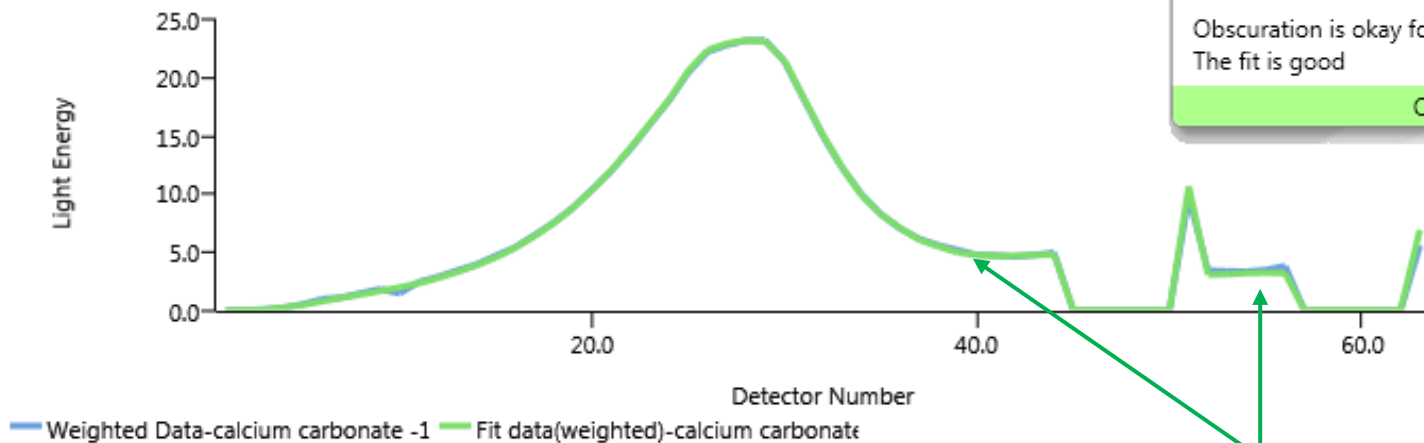
Weighted fit

Weighted residual = 0.48

Rec 23 - calcium carbonate 

Obscuration is okay for this particle size.
The fit is good

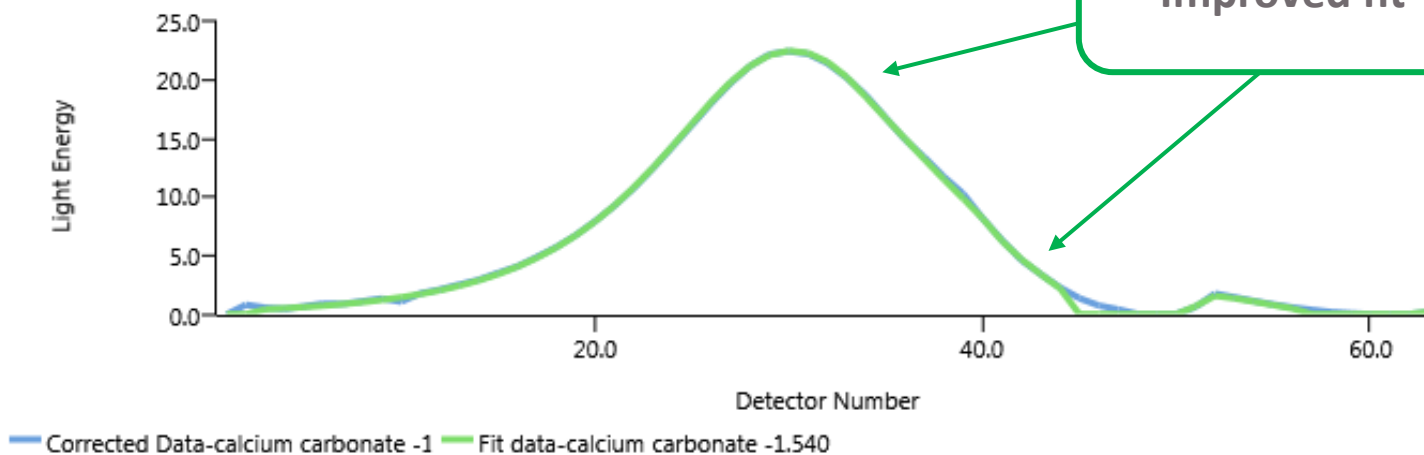
Overall data quality: Pass 



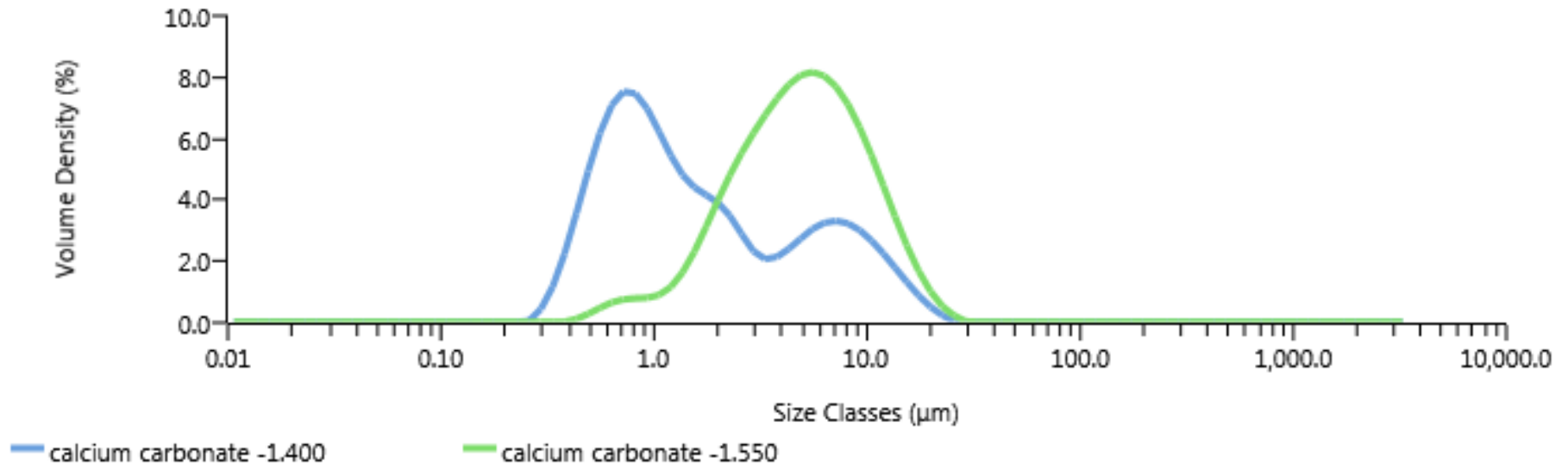
Improved fit

Un-Weighted fit

Residual = 0.57



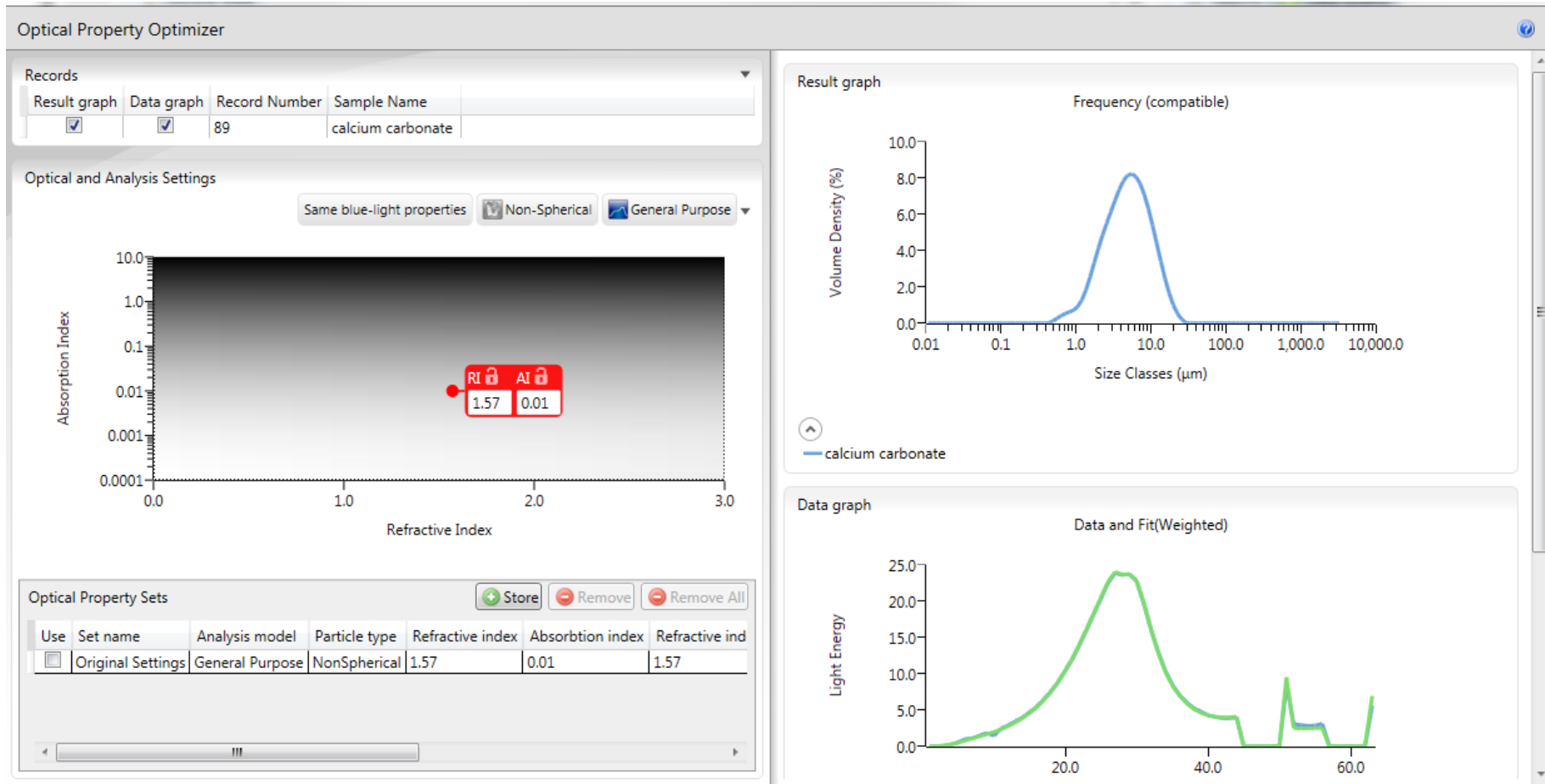
Looking at the results



- Sample is calcium carbonate
 - Reference RI is between 1.53 and 1.63

The optical property optimiser (OPO)

- Offers a quick way to adjust optical properties and assess the fit and result



Setting optical properties of mixtures

- Sensitivity to optical properties increases as particle size decreases.

Option 1 Consider the smallest

- Set the optical properties to suit the finest particles.

Option 2 Consider the average

- Calculate the volume-weighted average optical properties.
- Accept there may be errors in the fines.

Option 3 Consider using Fraunhofer

- Use the Fraunhofer Approximation.
- Accept there may be errors in the fines.

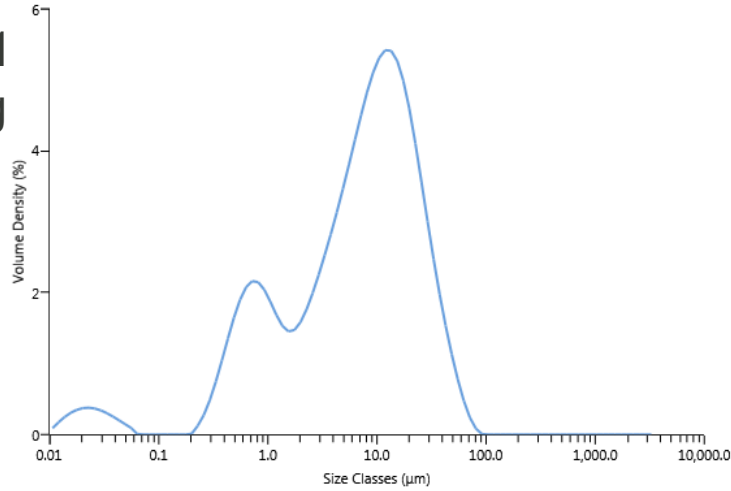
Plastisol – a real life example

- Plastisol uses include:
 - Slush moulding
 - Textile ink
 - Binder in Electric Solid Propellant (rocket fuel)
- Plastisol is a suspension of PVC particles in a liquid plasticizer:

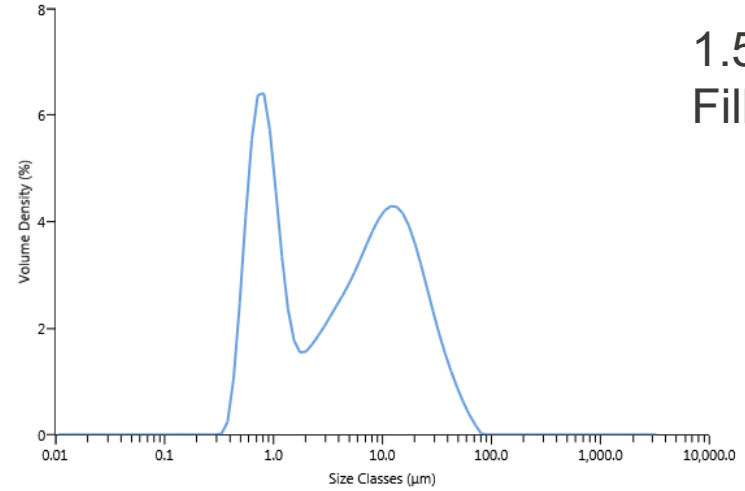
| | Refractive index | Absorption |
|--------------------|------------------|------------|
| Blowing Agent | 1.76 | 0.1 |
| Filler | 1.59 | 0.1 |
| Pigment | 2.51 | 0.01 |
| Polyvinyl chloride | 1.54 | 0.01 |

Plastisol – by optical properties

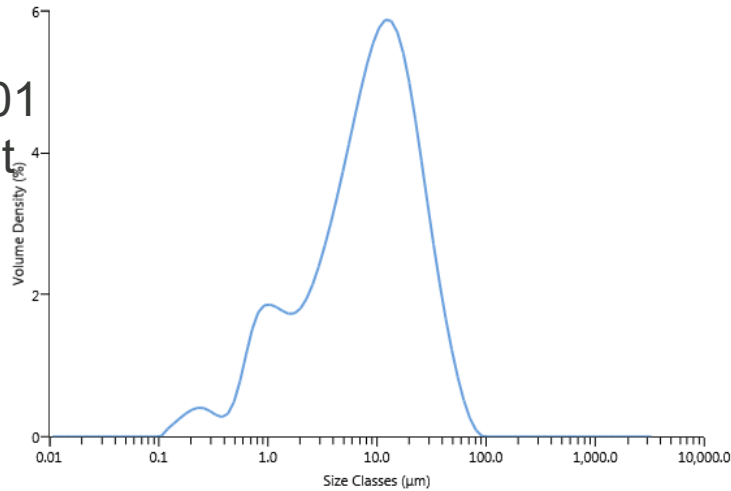
1.76/0.1
Blowing
agent



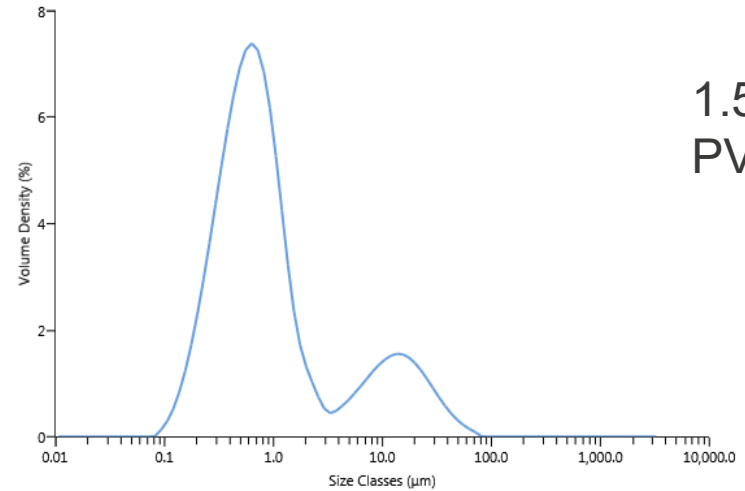
1.59/0.1
Filler



2.51/0.01
Pigment



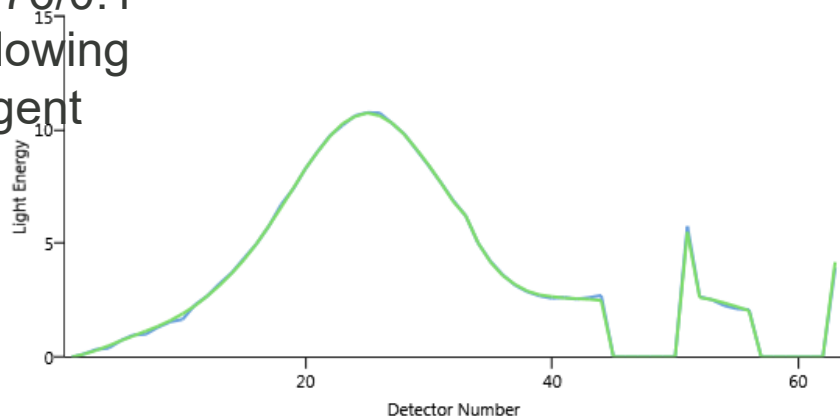
1.54/0.01
PVC



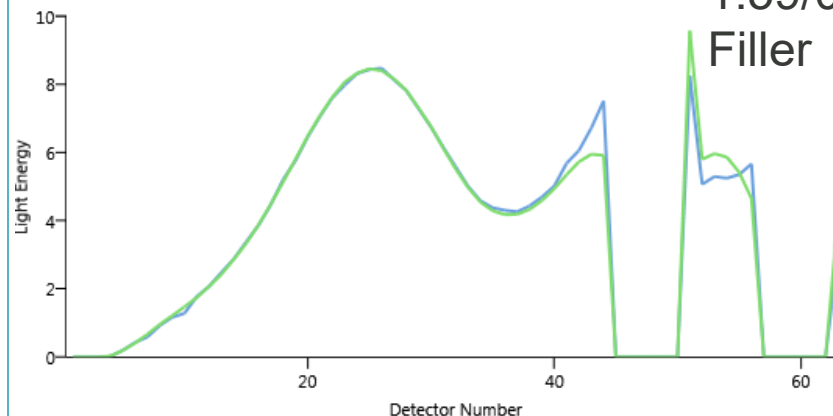


Plastisol – by fit and residual

1.76/0.1
Blowing
agent

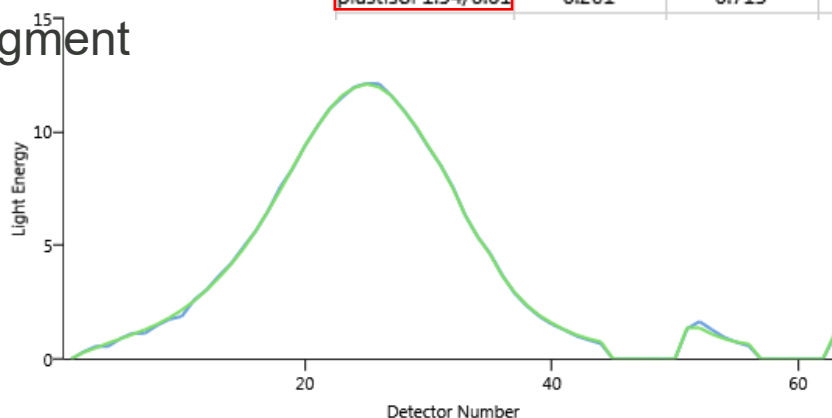


1.59/0.1
Filler

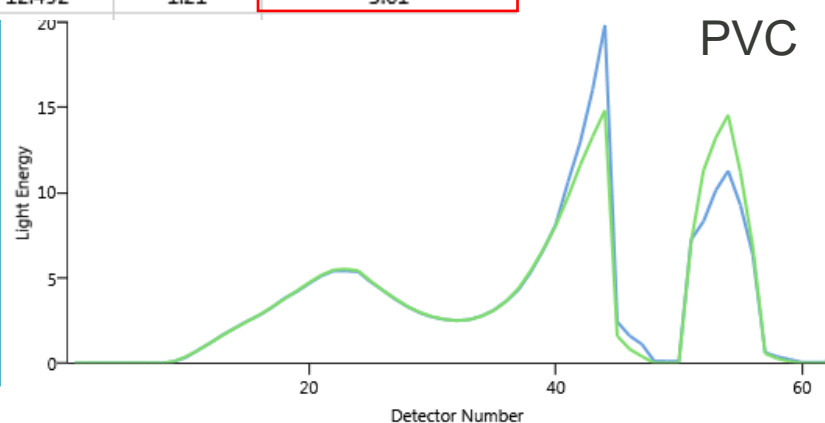


| Sample Name | Dx (10) (µm) | Dx (50) (µm) | Dx (90) (µm) | Residual (%) | Weighted Residual (%) |
|---------------------|--------------|--------------|--------------|--------------|-----------------------|
| plastisol 1.76/0.1 | 0.624 | 7.705 | 27.053 | 0.70 | 0.26 |
| plastisol 1.59/0.1 | 0.655 | 4.606 | 23.874 | 0.86 | 1.26 |
| plastisol 2.51/0.01 | 1.069 | 8.740 | 28.163 | 0.72 | 0.26 |
| plastisol 1.54/0.01 | 0.261 | 0.713 | 12.492 | 1.21 | 3.61 |

2.51/0.01
Pigment

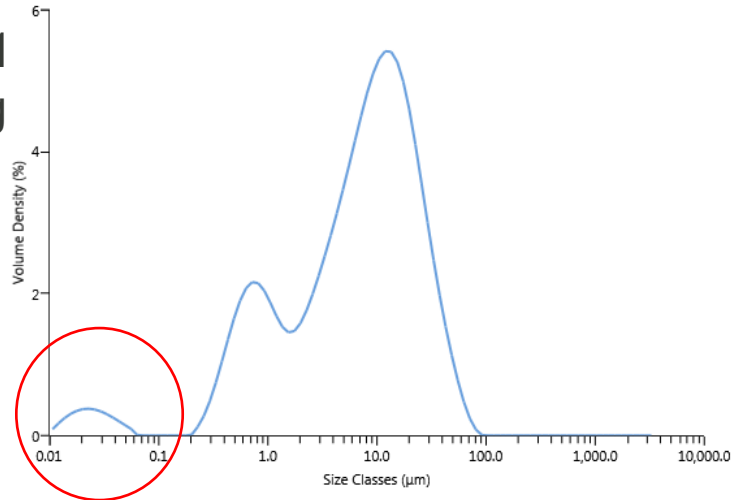


1.54/0.01
PVC

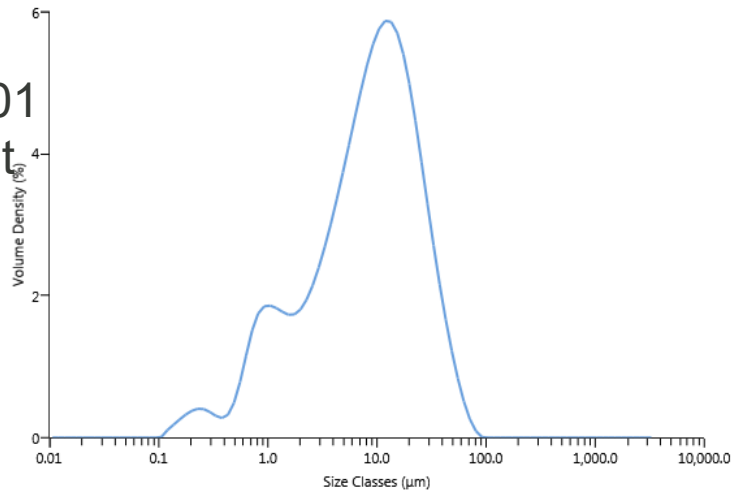


Plastisol – which optical properties to chose?

1.76/0.1
Blowing agent



2.51/0.01
Pigment

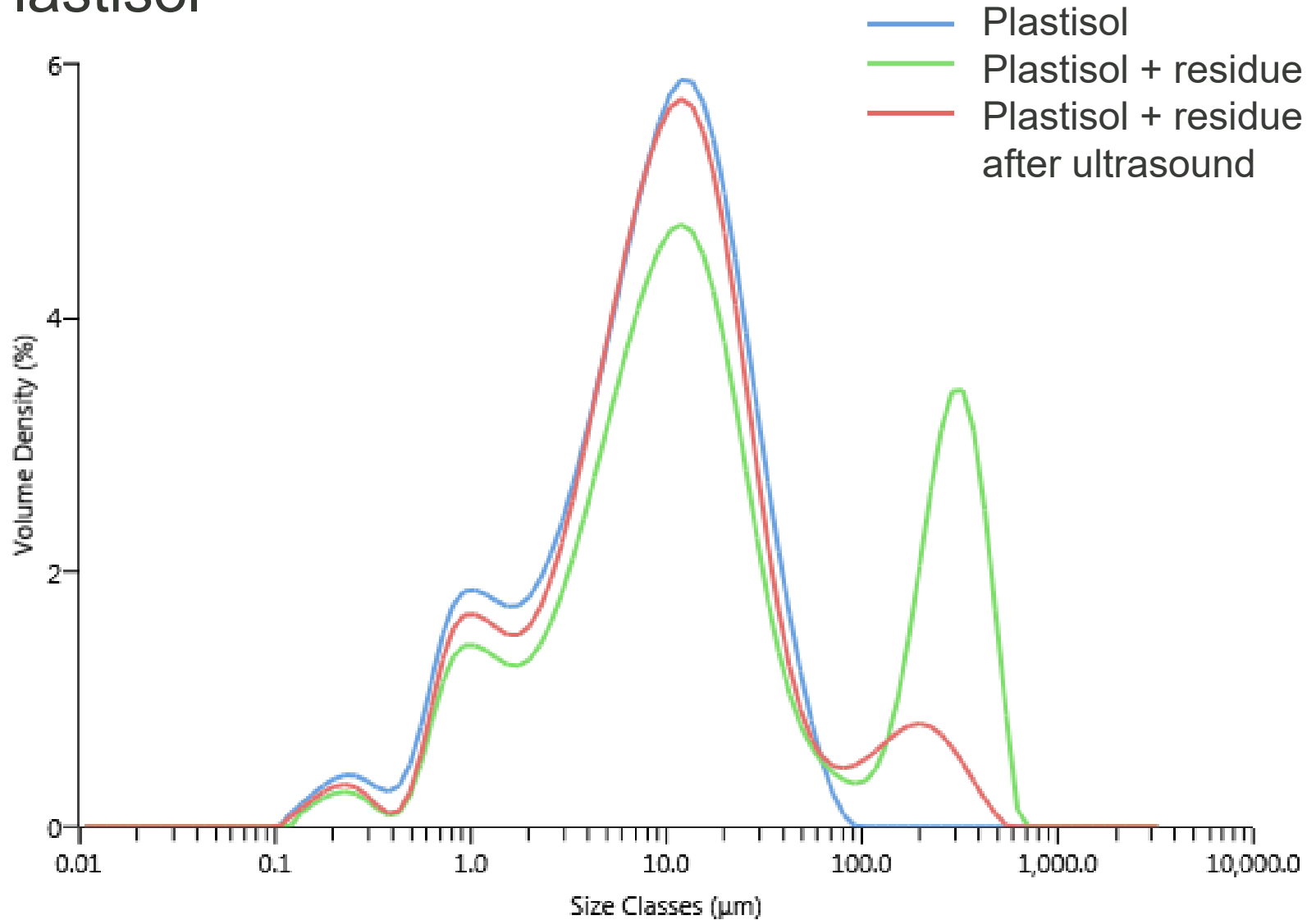


- Similar residuals
- Unresolved peak in 1.76/0.1
- Sample knowledge:
 - Blowing agent >10μm
 - Pigment <1μm
- Laser diffraction knowledge:
 - Sensitivity to optical properties increases as size decreases
- Select 2.51/0.01

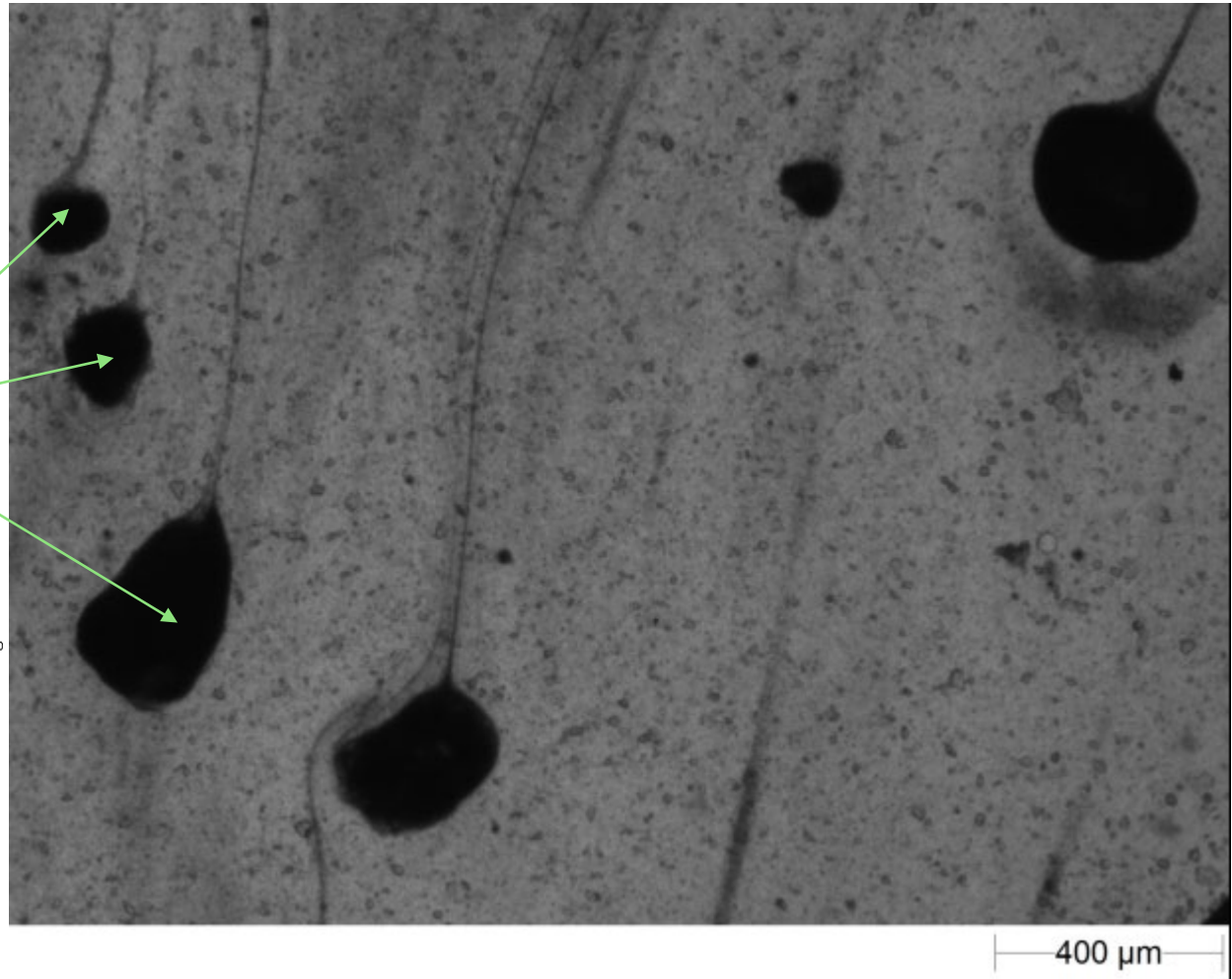
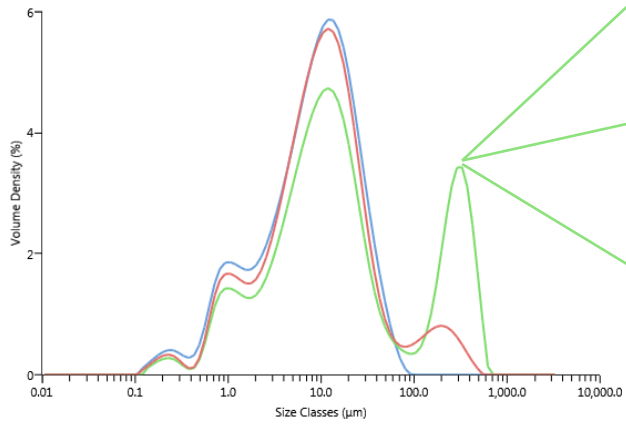


Malvern
Panalytical

Plastisol

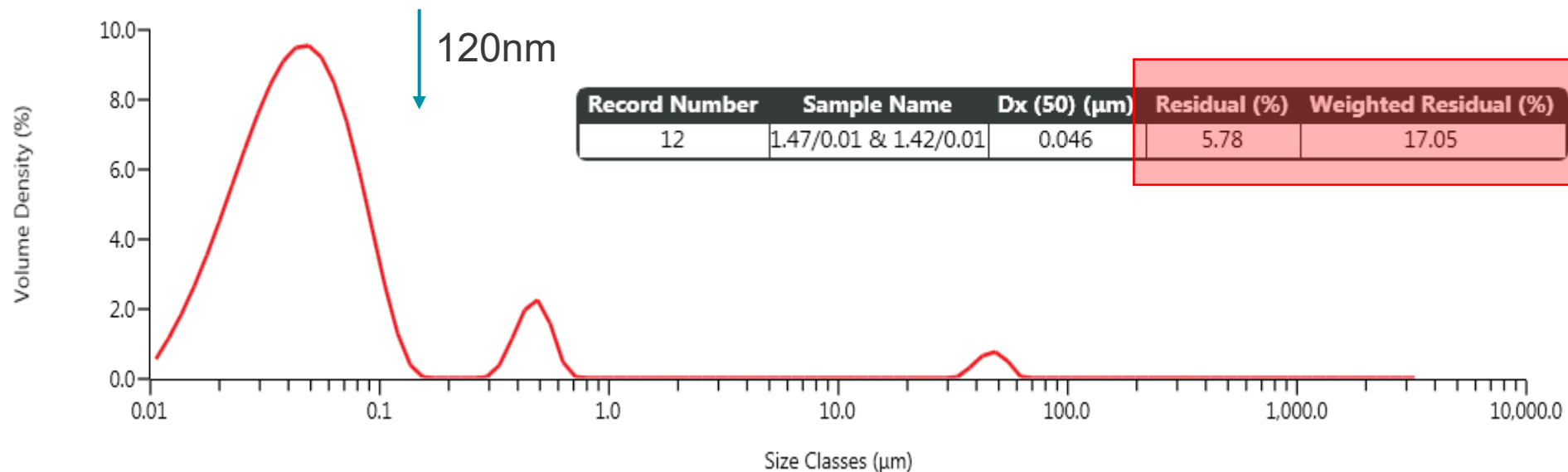


Plastisol



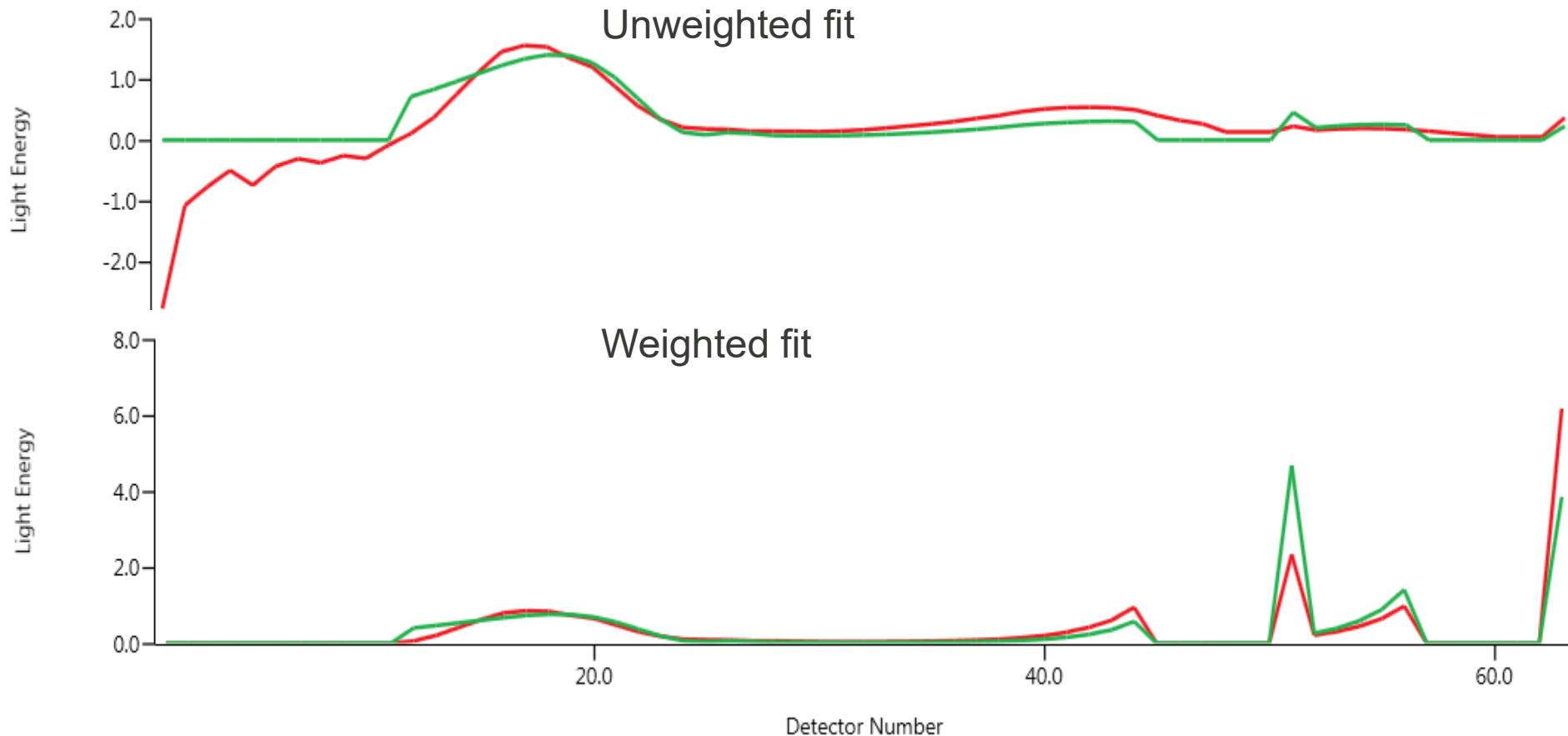
Pigment – a real life example of using the OPO

- Red ink
- Expected Dv50 120nm
- Red 1.47/0.01 & Blue 1.42/0.01 (reference)



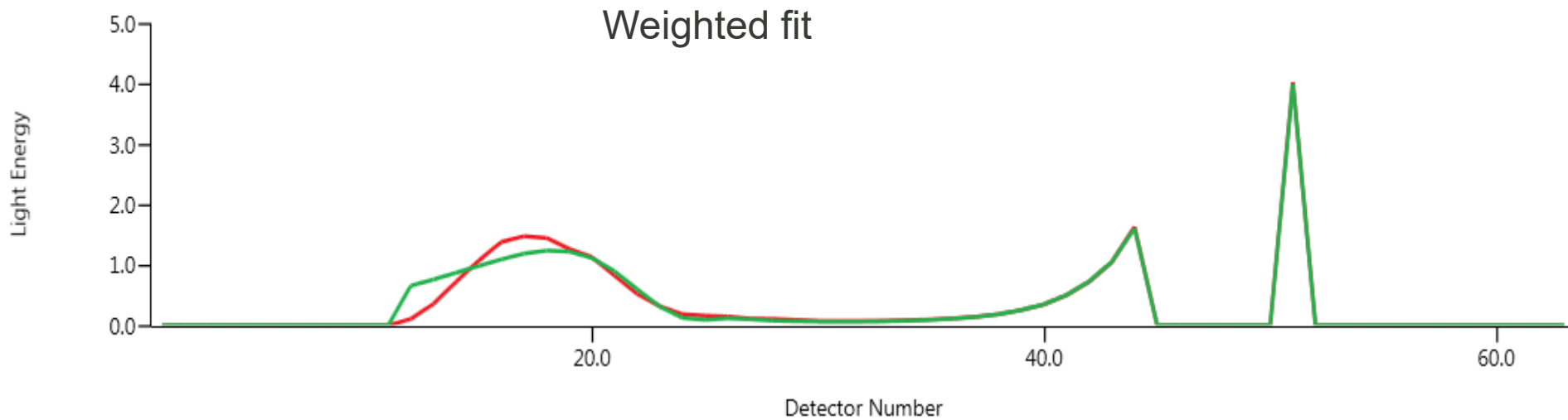
Looking at the fit report for the red ink

- Check fit report



Red ink

- Check fit – **red** light only



| Record Number | Sample Name | Dx (50) (μm) | Residual (%) | Weighted Residual (%) |
|---------------|-------------|---------------------------|--------------|-----------------------|
| 3 | Red ink | 0.179 | 5.80 | 4.09 |

OPO – red ink

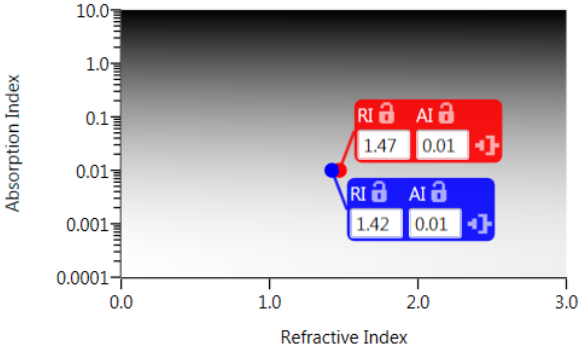
Optical Property Optimizer

Records

| Result graph | Data graph | Record Number | Sample Name |
|-------------------------------------|-------------------------------------|---------------|-----------------------|
| <input checked="" type="checkbox"/> | <input checked="" type="checkbox"/> | 12 | 1.47/0.01 & 1.42/0.01 |

Optical and Analysis Settings

Different blue-light properties Non-Spherical Narrow Modes



Absorption Index

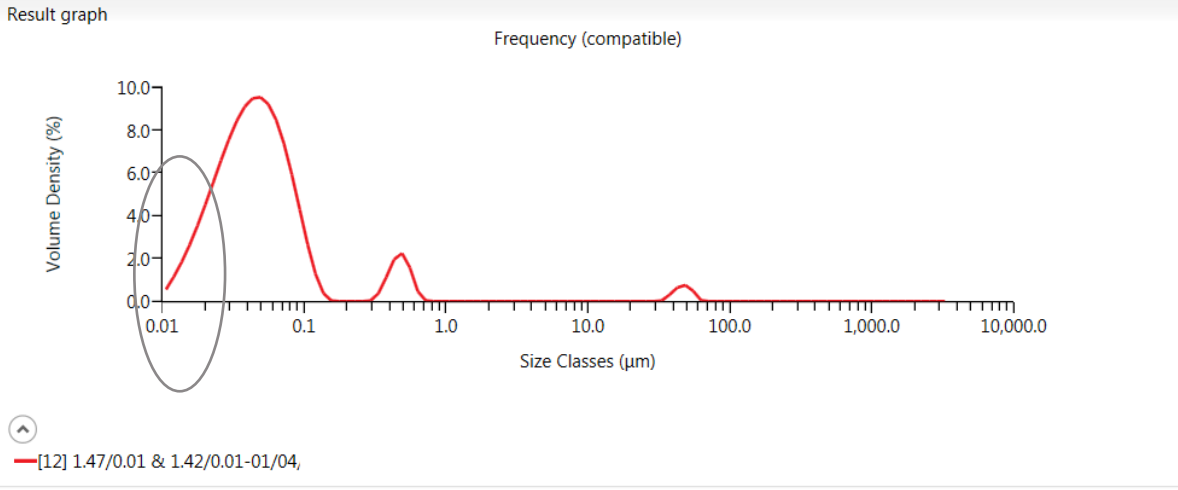
Refractive Index

Optical Property Sets

| Use | Set name | Analysis model | Particle type | Refractive index | Ab: |
|--------------------------|-------------------|----------------|---------------|------------------|------|
| <input type="checkbox"/> | Original Settings | Narrow Modes | NonSpherical | 1.47 | 0.01 |

Result graph

Frequency (compatible)



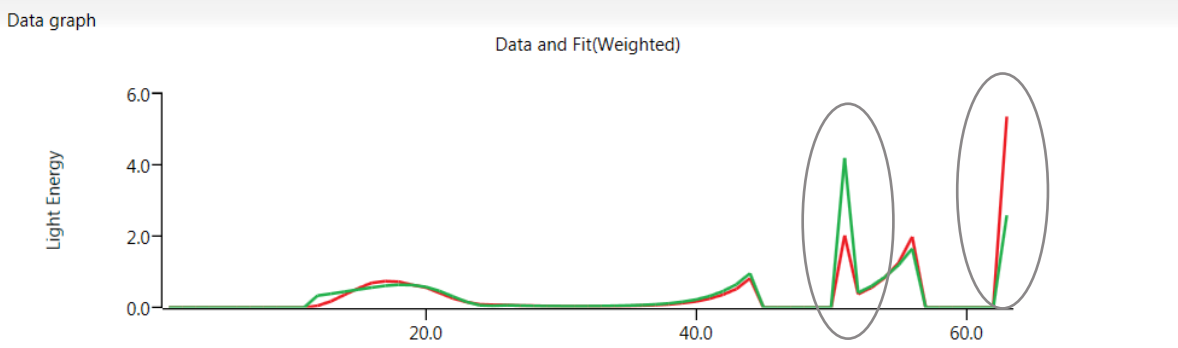
Volume Density (%)

Size Classes (µm)

— [12] 1.47/0.01 & 1.42/0.01-01/04,

Data graph

Data and Fit(Weighted)



Light Energy

| Record Number | Sample Name | Dx (50) (µm) | Residual (%) | Weighted Residual (%) |
|---------------|-----------------------|--------------|--------------|-----------------------|
| 12 | 1.47/0.01 & 1.42/0.01 | 0.046 | 5.78 | 17.05 |

OPO – red ink

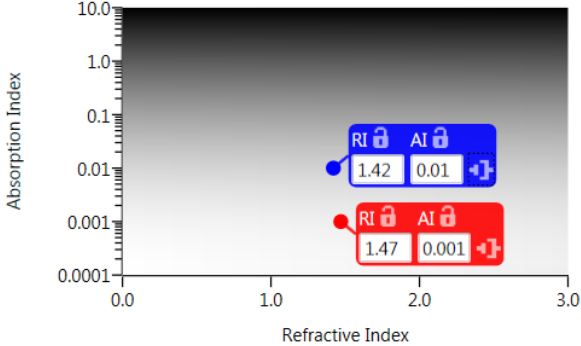
Optical Property Optimizer

Records

| Result graph | Data graph | Record Number | Sample Name |
|-------------------------------------|-------------------------------------|---------------|-----------------------|
| <input checked="" type="checkbox"/> | <input checked="" type="checkbox"/> | 12 | 1.47/0.01 & 1.42/0.01 |

Optical and Analysis Settings

Different blue-light properties Non-Spherical Narrow Modes



Absorption Index vs Refractive Index plot. The y-axis is logarithmic from 0.0001 to 10.0. The x-axis is linear from 0.0 to 3.0. Two data points are shown: a blue point at RI=1.42, AI=0.01 and a red point at RI=1.47, AI=0.001.

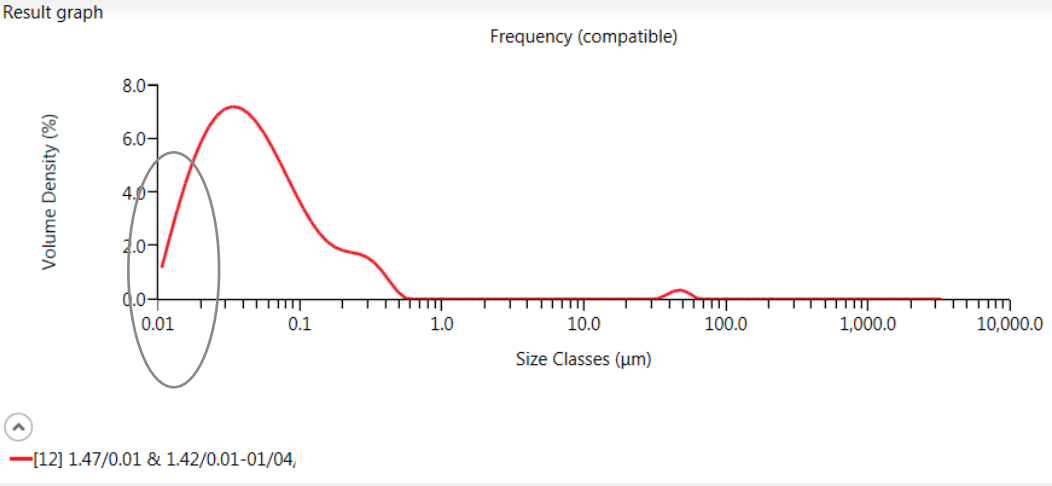
Optical Property Sets

Store Remove Remove All

| Use | Set name | Analysis model | Particle type | Refractive index | Ab: |
|-------------------------------------|-------------------|----------------|---------------|------------------|------|
| <input checked="" type="checkbox"/> | Original Settings | Narrow Modes | NonSpherical | 1.47 | 0.01 |

Result graph

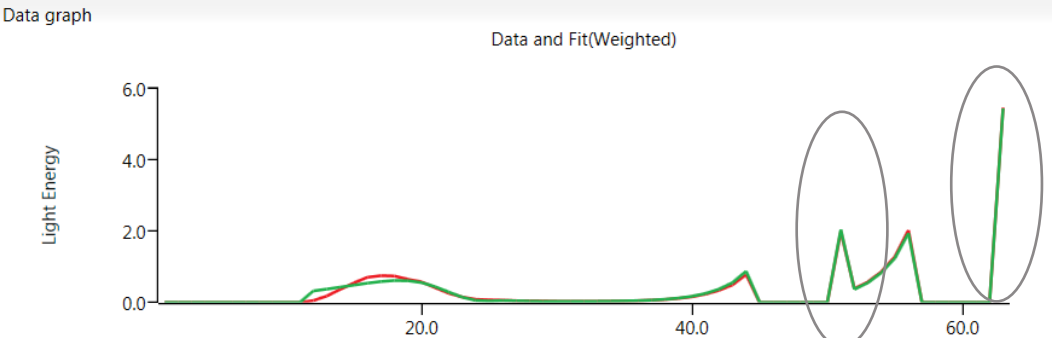
Frequency (compatible)



Volume Density (%) vs Size Classes (µm) plot. The y-axis is linear from 0.0 to 8.0. The x-axis is logarithmic from 0.01 to 10,000.0. A red curve shows a peak at approximately 0.05 µm. A grey oval highlights the peak region.

Data graph

Data and Fit(Weighted)



Light Energy vs Wavelength plot. The y-axis is linear from 0.0 to 6.0. The x-axis is linear from 0.0 to 60.0. A green curve shows two peaks at approximately 500 nm and 650 nm. Grey ovals highlight these peaks.

| Record Number | Sample Name | Dx (50) (µm) | Residual (%) | Weighted Residual (%) |
|---------------|------------------------|--------------|--------------|-----------------------|
| 13 | 1.47/0.001 & 1.42/0.01 | 0.043 | 5.50 | 2.26 |

OPO – red ink

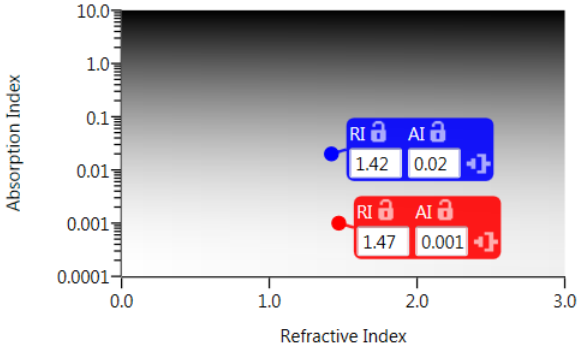
Optical Property Optimizer

Records

| Result graph | Data graph | Record Number | Sample Name |
|-------------------------------------|-------------------------------------|---------------|------------------------|
| <input checked="" type="checkbox"/> | <input checked="" type="checkbox"/> | 13 | 1.47/0.001 & 1.42/0.01 |

Optical and Analysis Settings

Different blue-light properties Non-Spherical Narrow Modes



Absorption Index

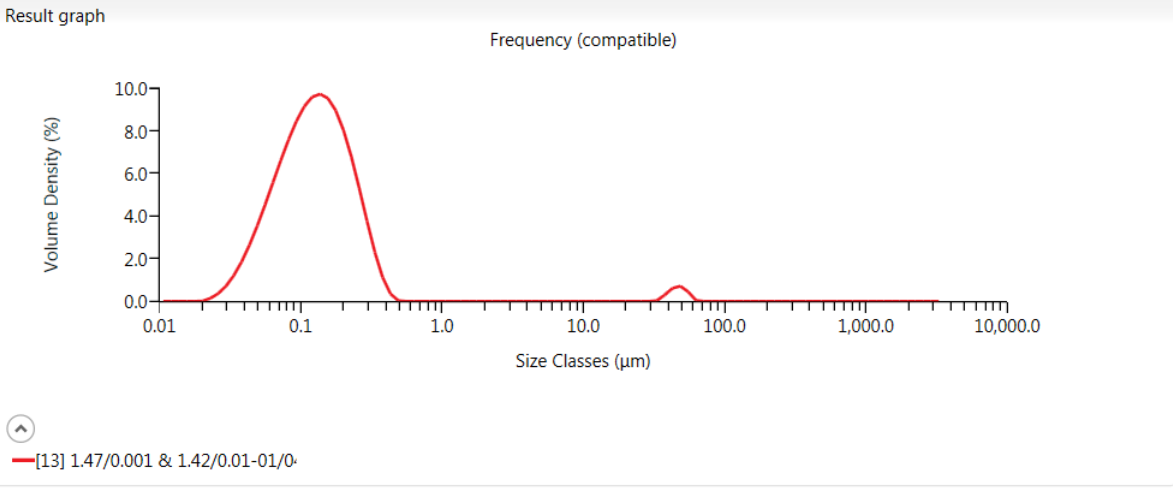
Refractive Index

Optical Property Sets

| Use | Set name | Analysis model | Particle type | Refractive index | Abs |
|-------------------------------------|-------------------|----------------|---------------|------------------|------|
| <input type="checkbox"/> | intermediate | Narrow Modes | NonSpherical | 1.47 | 0.00 |
| <input type="checkbox"/> | Original Settings | Narrow Modes | NonSpherical | 1.47 | 0.00 |
| <input checked="" type="checkbox"/> | final | Narrow Modes | NonSpherical | 1.47 | 0.00 |

Result graph

Frequency (compatible)



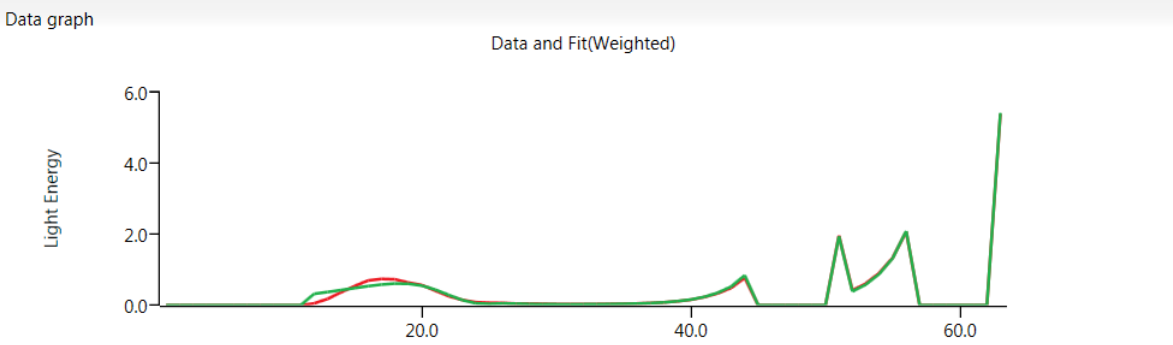
Volume Density (%)

Size Classes (µm)

[13] 1.47/0.001 & 1.42/0.01-01/0-

Data graph

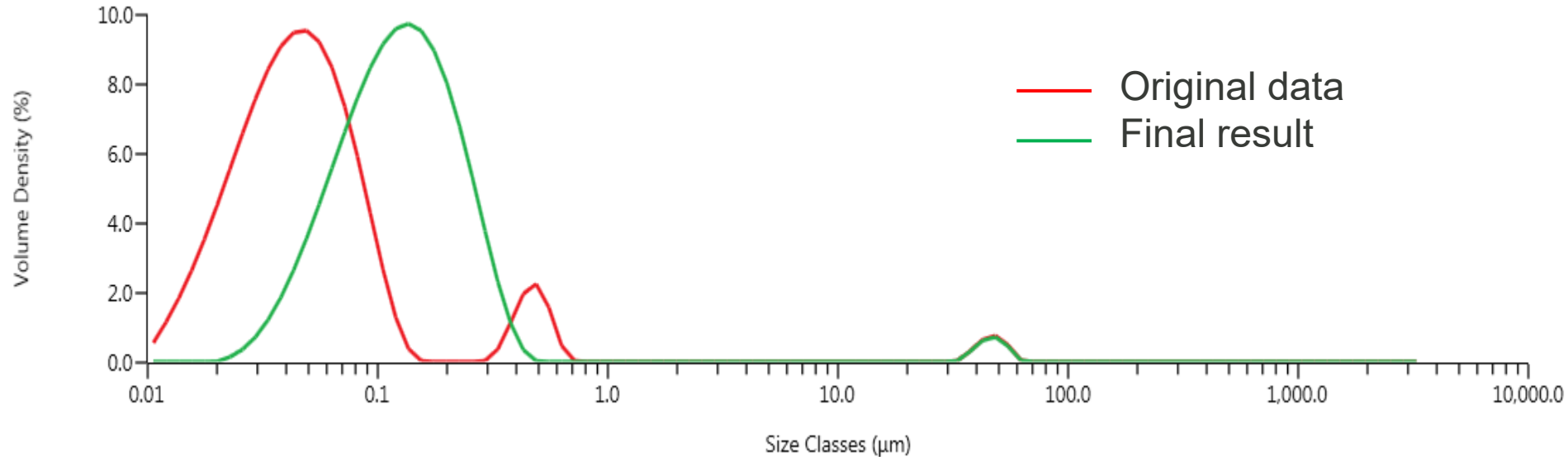
Data and Fit(Weighted)



Light Energy

| Record Number | Sample Name | Dx (50) (µm) | Residual (%) | Weighted Residual (%) |
|---------------|------------------------|--------------|--------------|-----------------------|
| 15 | 1.47/0.001 & 1.42/0.02 | 0.125 | 5.50 | 2.14 |

Corrected red ink result



| Record Number | Sample Name | Dx (50) (µm) | Residual (%) | Weighted Residual (%) |
|---------------|------------------------|--------------|--------------|-----------------------|
| 12 | 1.47/0.01 & 1.42/0.01 | 0.046 | 5.78 | 17.05 |
| 15 | 1.47/0.001 & 1.42/0.02 | 0.125 | 5.50 | 2.14 |

- Dv50 now 125nm
- Close to expected value of 120nm



General Purpose

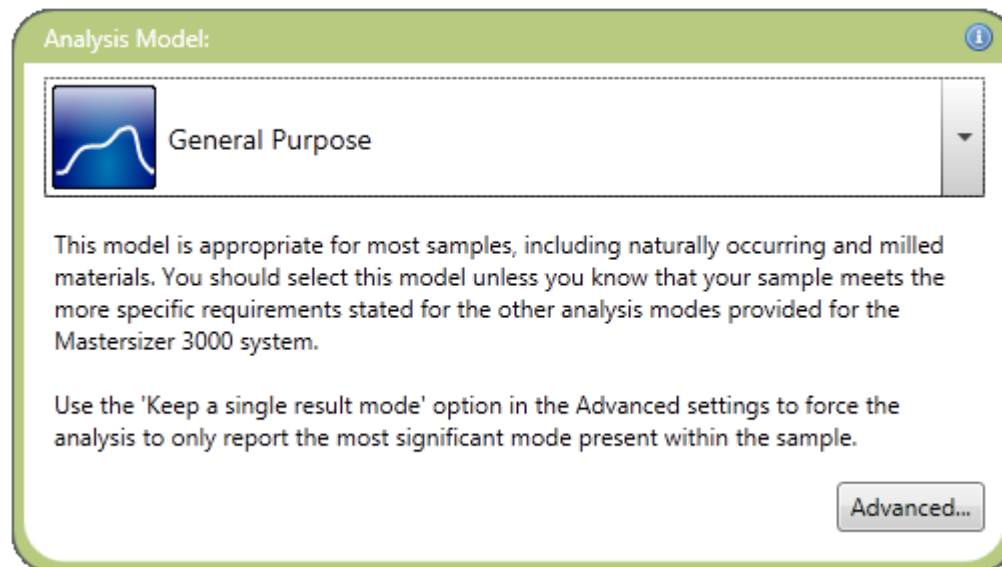
This model is appropriate for most samples, including solids. You should select this model unless you know of specific requirements stated for the other analysis models on the Versizer 3000 system.

Select the 'Keep a single result mode' option in the Advanced Analysis to only report the most significant mode present.

Analysis models

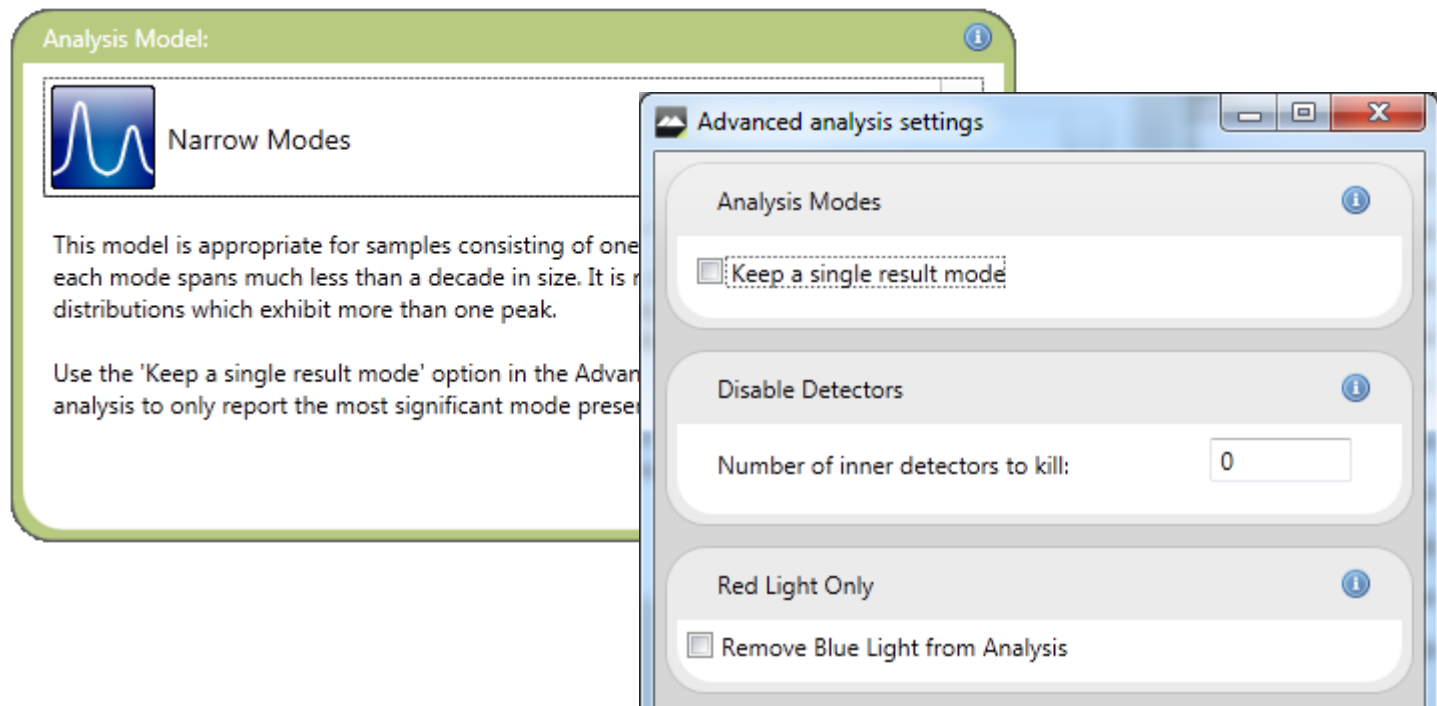
Analysis Models

- Using a particular analysis model enables the software to better interpret the light scattering data.
- General purpose
 - Suitable for the majority of samples



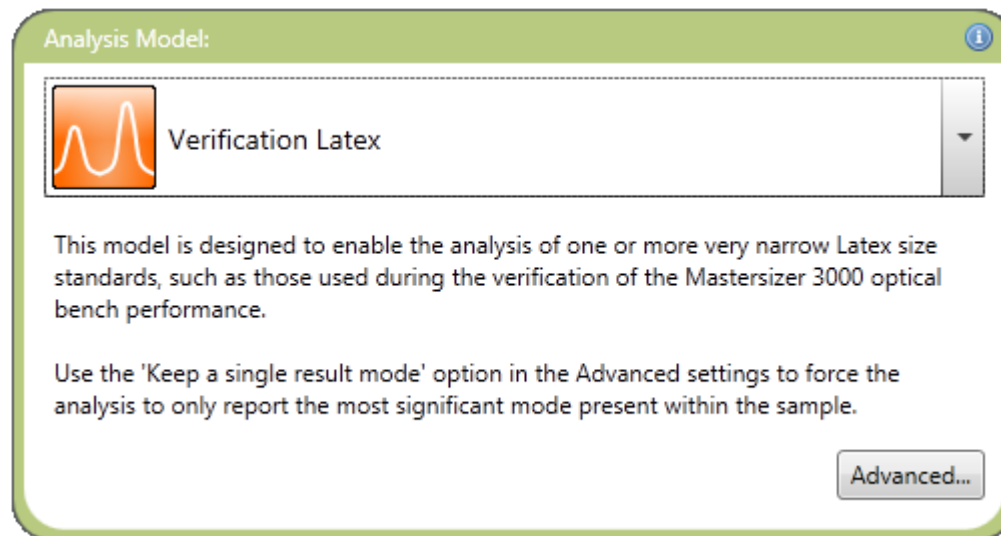
Analysis Models

- Using a particular analysis model enables the software to better interpret the light scattering data.
- Narrow modes
 - Suitable for samples with modes spanning much less than a decade in size



Analysis Models


- Using a particular analysis model enables the software to better interpret the light scattering data.
- Verification latex
 - Suitable for use with very narrow latex size standards



Particle shape, irregular or spherical?

- Irregularly shaped small particles ($<1\mu\text{m}$) depolarise light more strongly in one direction
- The Non-Spherical option enables this high angle scattering to be correctly interpreted.
- Non-Spherical is the default option as the majority of materials are irregular

Particle Type ?

 Non-Spherical


This particle type is applicable for particles which are irregular in shape, or have a rough surface structure. For example, it should be selected for milled or crushed materials.

This type uses Mie Theory, and therefore requires input of the optical properties of your sample in order to calculate a particle size distribution. The advantage of this is that it provides the possibility of obtaining accurate size distributions for all particle sizes.

Particle shape, irregular or spherical?

- Spherical samples are rarer but will include glass beads, latex spheres and emulsions.

Particle Type ⓘ

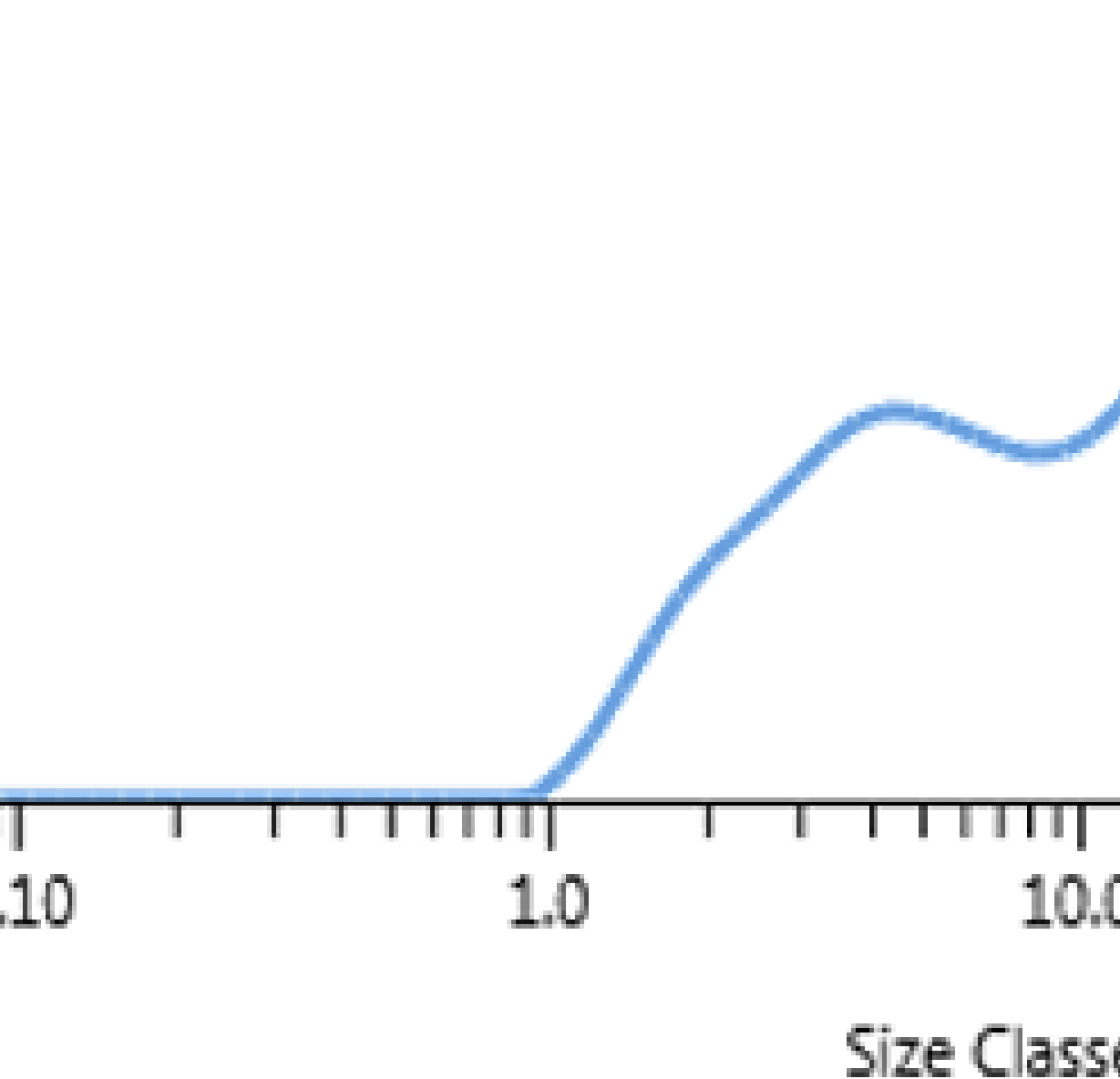
 Spherical

This particle type is applicable for particles which are perfectly spherical in shape. For example, it should be selected for polymer latex samples or for emulsions.

This type uses Mie Theory, and therefore requires input of the optical properties of your sample in order to calculate a particle size distribution. The advantage of this is that it provides the possibility of obtaining accurate size distributions for all particle sizes.

Fine powder mode

- This removes the noise on the inner detectors caused by temperature gradients in dry powder feeder measurements
 - Or from using ultrasound in organic solvents
- It is only used in measurements of samples which are finer than 600 microns



Understanding
the size
distribution



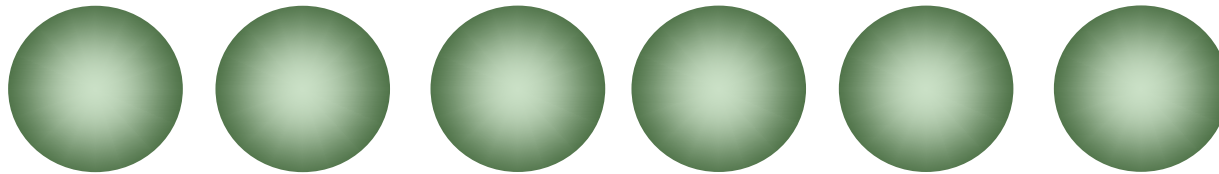
Understanding the size distribution



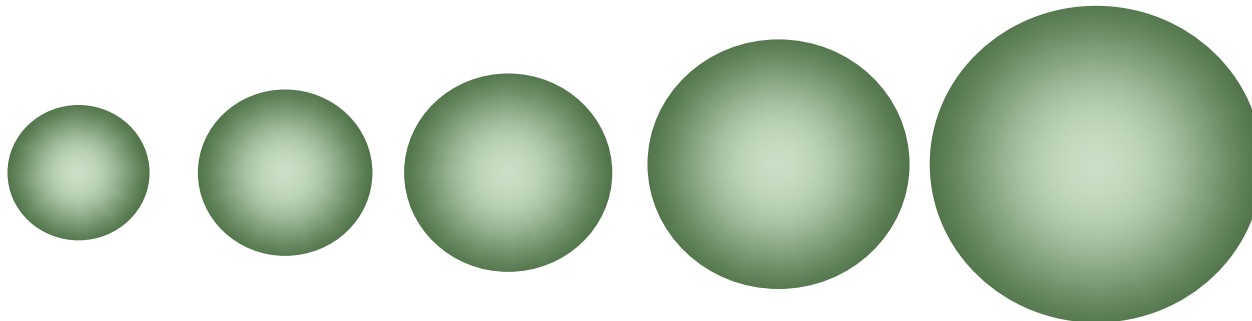
- The Mastersizer 3000 system is designed such that equal volumes of particles of different sizes produce a similar measured scattering intensity.
- The size distribution is reported as a volume distribution as this best reflects the sensitivity of the system.
- What does this mean in practice?

Particle size distributions

- If a sample contains particles that are all the same size it is described as monodisperse

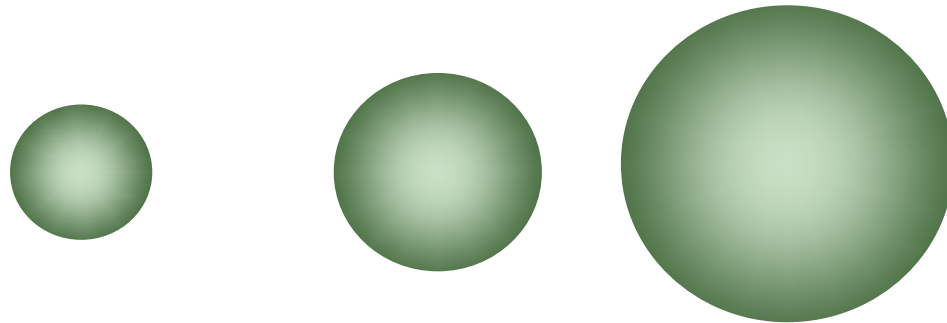


- Most real world samples will contain a distribution of particle sizes and are described as polydisperse



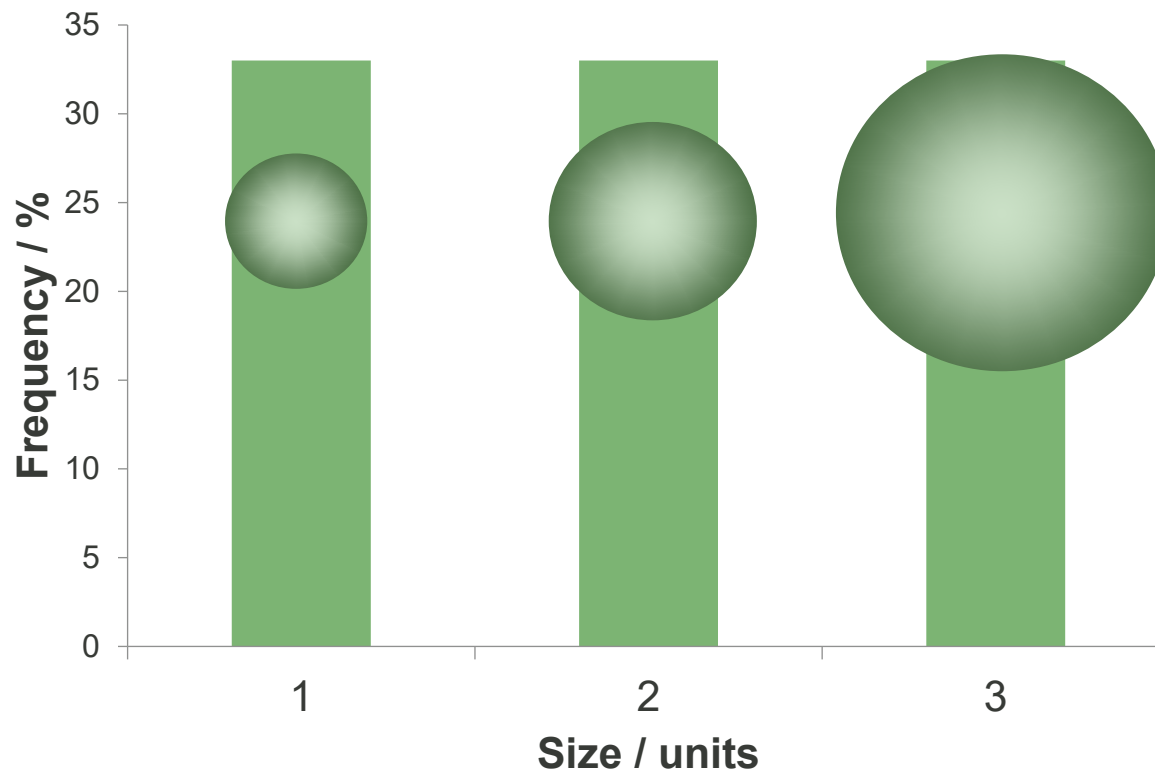
Understanding distribution weighting

- If we have a sample made up of 3 particles
 - With diameters of 1, 2 and 3 units.



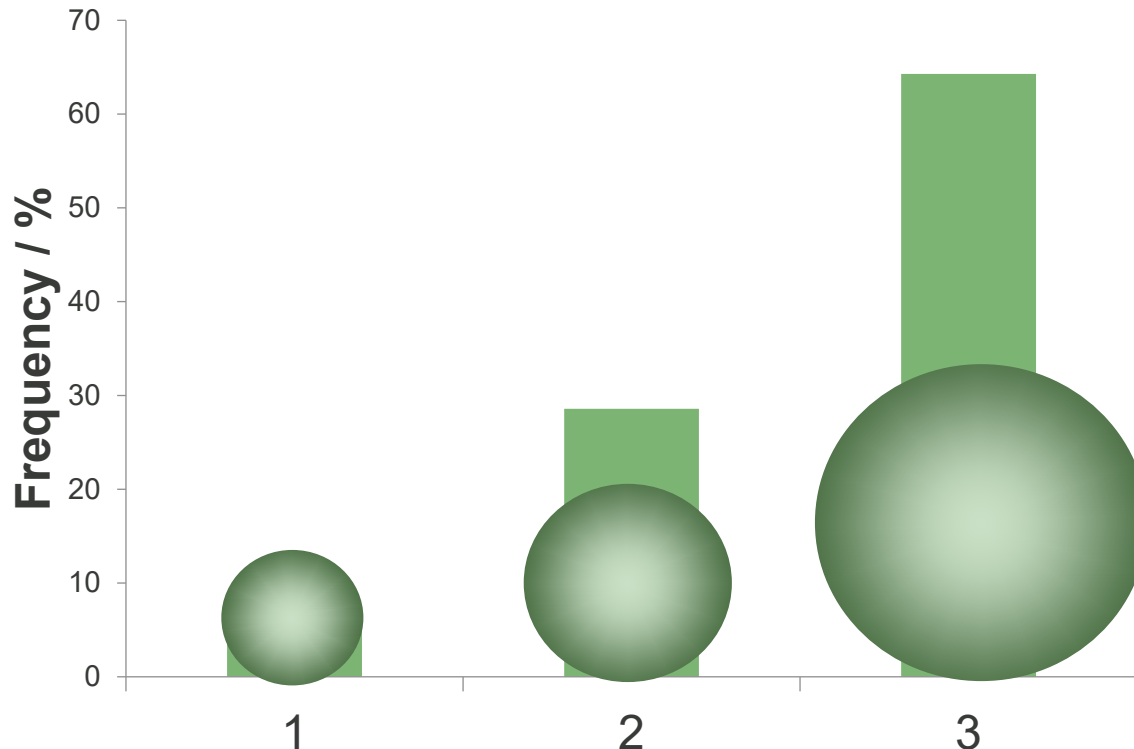
Understanding distribution weighting

- If we measure the sample by a counting technique (such as microscopy) we get a number weighted distribution...
 - And each particle contributes equally

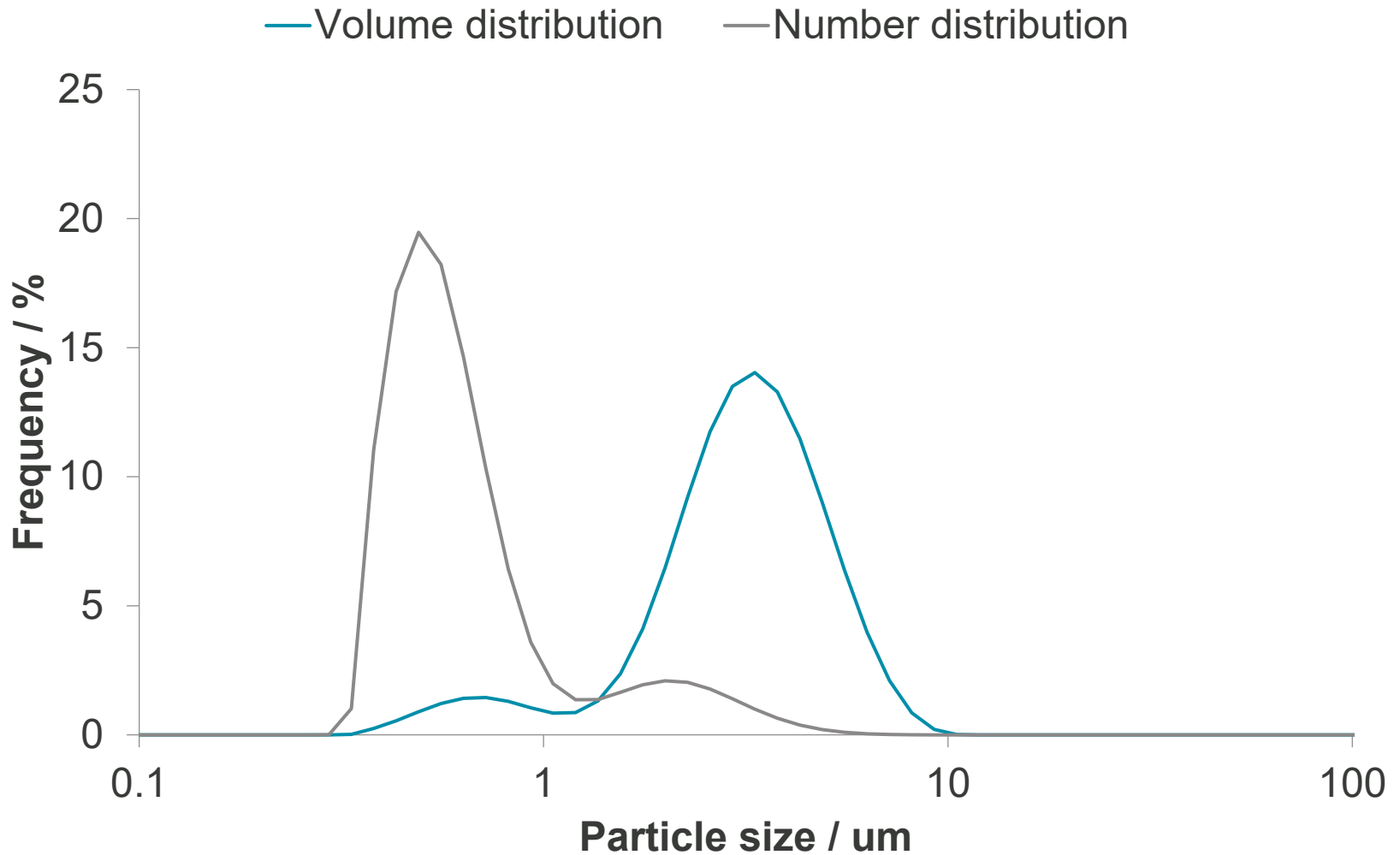


Understanding distribution weighting

- If we measure the sample by laser diffraction we get a distribution weighted by volume
 - And the largest particle features most prominently

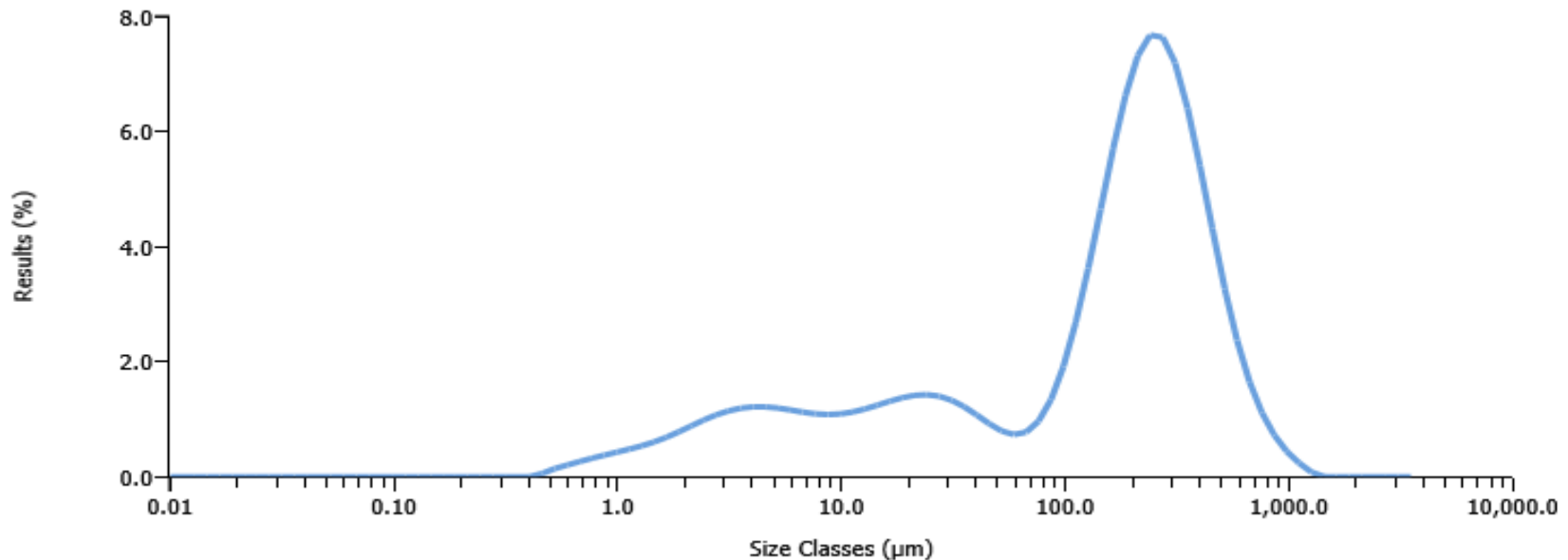


Real example of number and volume distributions



Representing the size distribution

- The particle size distributions are plotted on a logarithmic axis

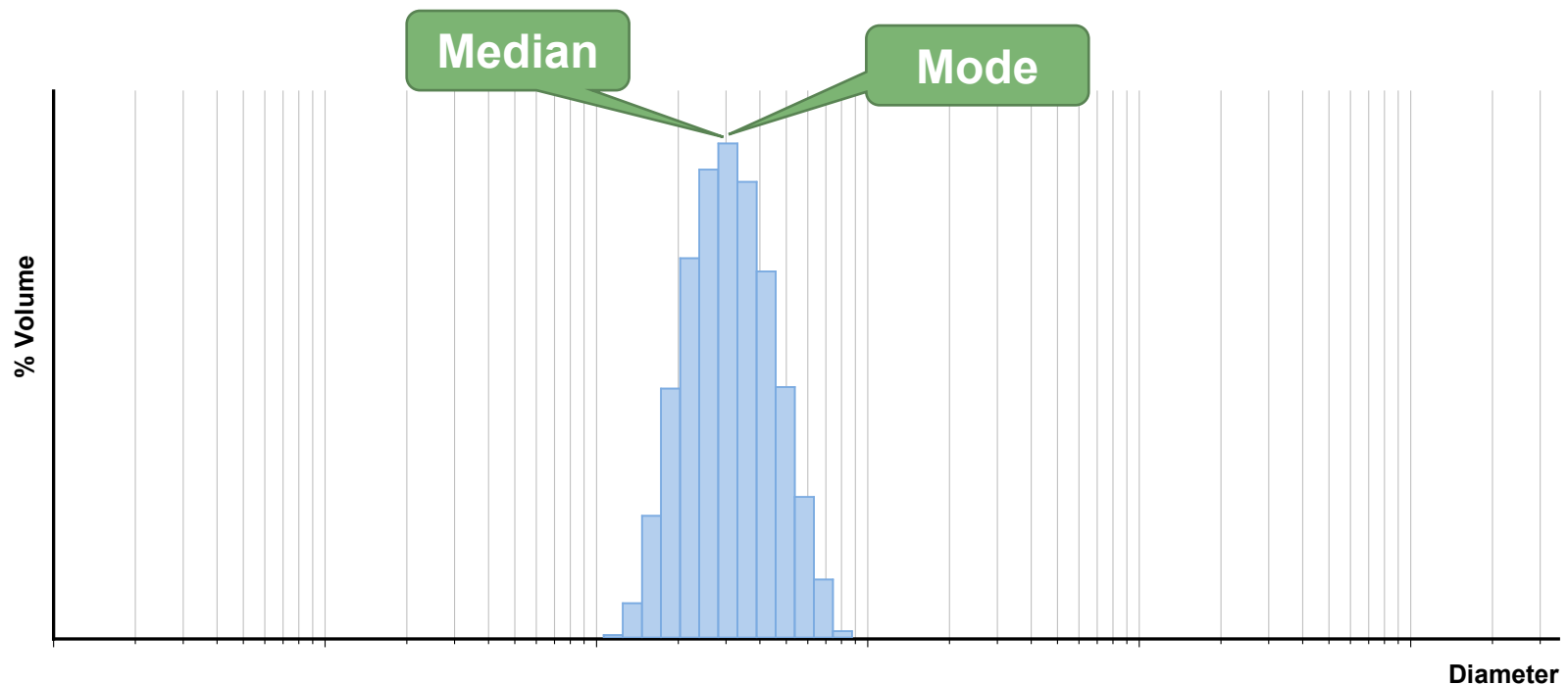


| Result | |
|---|----------------------------|
| Concentration 0.0834 % | Span 0.910 |
| Uniformity 0.277 | Result Units Volume |
| Specific Surface Area 104.6 m ² /kg | Dv (10) 38.670 µm |
| D [3,2] 57.344 µm | Dv (50) 61.177 µm |
| D [4,3] 64.046 µm | Dv (90) 94.343 µm |

Particle size distribution statistics: Median and Mode

- Median = midpoint of the distribution
- Mode = most commonly occurring size class

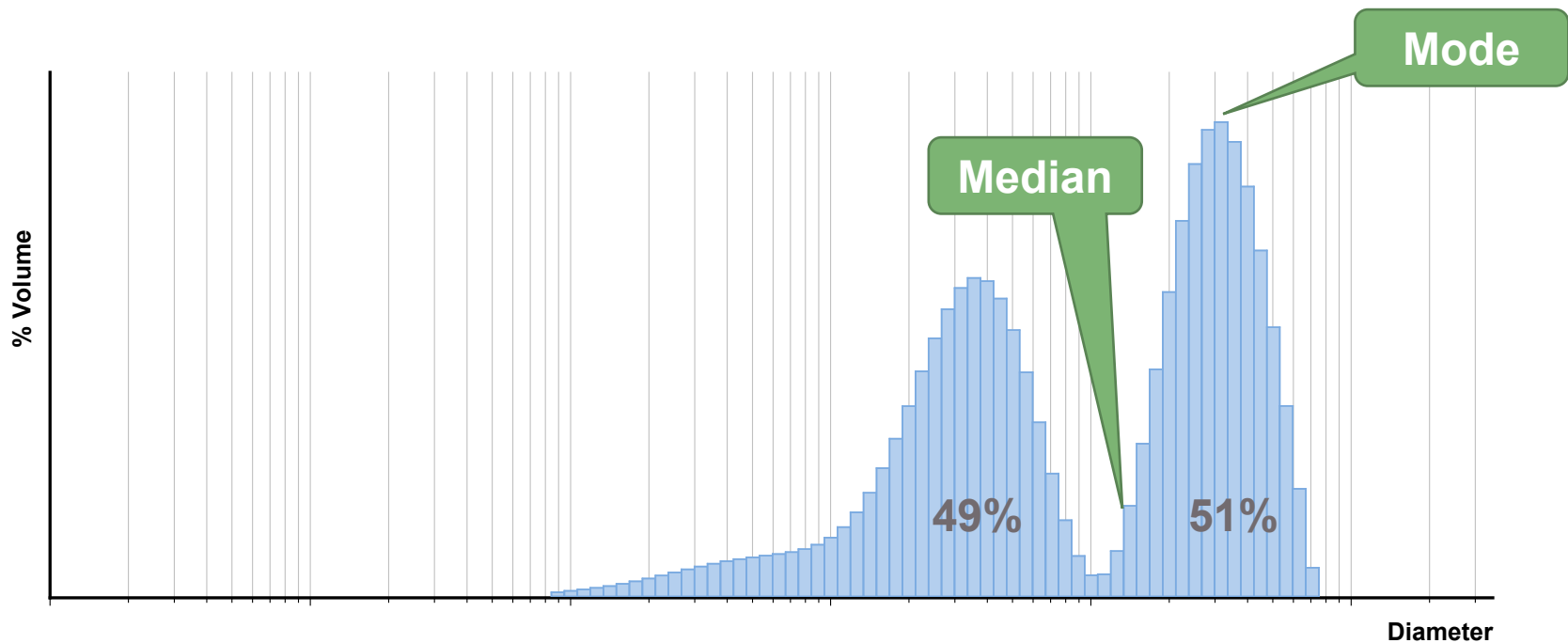
Gaussian Distribution



PSD Statistics: Median and Mode

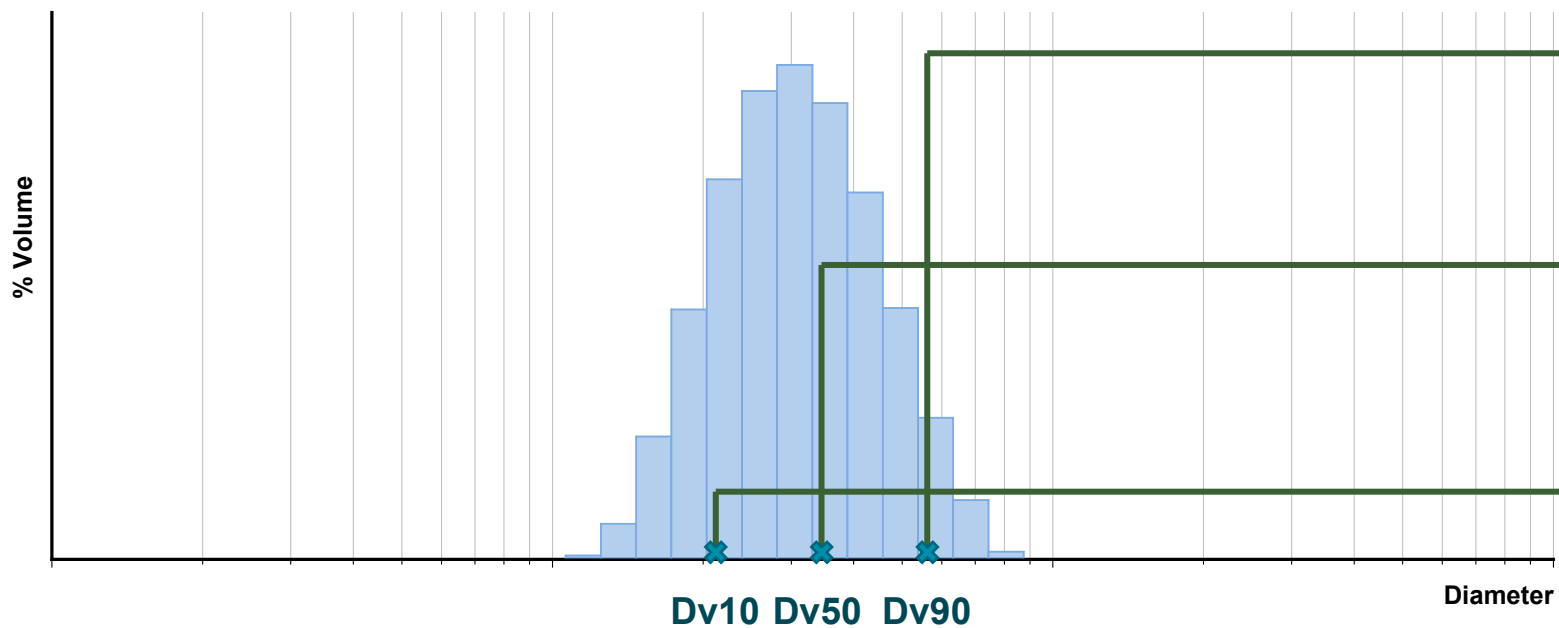
- If the distribution shape is more complex then these parameters will diverge

Bimodal Distribution



PSD Statistics: Percentiles

- Percentiles are the size below which there is a certain volume of the sample
- Taken from the cumulative distribution



| Record Number | Dx (10) (μm) | Dx (50) (μm) | Dx (90) (μm) |
|---------------|---------------------------|---------------------------|---------------------------|
| 43 | 2.08 | 3.32 | 5.43 |

PSD Statistics: Percentiles

- Percentiles can be referred to in several ways
 - The 50th percentile
 - Dv50
 - D[v,0.5]
 - $x_{0.5}$ (in the ISO standard)
- The 'v' signifies that the percentile is from a volume distribution
 - Dx50 is used when the result may be volume or number
- The Dv10 and Dv90 are also reported by default
- Other percentiles can also be reported

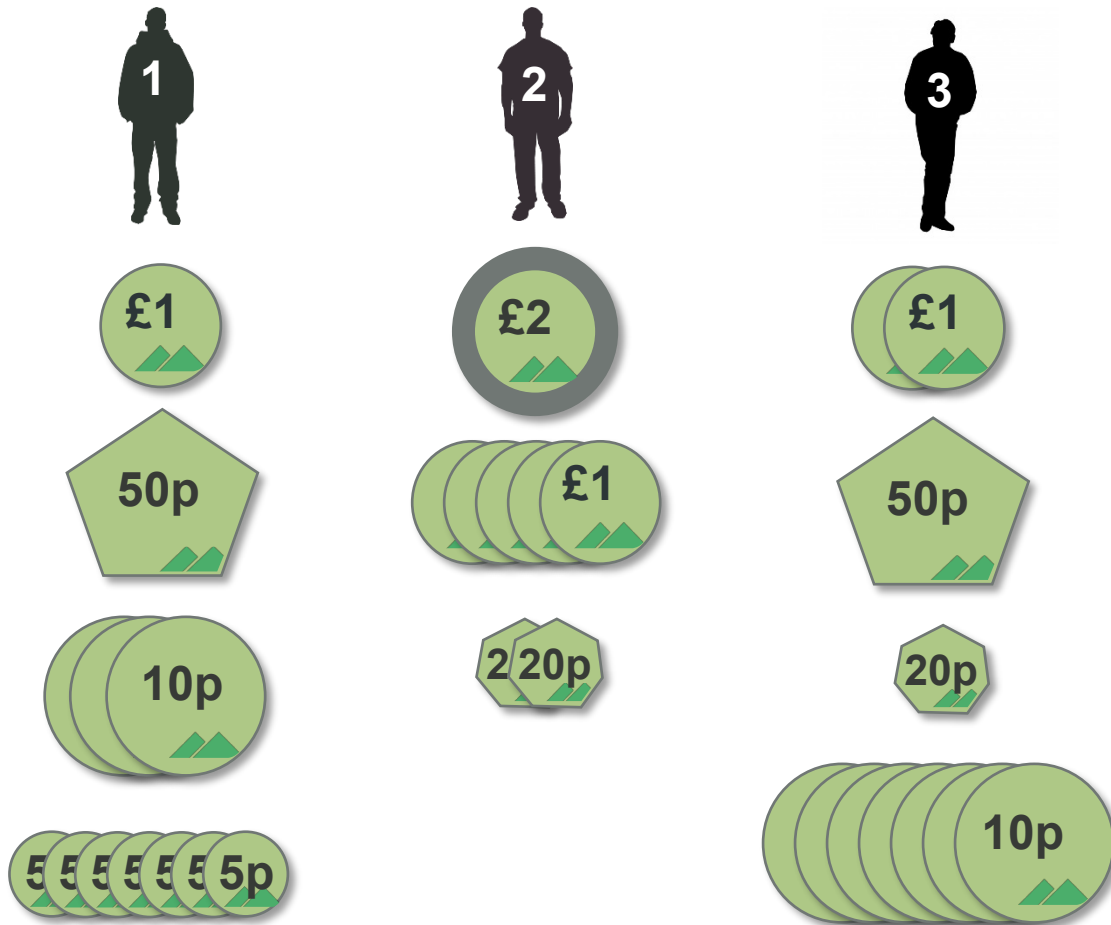
PSD Statistics: Mean particle sizes

- The most familiar mean is the arithmetic mean

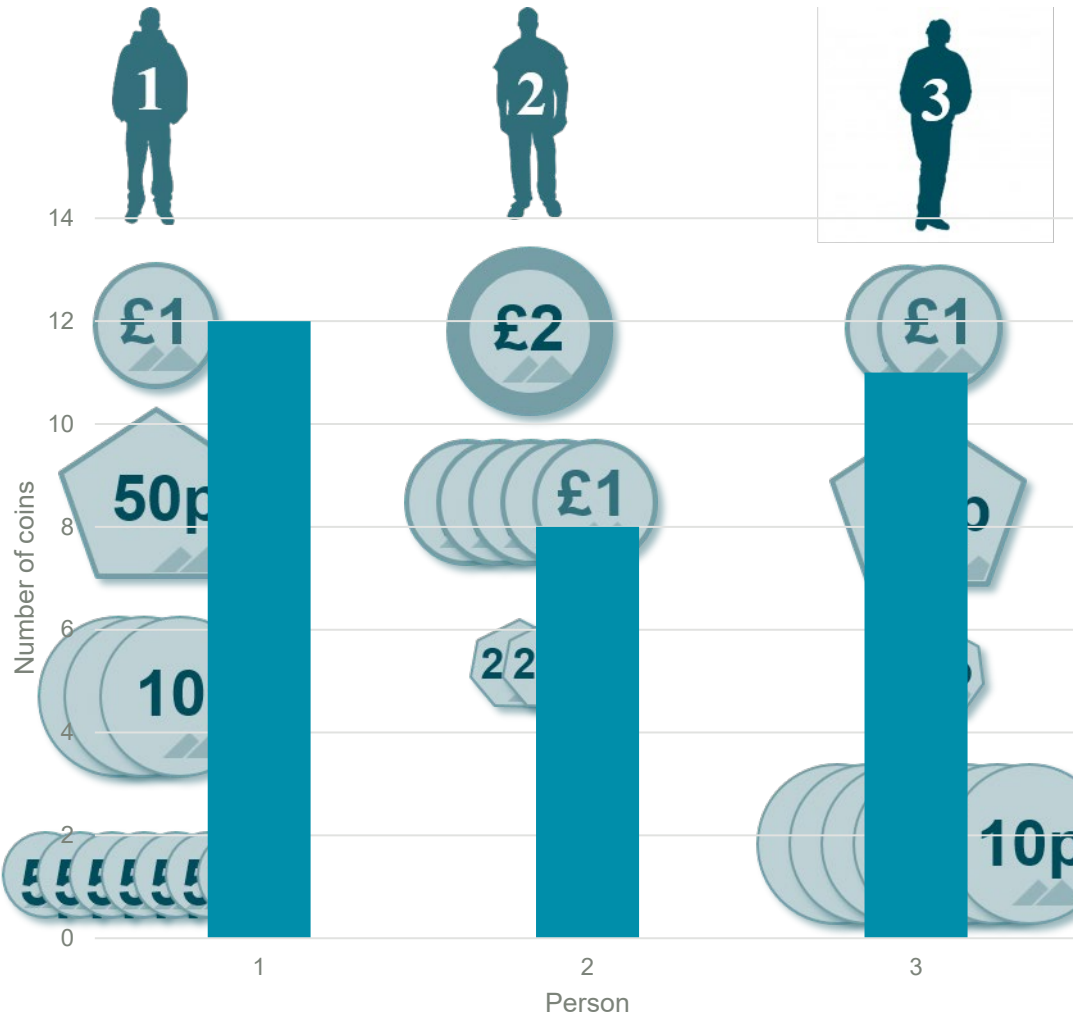
$$X_{nl} = D[1,0] = \frac{1^1 + 2^1 + 3^1}{1^0 + 2^0 + 3^0} = \frac{1 + 2 + 3}{3} = 2$$

- Laser diffraction does not use the arithmetic mean
 - As it measures the volume of particles, not the number
- Laser diffraction reports the volume weighted mean
 - $D[4,3]$
- And surface area weighted mean
 - $D[3,2]$

Different types of average

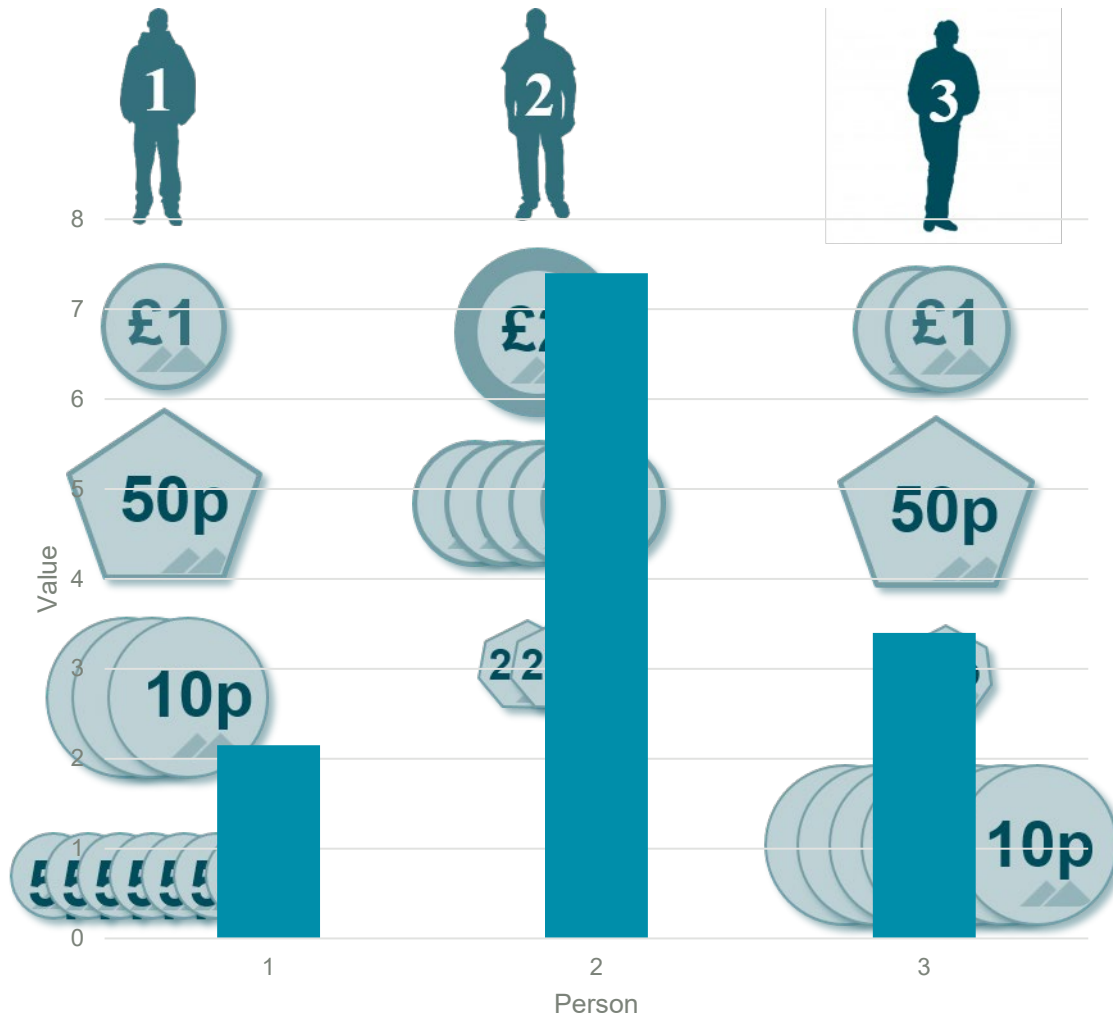


Different types of average



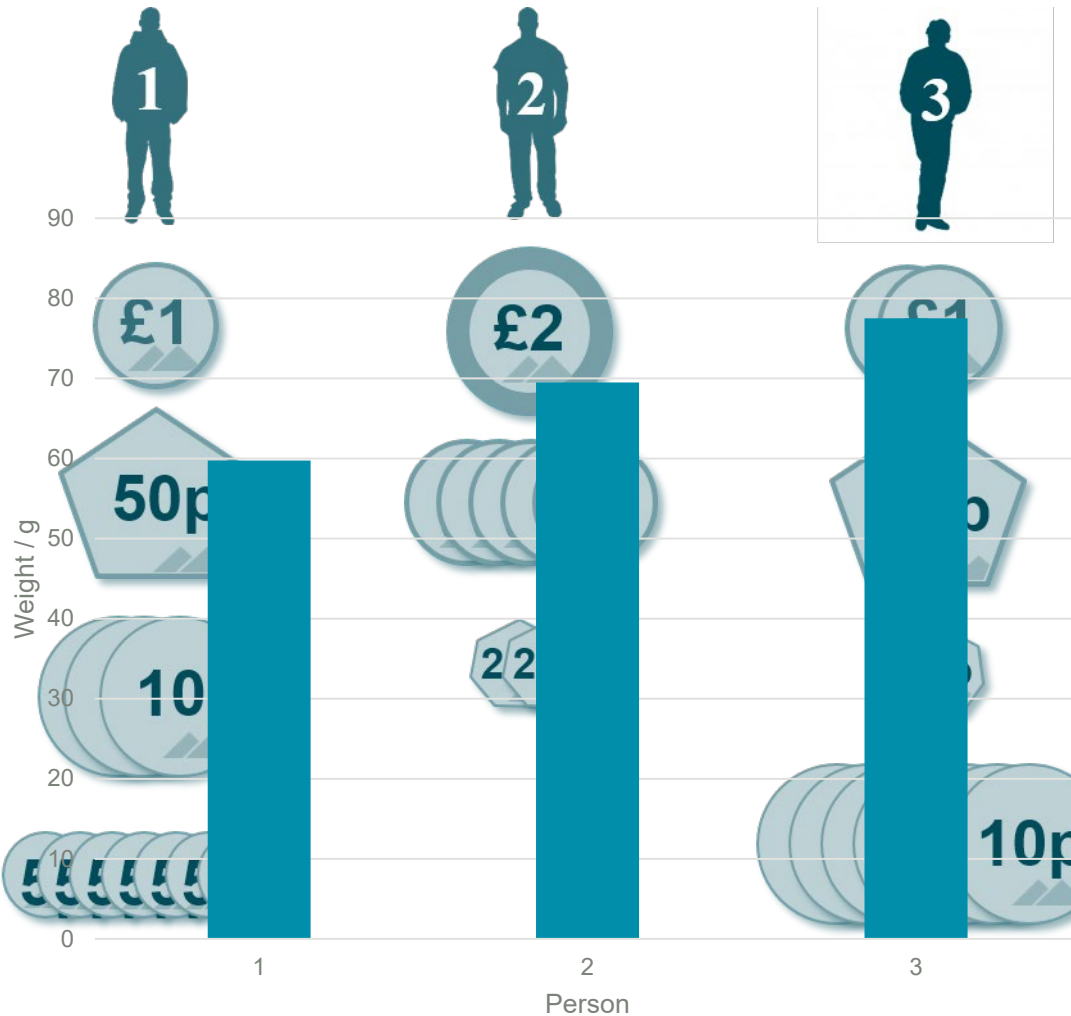
| Type | Average |
|--------|---------|
| Number | 10.33 |

Different types of average



| Type | Average |
|--------|---------|
| Number | 10.33 |
| Value | 4.31 |

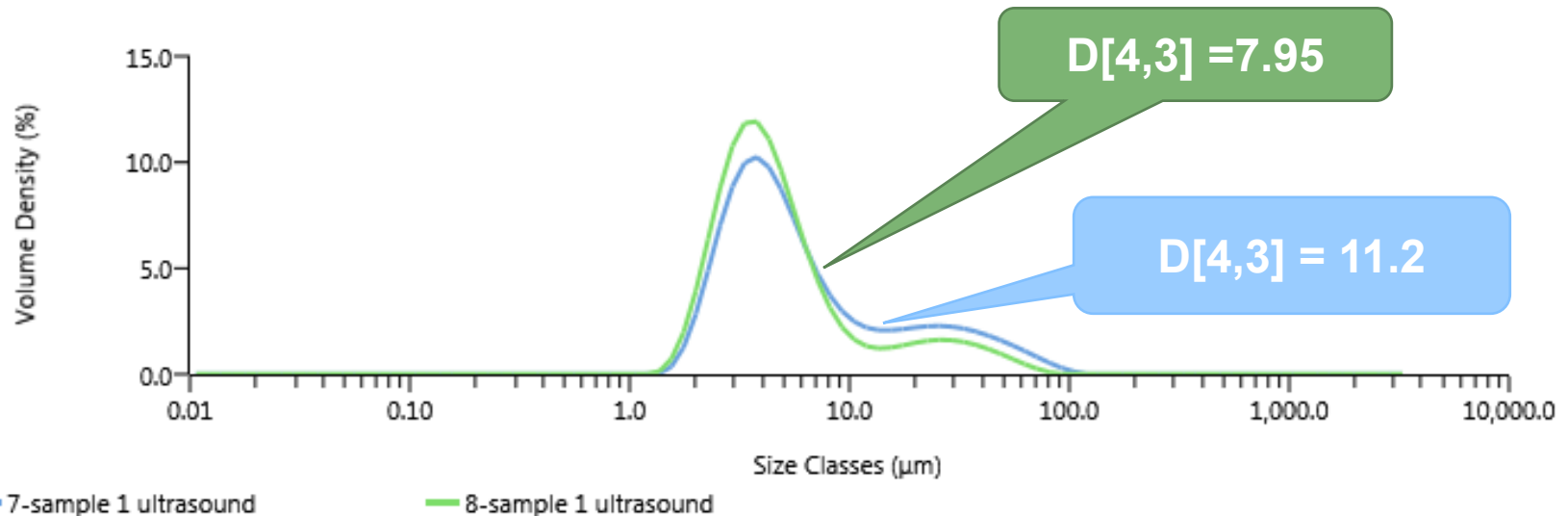
Different types of average



| Type | Average |
|--------|---------|
| Number | 10.33 |
| Value | 4.31 |
| Weight | 68.9 |

PSD Statistics: Volume weighted mean

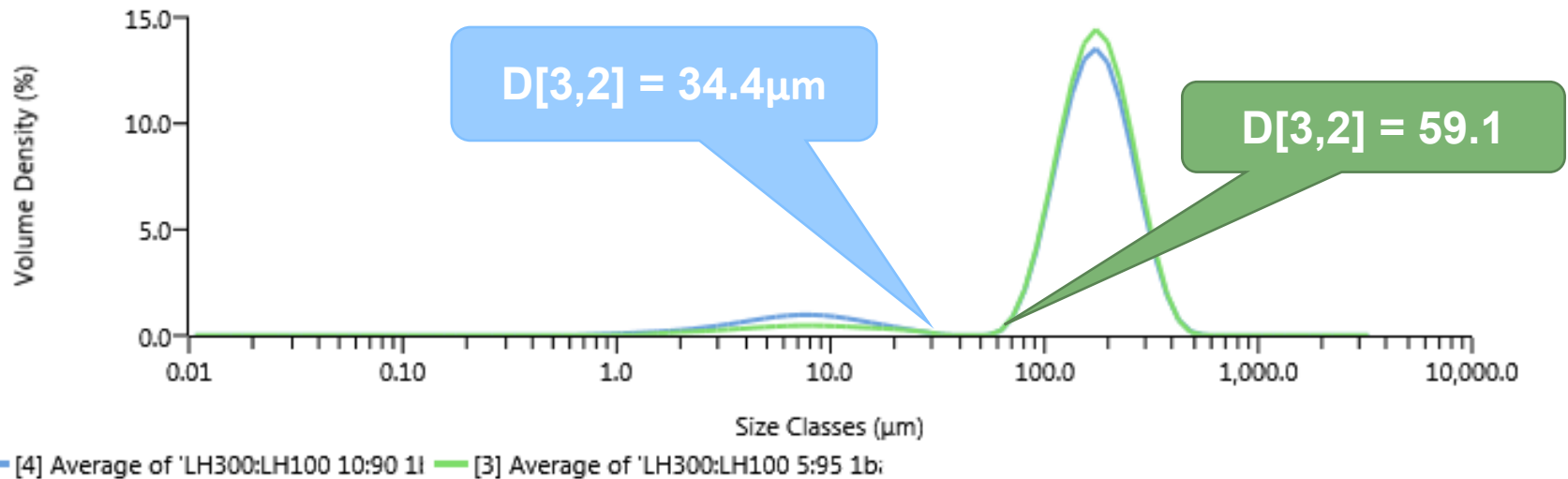
- D[4,3] is sensitive to changes in the coarse particle fraction
 - Useful for monitoring milling or dispersion



| Record Number | Sample Name | Dx (10) (μm) | Dx (50) (μm) | Dx (90) (μm) | D [3,2] (μm) | D [4,3] (μm) |
|---------------|---------------------|---------------------------|---------------------------|---------------------------|---------------------------|---------------------------|
| 7 | sample 1 ultrasound | 2.55 | 4.92 | 30.6 | 4.78 | 11.2 |
| 8 | sample 1 ultrasound | 2.38 | 4.21 | 20.0 | 4.11 | 7.95 |

PSD Statistics: Surface area weighted mean

- D[3,2] is sensitive to changes in the fine particle fraction
 - Useful when surface area is important

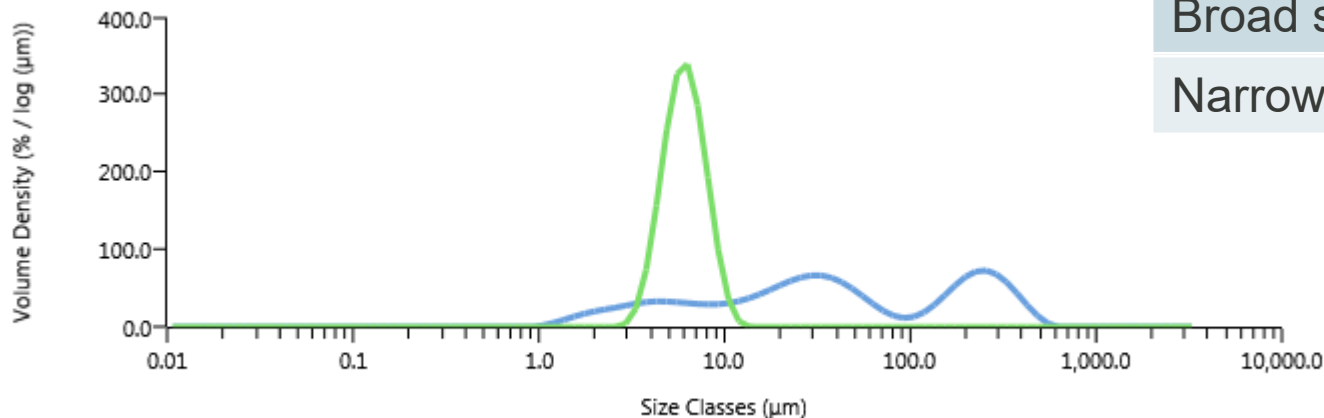


| Record Number | Sample Name | Dx (10) (µm) | Dx (50) (µm) | Dx (90) (µm) | D [3,2] (µm) | D [4,3] (µm) |
|---------------|-------------------------------------|--------------|--------------|--------------|--------------|--------------|
| 4 | Average of 'LH300:LH100 10:90 1bar' | 14.9 | 162 | 275 | 34.4 | 165 |
| 3 | Average of 'LH300:LH100 5:95 1bar' | 90.3 | 168 | 279 | 59.1 | 175 |

PSD Statistics: Other parameters

- Span is a measure of distribution width

$$\text{Span} = \frac{(Dv90 - Dv10)}{Dv50}$$



— Average of 'broad sample 10s' — Average of 'Narrow sample'

| | Span |
|---------------|-------|
| Broad sample | 8.619 |
| Narrow sample | 0.695 |

PSD Statistics: Other parameters

- The Uniformity describes distribution spread (has a lower value for narrower distributions)

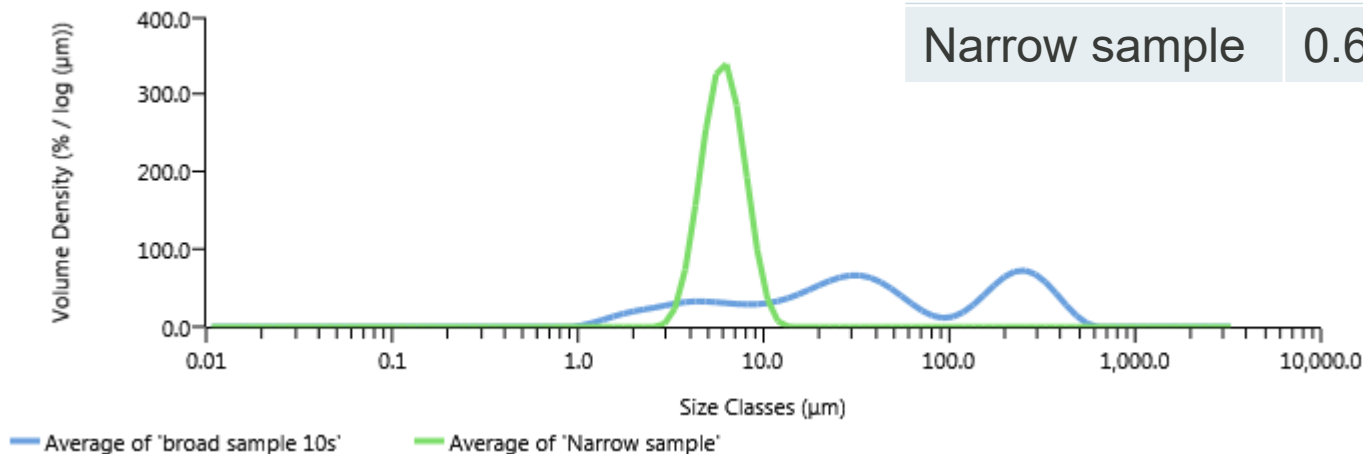
$$\text{Uniformity} = \frac{\sum V_i |Dv50 - D_i|}{Dv50 \sum V_i}$$

$Dv50$ = median

D_i = diameter of size class

V_i = volume in size class

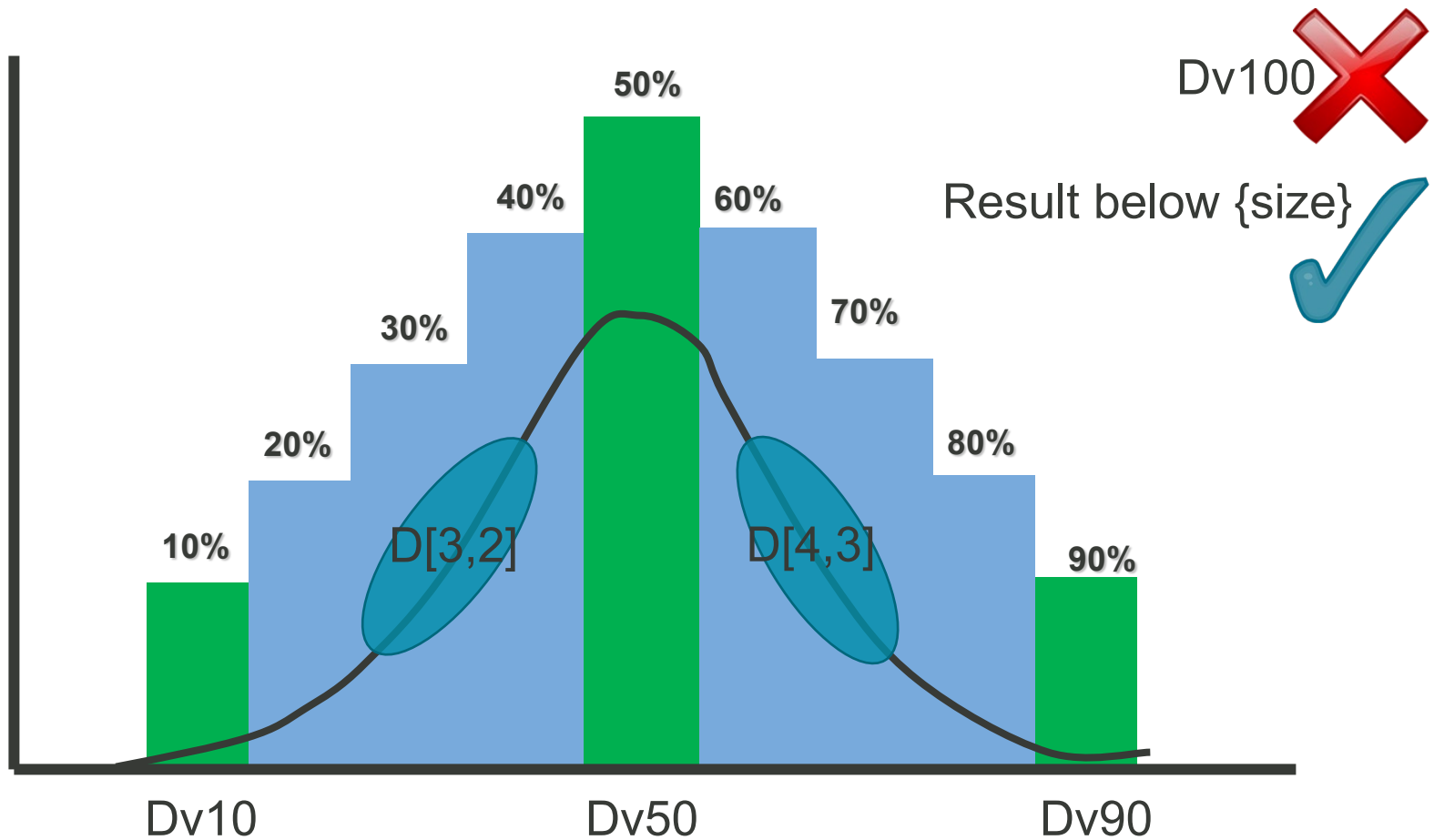
| | Span | Uniformity |
|---------------|-------|------------|
| Broad sample | 8.619 | 2.525 |
| Narrow sample | 0.695 | 0.213 |



PSD Statistics: Other parameters

- **Specific surface area**
 - Total surface area of particles in the sample divided by their total weight
 - Requires the density of the material to be input on the Material page of the SOP/record (default value is 1g/cc)
- **Concentration**
 - Volume concentration (%) is calculated using the Beer-Lambert law

Chose a parameter that is sensitive to the changes you wish to monitor

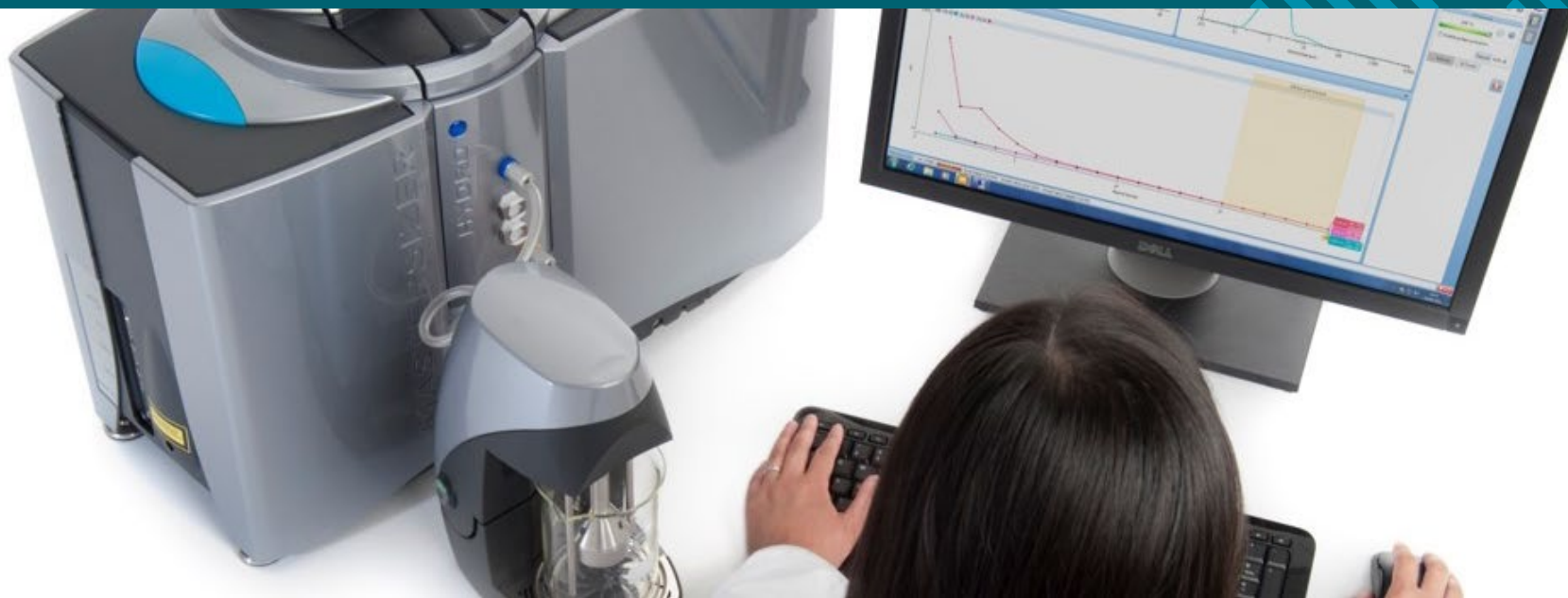




**Malvern
Panalytical**

Mastersizer 3000 Part 3:

Method development for wet or liquid dispersions



The purpose of method development

- A laser diffraction measurement requires

‘a representative sample, dispersed at an adequate concentration in a suitable liquid or gas’

<USP429>

- Method development must define appropriate
 - Sampling
 - Dispersion
 - Measurement conditions

What is the biggest source of error?

Fine particles

Coarse particles

Choice of dispersant

Sampling

Sonication
(dispersion)

Measurement
time

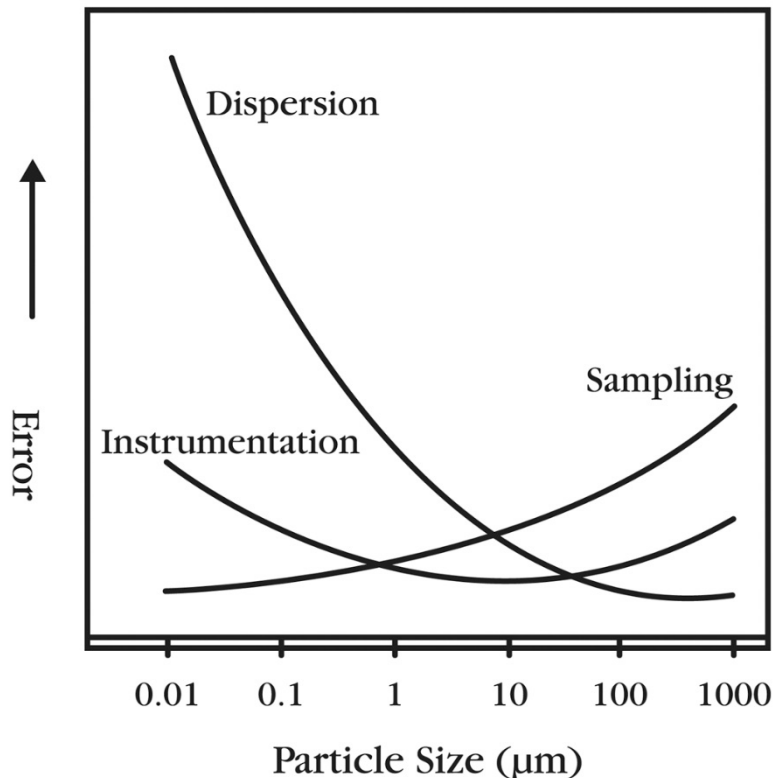
Optical
properties

Dispersant

Sampling

Optical
properties

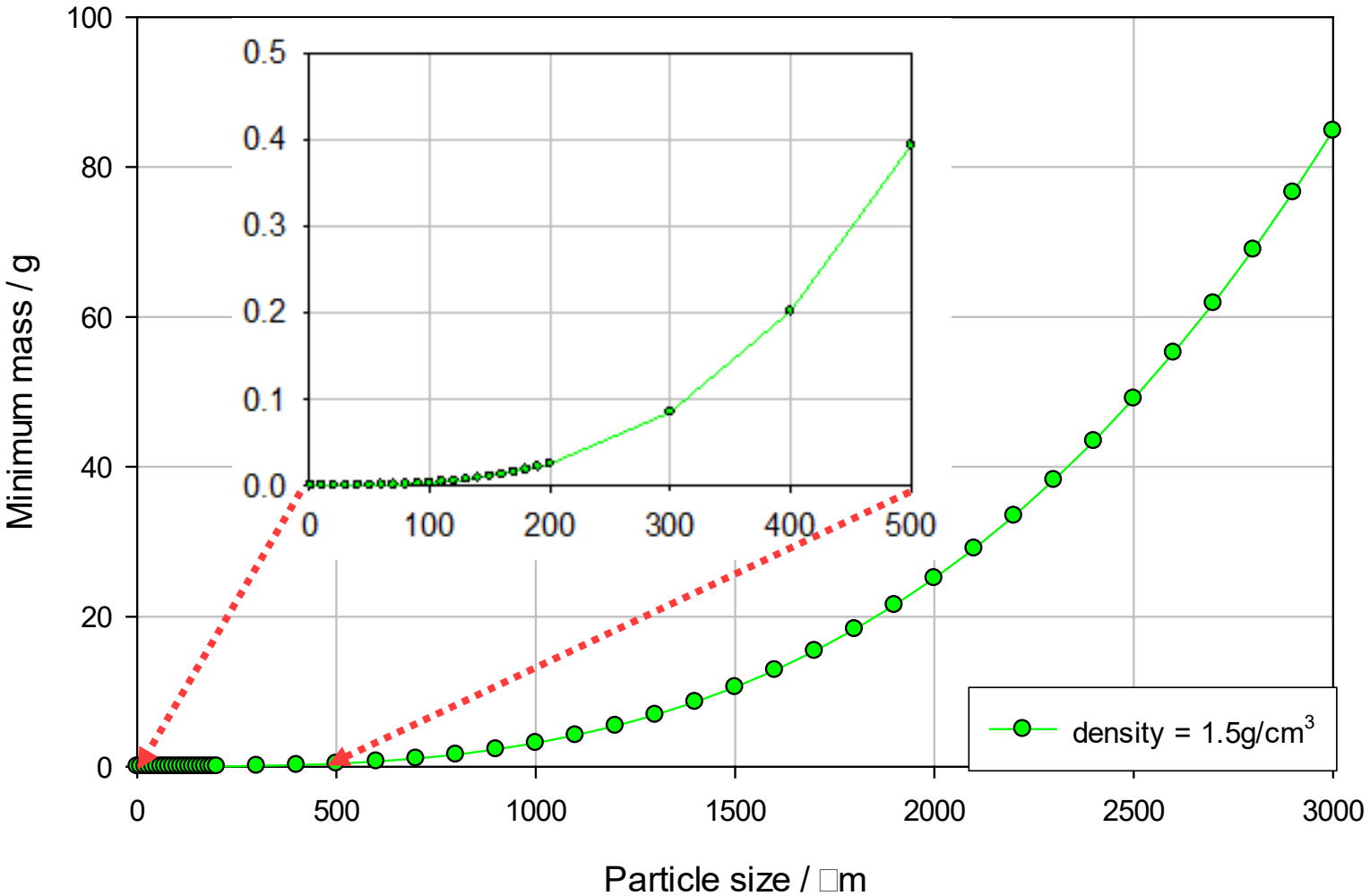
What do we need to control when making measurements?



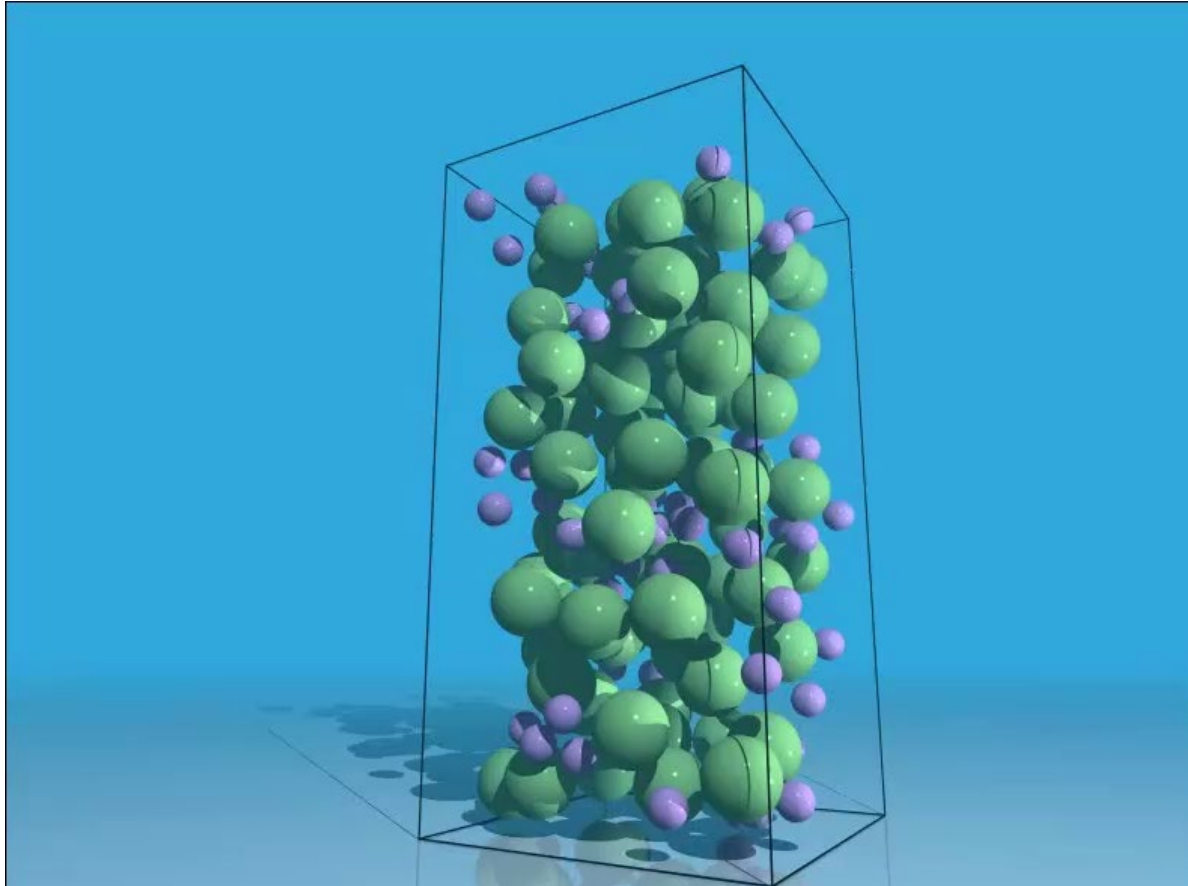
“Novices in the size measurement field must understand that most errors in size measurement arise through poor **sampling** and **dispersion** and not through instrument inadequacies.”

T. Allen, *Advances in Ceramics, Vol 21: Ceramic Powder Science*, page 721, The American Ceramic Society Inc. (1987)

How much sample do I need to measure?

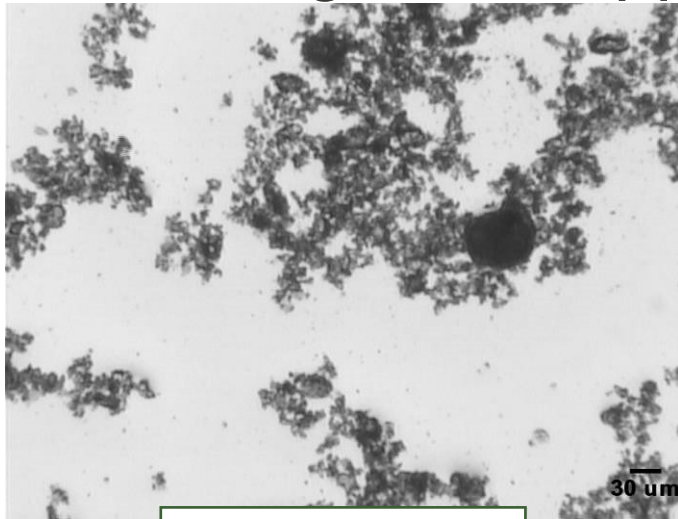


What happens to particles in transit?

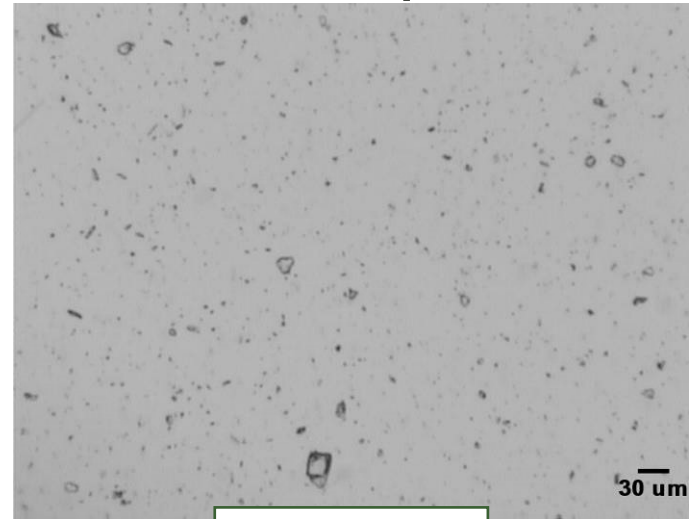
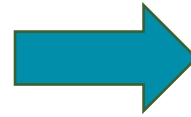


Courtesy of A.J. Morris, M. Glover and M. Probert
© Malvern Panalytical 2017

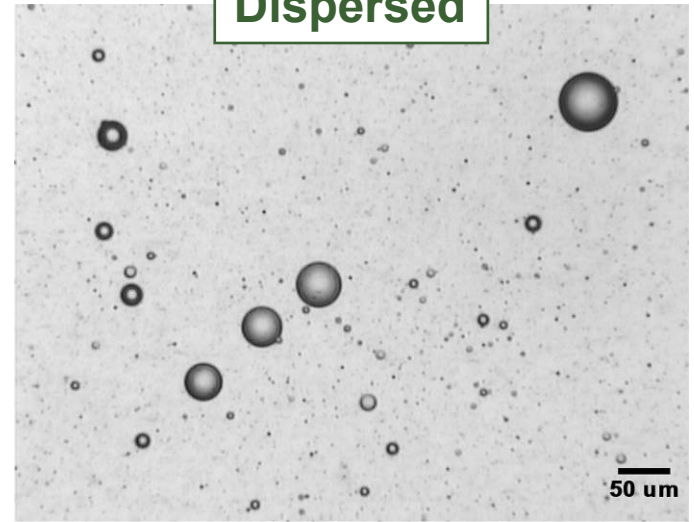
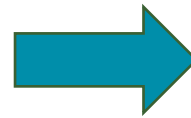
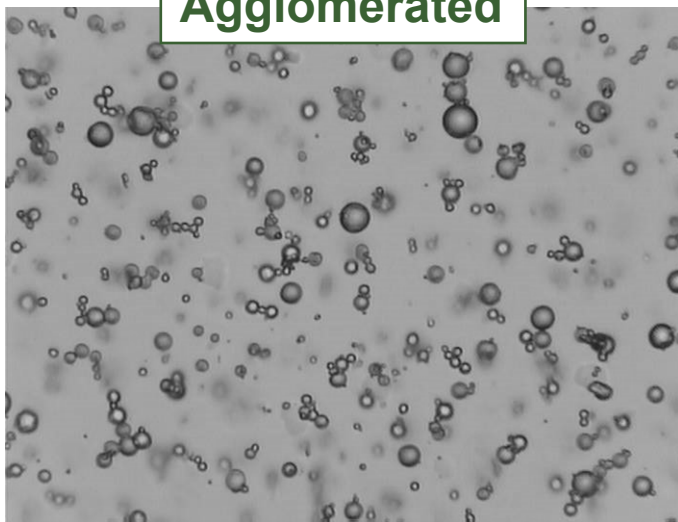
Measuring in the appropriate state of dispersion



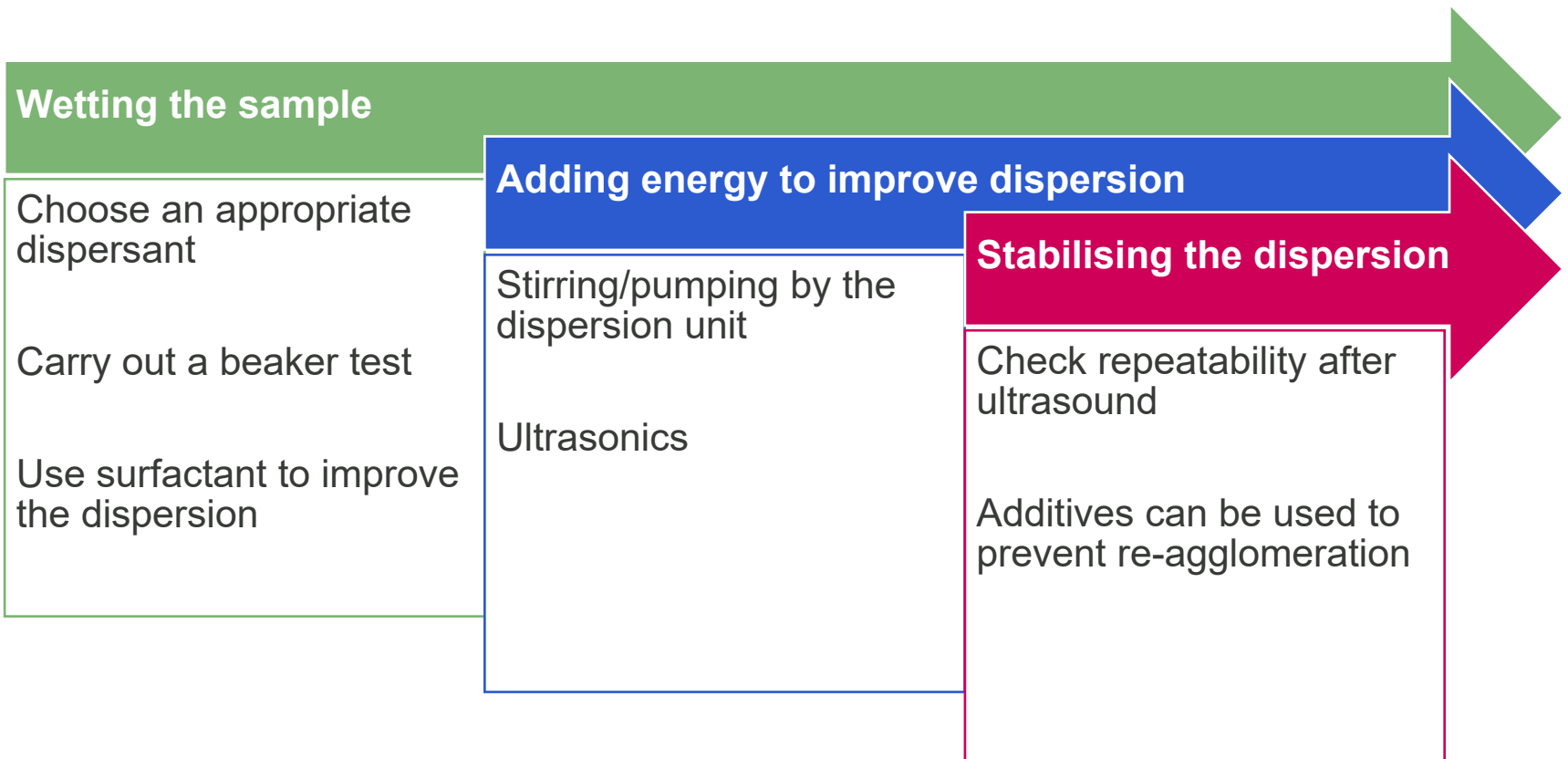
Agglomerated



Dispersed



The wet dispersion process

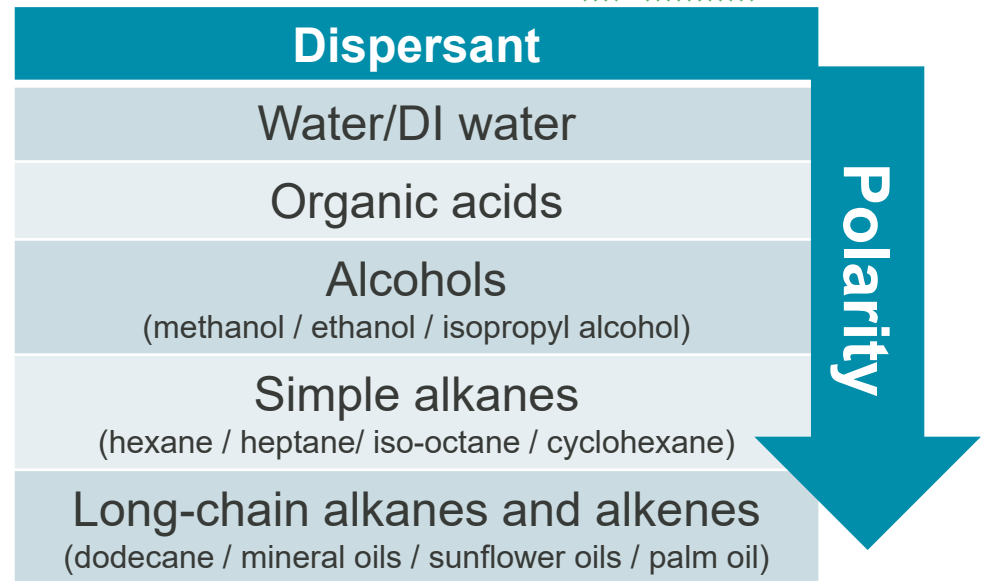


Wetting the sample

Choose an appropriate dispersant

Carry out a beaker test

Use surfactants to improve wetting

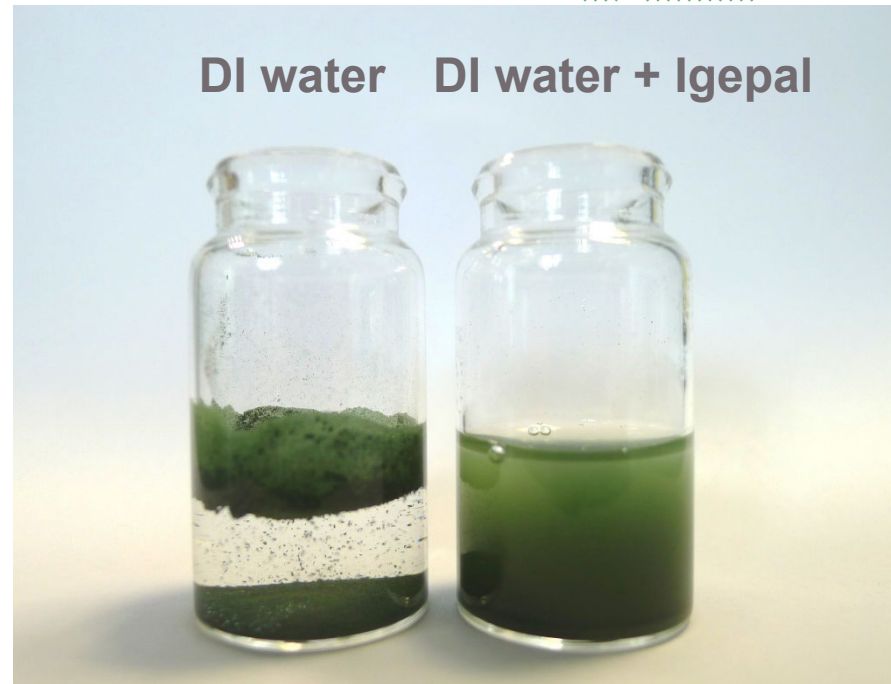


Wetting the sample

Choose an appropriate dispersant

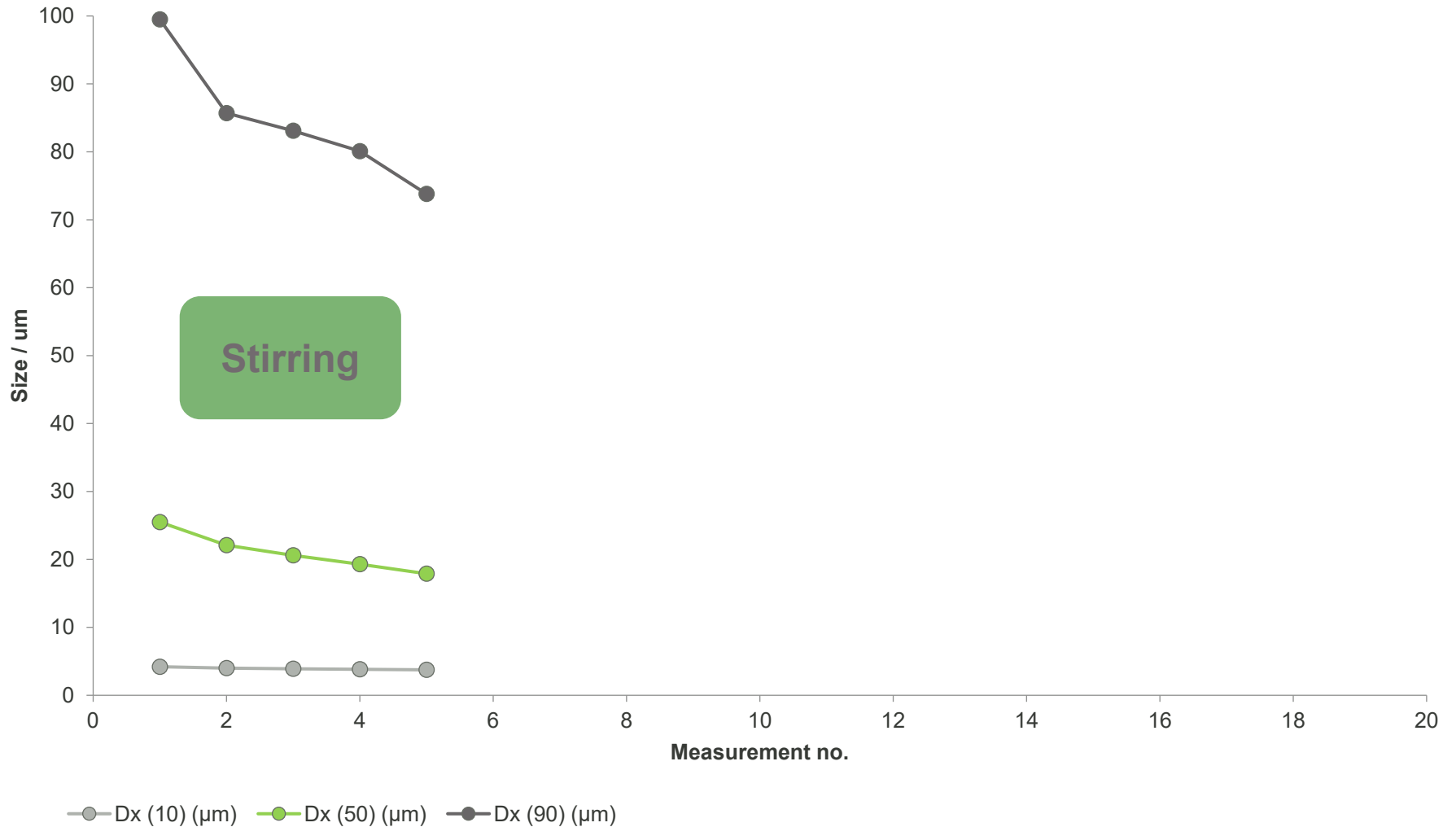
Carry out a beaker test

Use surfactants to improve wetting

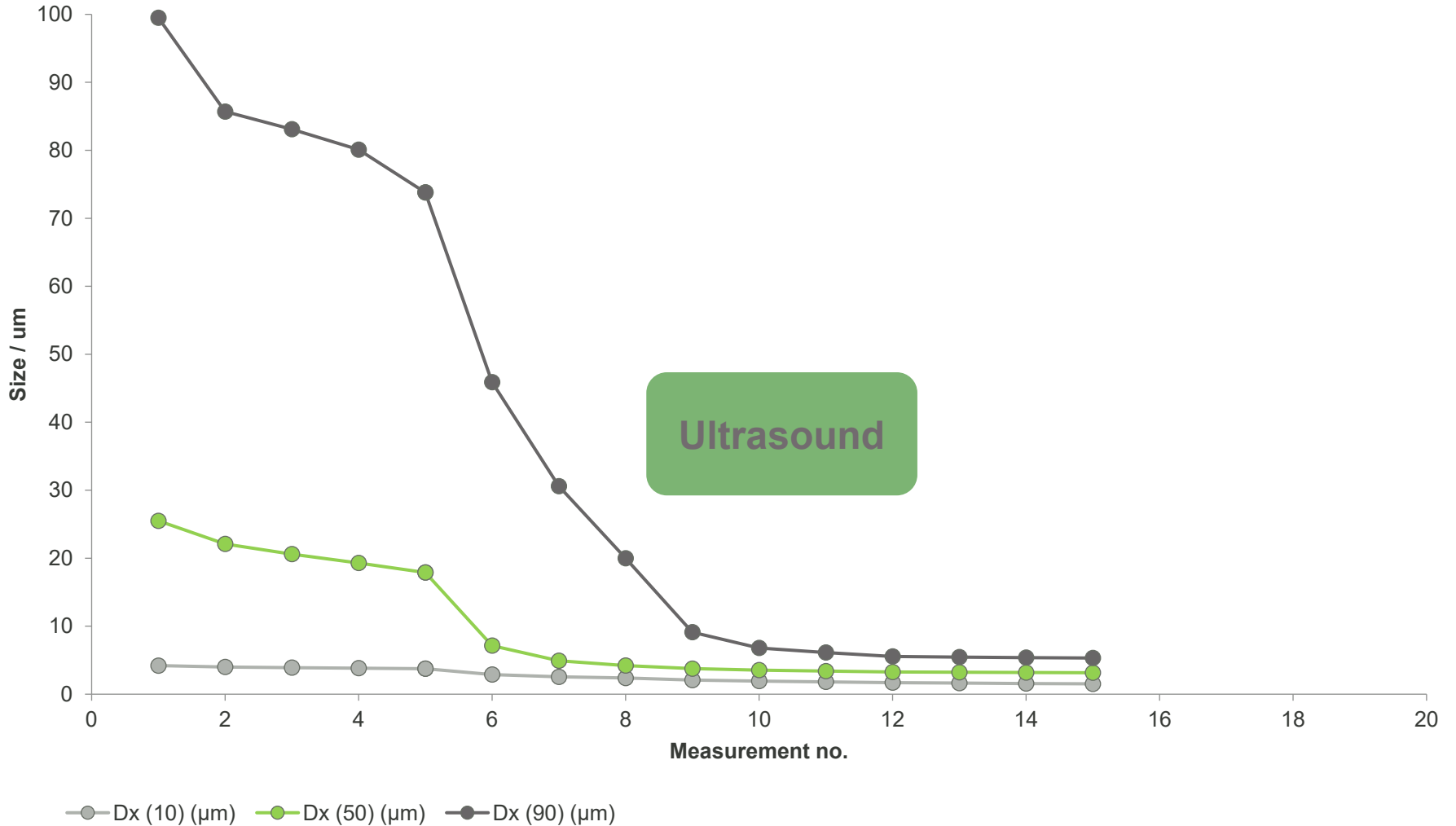


| Stabilization | Examples |
|---------------|---|
| Steric | Igepal CA-360, Tween 20/80, Span 20/80 |
| Electrosteric | Anionic: SDS (sodium dodecylsulfate), AOT (sodium-bis-2-ethylhexylsulfosuccinate) |
| | Cationic: CTAB (cetyltrimethylammonium bromide) |

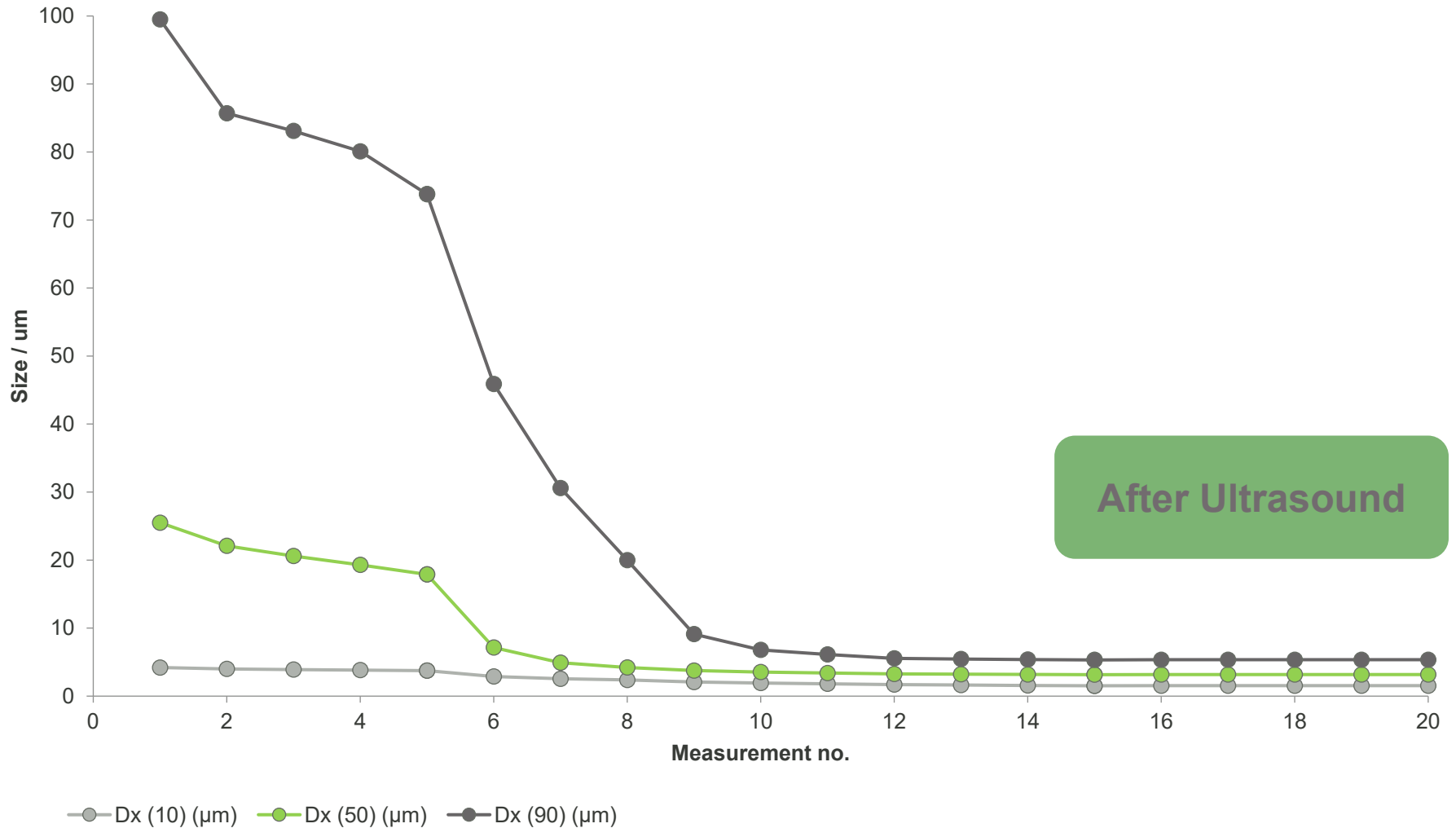
Adding energy to improve dispersion



Adding energy to improve dispersion

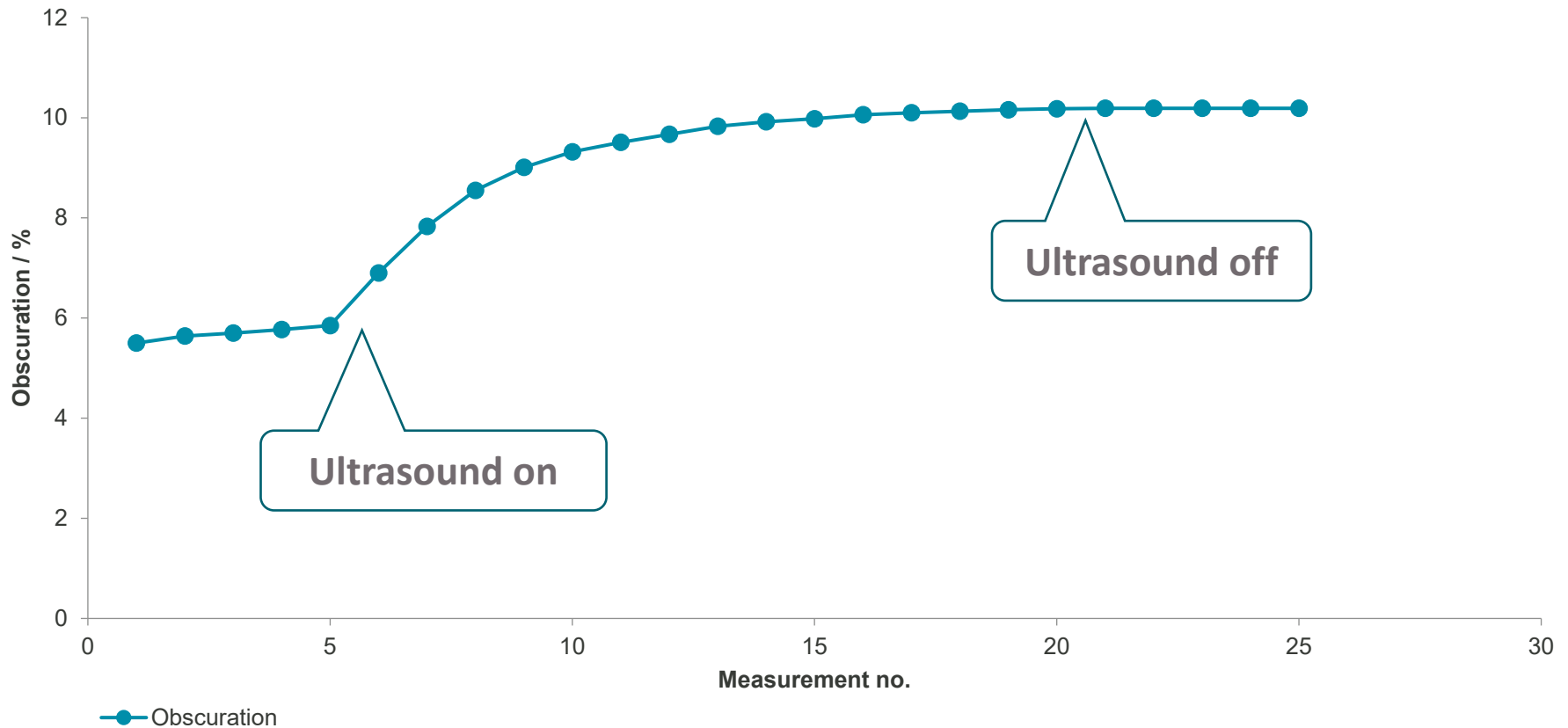


Adding energy to improve dispersion



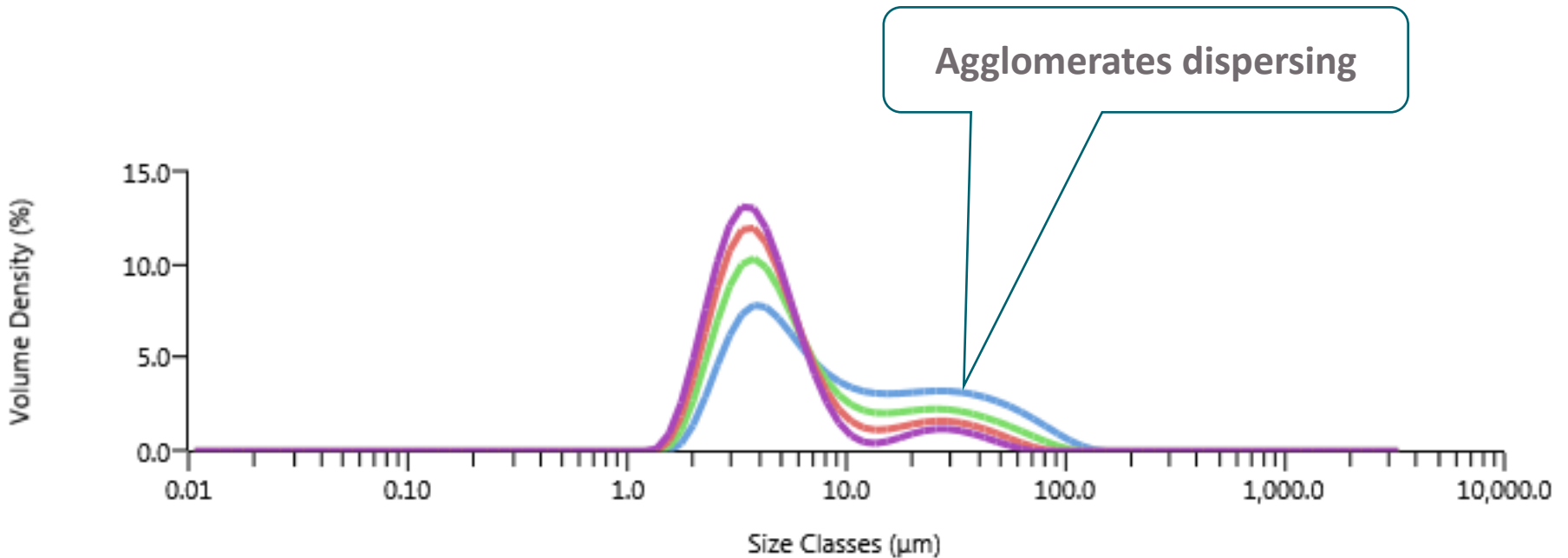
Identifying dispersion: obscuration

- Obscuration increases as agglomerates disperse



Dispersion trend: Particle size distribution

- Overlay the results of an ultrasound titration
 - Should show gradual dispersion



Following trends on the records view: dispersion

Record View

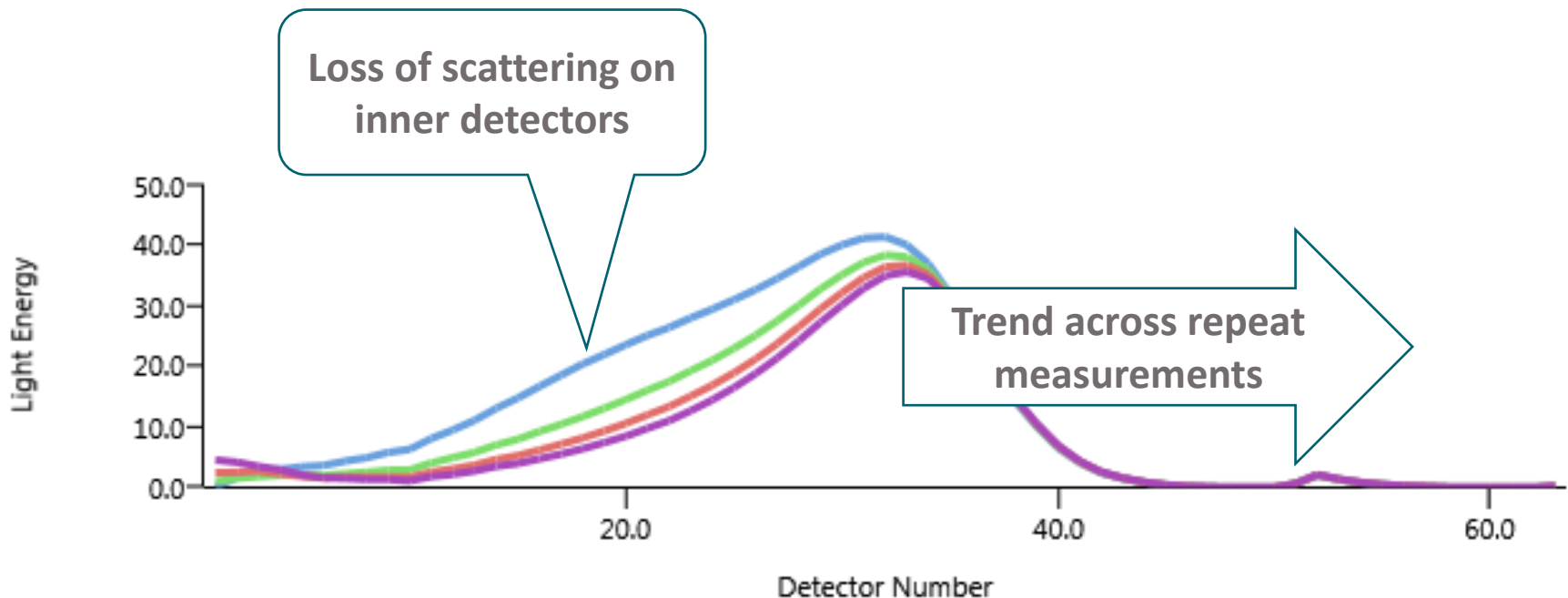
Measurement file 40

Drag column header here to group by that column

| Record Number | Sample Name | Laser Obscuration (%) | Dx (10) (µm) | Dx (50) (µm) | Dx (90) (µm) |
|---------------|-----------------------|-----------------------|--------------|--------------|--------------|
| 1 | Toner | 5.50 | 4.84 | 20.2 | 325 |
| 2 | Toner | 5.64 | 4.86 | 21.2 | 341 |
| 3 | Toner | 5.70 | 4.83 | 19.6 | 329 |
| 4 | Toner | 5.77 | 4.81 | 18.7 | 337 |
| 5 | Toner | 5.85 | 4.86 | 21.7 | 349 |
| 6 | Toner with ultrasound | 6.90 | 4.60 | 11.2 | 295 |
| 7 | Toner with ultrasound | 7.83 | 4.32 | 8.35 | 211 |
| 8 | Toner with ultrasound | 8.55 | 4.18 | 7.44 | 65.3 |
| 9 | Toner with ultrasound | 9.01 | 4.09 | 6.95 | 29.6 |
| 10 | Toner with ultrasound | 9.32 | 4.02 | 6.60 | 12.4 |
| 11 | Toner with ultrasound | 9.51 | 4.17 | 6.48 | 10.9 |
| 12 | Toner with ultrasound | 9.67 | 4.21 | 6.38 | 10.2 |
| 13 | Toner with ultrasound | 9.83 | 4.22 | 6.30 | 9.68 |
| 14 | Toner with ultrasound | 9.92 | 4.20 | 6.24 | 9.49 |
| 15 | Toner with ultrasound | 9.98 | 4.15 | 6.16 | 9.19 |
| 16 | Toner with ultrasound | 10.06 | 4.17 | 6.14 | 9.05 |
| 17 | Toner with ultrasound | 10.10 | 4.22 | 6.12 | 8.82 |
| 18 | Toner with ultrasound | 10.13 | 4.24 | 6.10 | 8.67 |
| 19 | Toner with ultrasound | 10.16 | 4.27 | 6.08 | 8.60 |
| 20 | Toner with ultrasound | 10.18 | 4.28 | 6.07 | 8.51 |

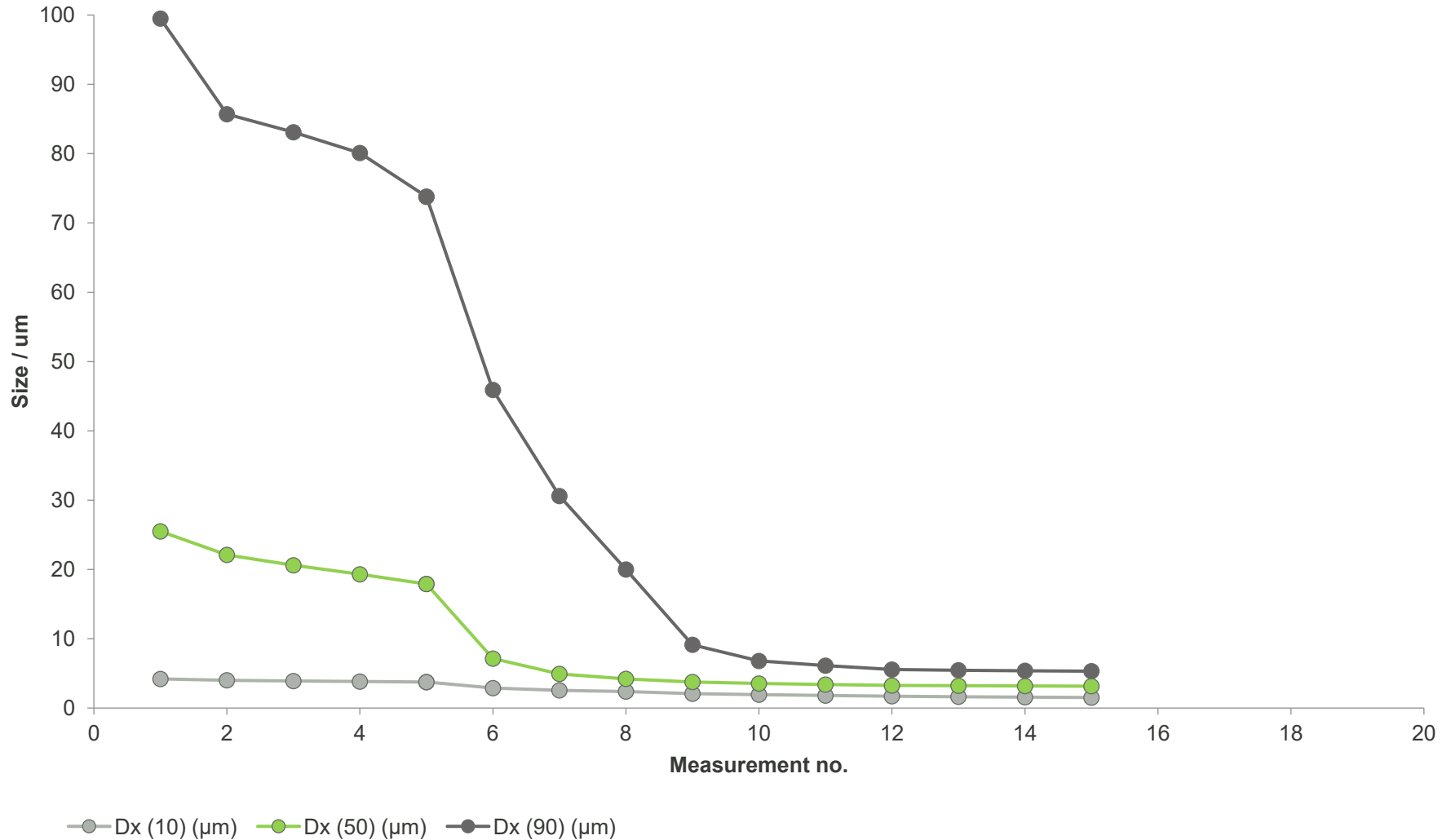
Dispersion trend: scattering data

- During dispersion, as the particles get smaller
 - Scattering on inner detectors decreases
 - Peak shifts to higher angle detectors

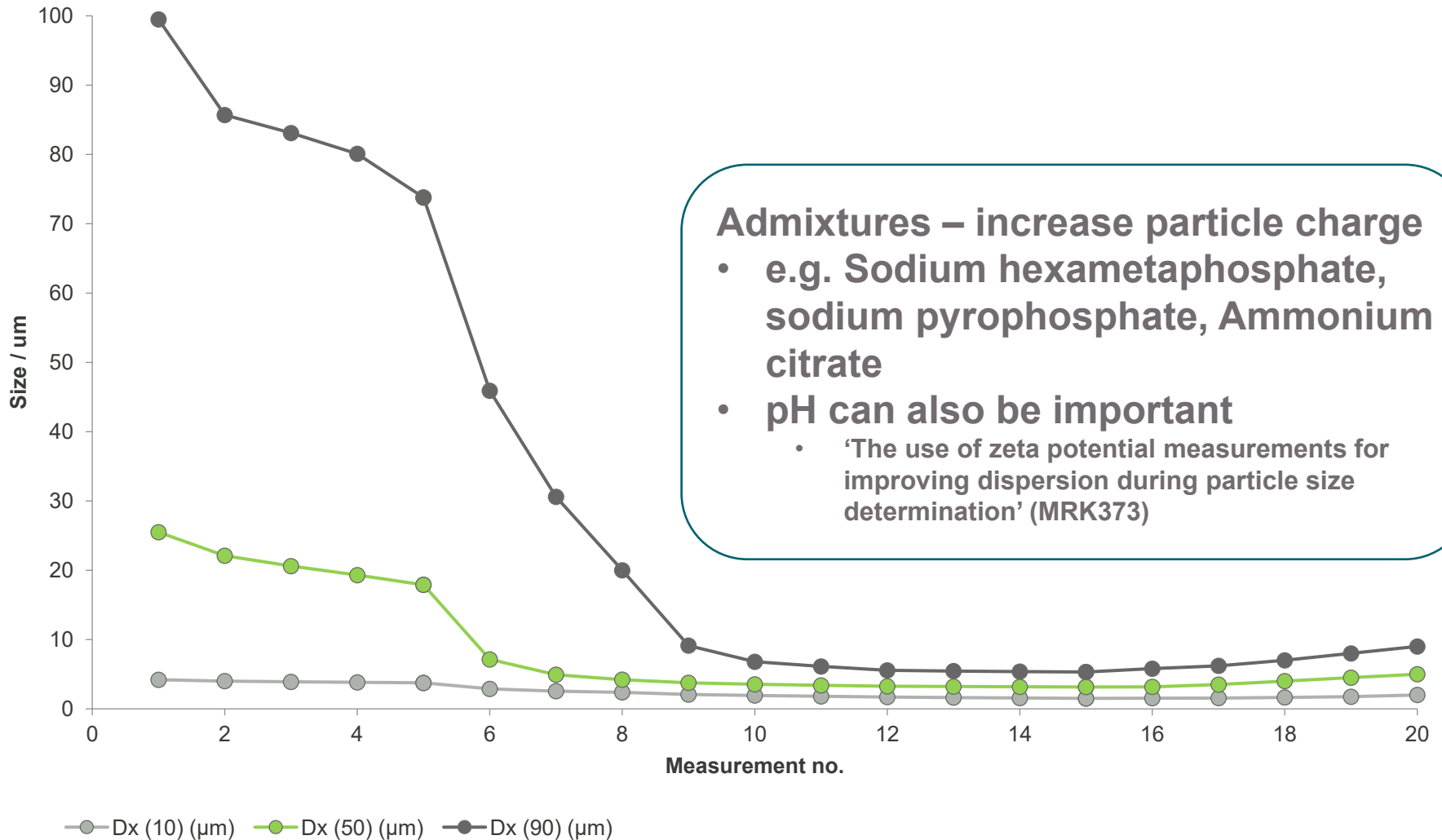


Reminder: detector number increases with angle
Scattering from larger particles falls on low angle detectors

Stabilising the dispersion



Stabilising the dispersion



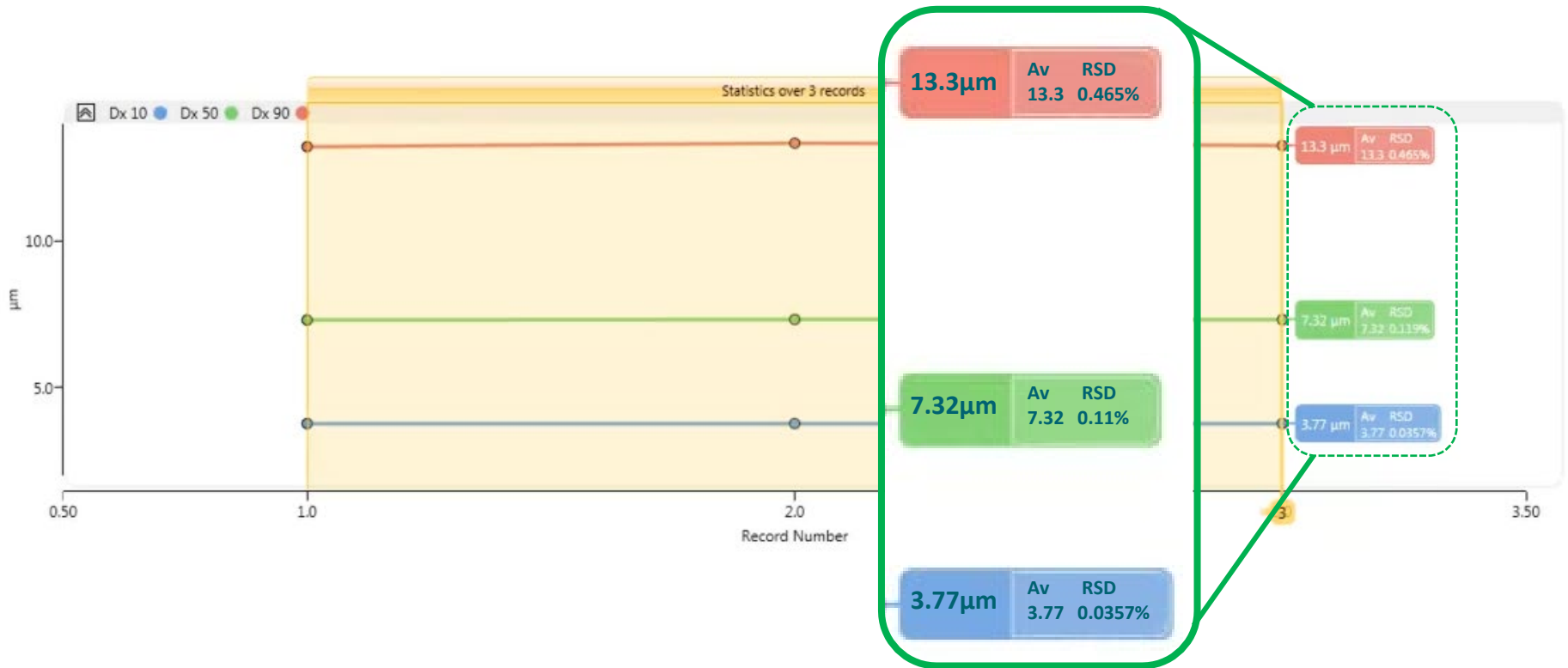
How stable should the results be?

- ISO13320-1: Section 6.4
 - Dv50 - 5 different readings: COV < 3%
 - Dv10 and Dv90: COV < 5%
 - “Below 10µm, these maximum values should be doubled.”
 - **Coefficient of variance = relative standard deviation (RSD)**
- In ideal conditions
 - 0.5% COV on parameters >1µm
 - 1% COV on parameters <1µm

$$\%COV = \frac{\textit{standard deviation}}{\textit{average}} \times 100$$

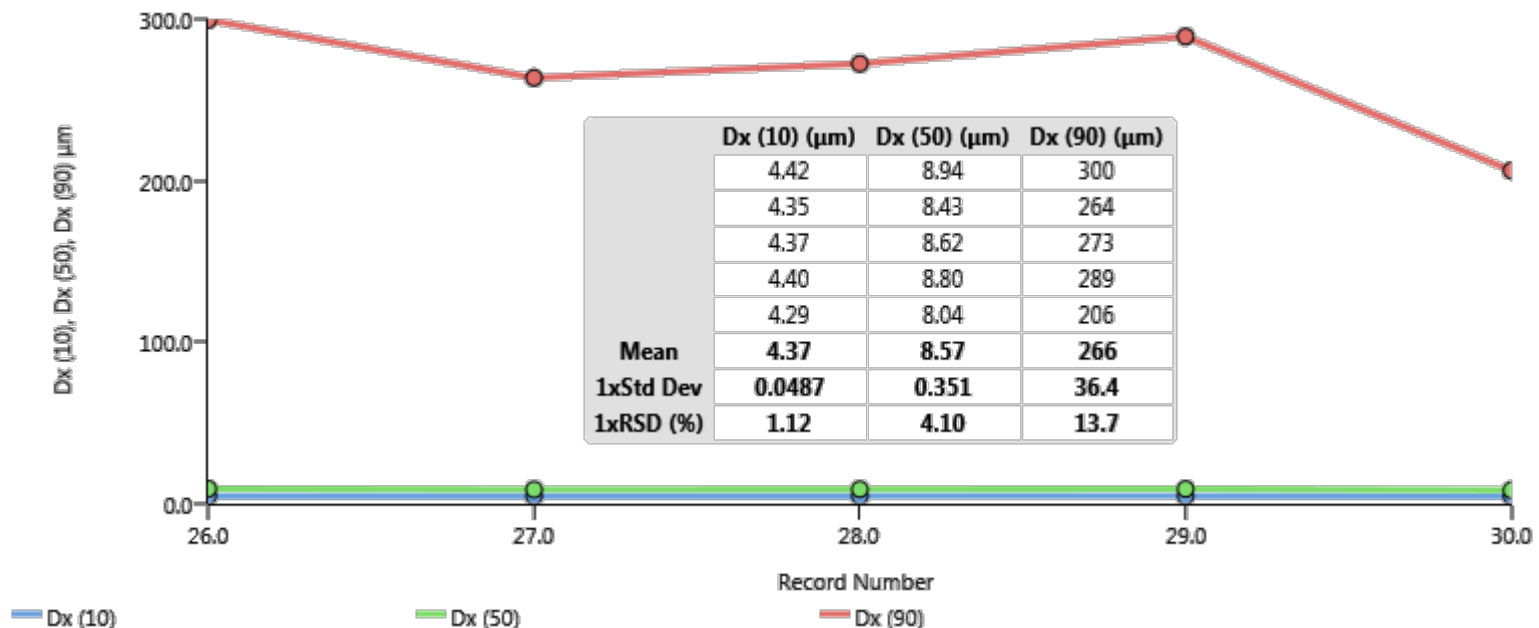
Checking the stability of the results

- The live trend shows the variability of the results
 - RSDs should be within ISO limits



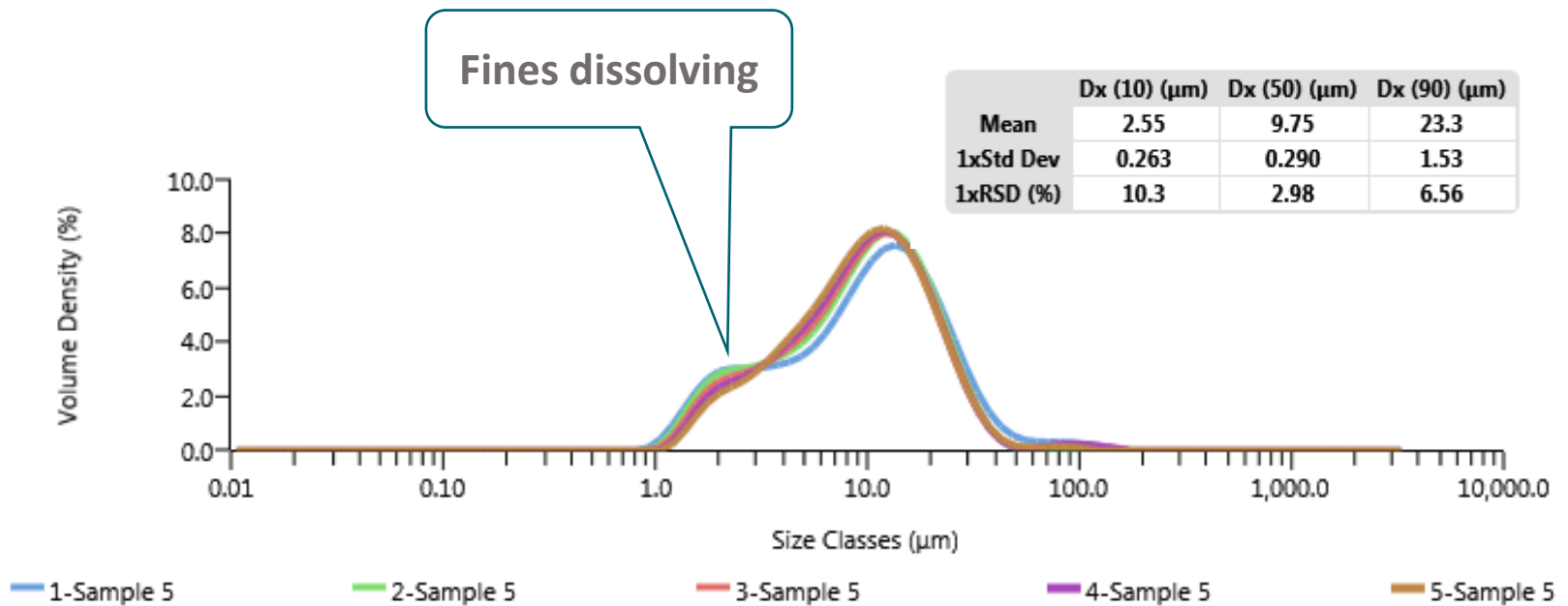
What if the results are not stable?

- If the results are not stable the sample could be:
 - Dissolving
 - Agglomerating
 - Breaking due to excessive ultrasound
 - Incomplete dispersion



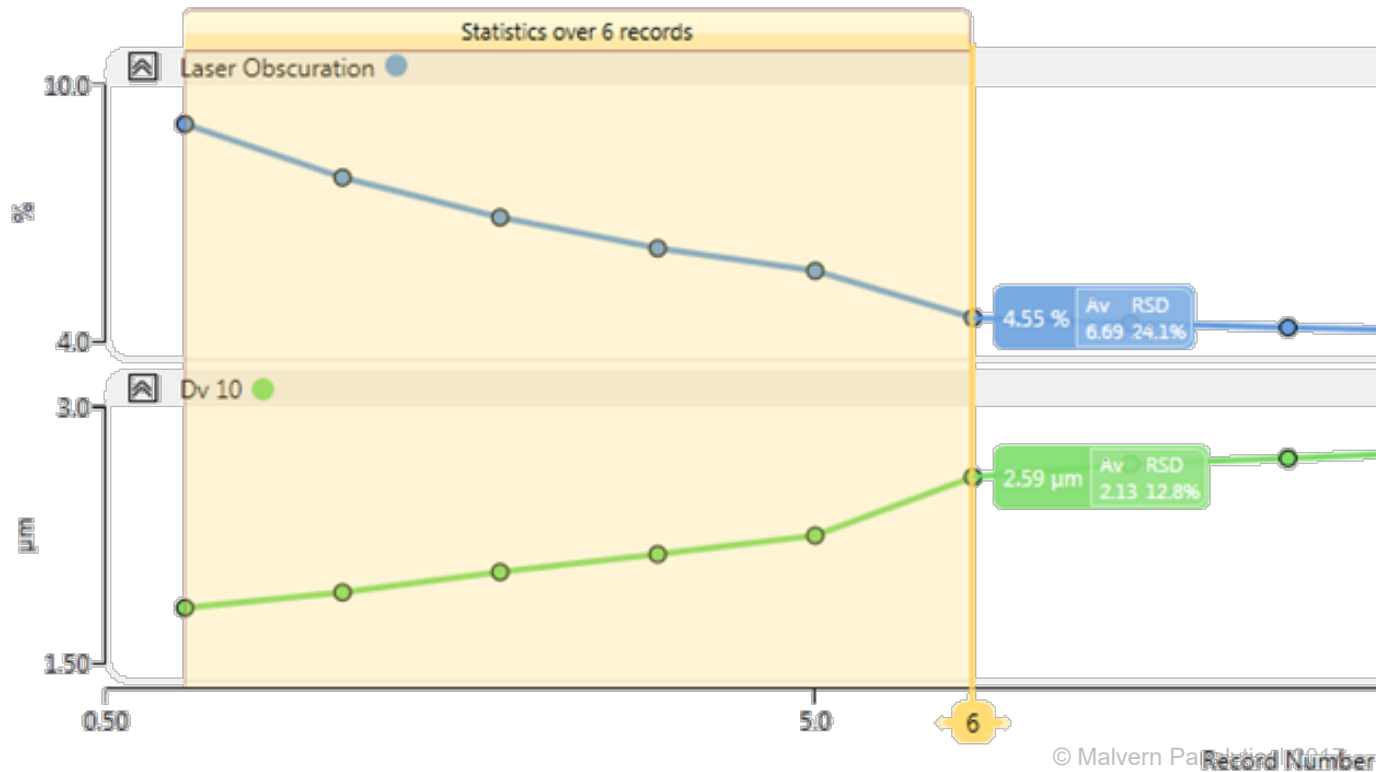
Identifying sample dissolution

- If the particles are dissolving
 - The obscuration will decrease
 - The fine particles will get smaller and then disappear
 - The Dv10 will increase
 - The coarse fraction will become more dominant



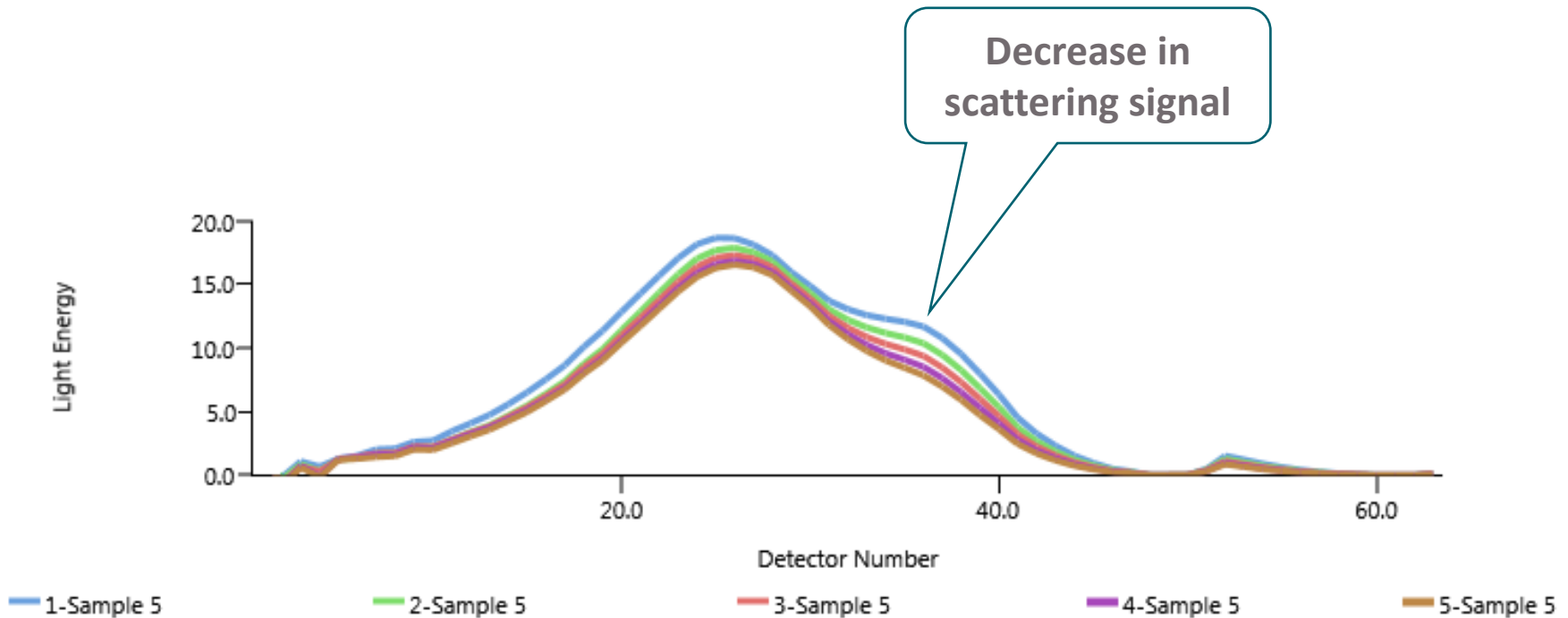
Identifying sample dissolution: Trend view

- Plotting the obscuration and Dv10 on the live trend can help to identify dissolution
 - Decreasing obscuration
 - Increasing Dv10



Identifying sample dissolution: Scattering data

- Scattering signal level decreases as sample dissolves
 - Fewer particles scatter less light



Remedies for sample dissolution

- Try other dispersants
 - As listed on slide 6
- Select the dispersant with the least dissolution
 - Determined from repeatability
- If unavoidable, measurements can be made with slight dissolution
 - Sample prep and measurement time must be controlled
- As a last resort, saturated solutions can be used
 - Must be used carefully
 - Temperature control etc.

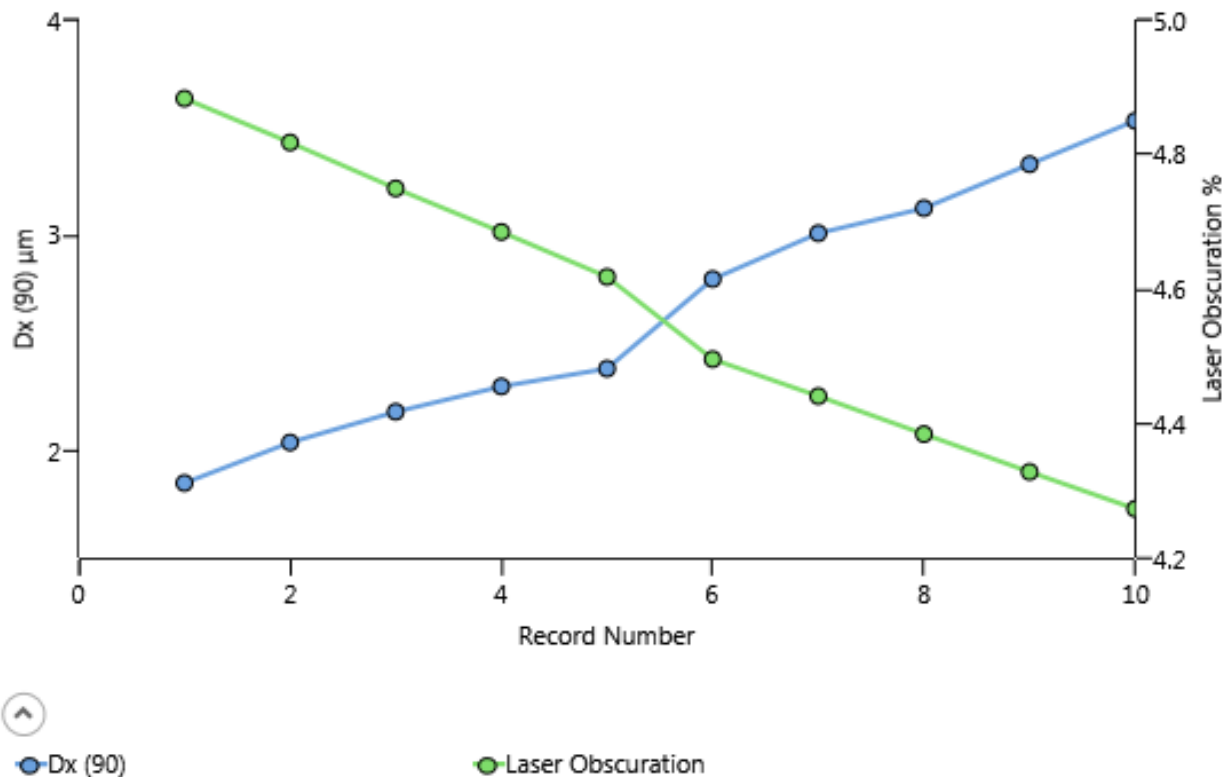
Identifying sample agglomeration

- If the sample is agglomerating
 - A tail of large particles will appear in the distribution
 - The obscuration will decrease gradually
 - As individual particles become agglomerates
 - Using ultrasound may remove the agglomerates
 - If the sample is not chemically stable then ultrasound can cause agglomeration



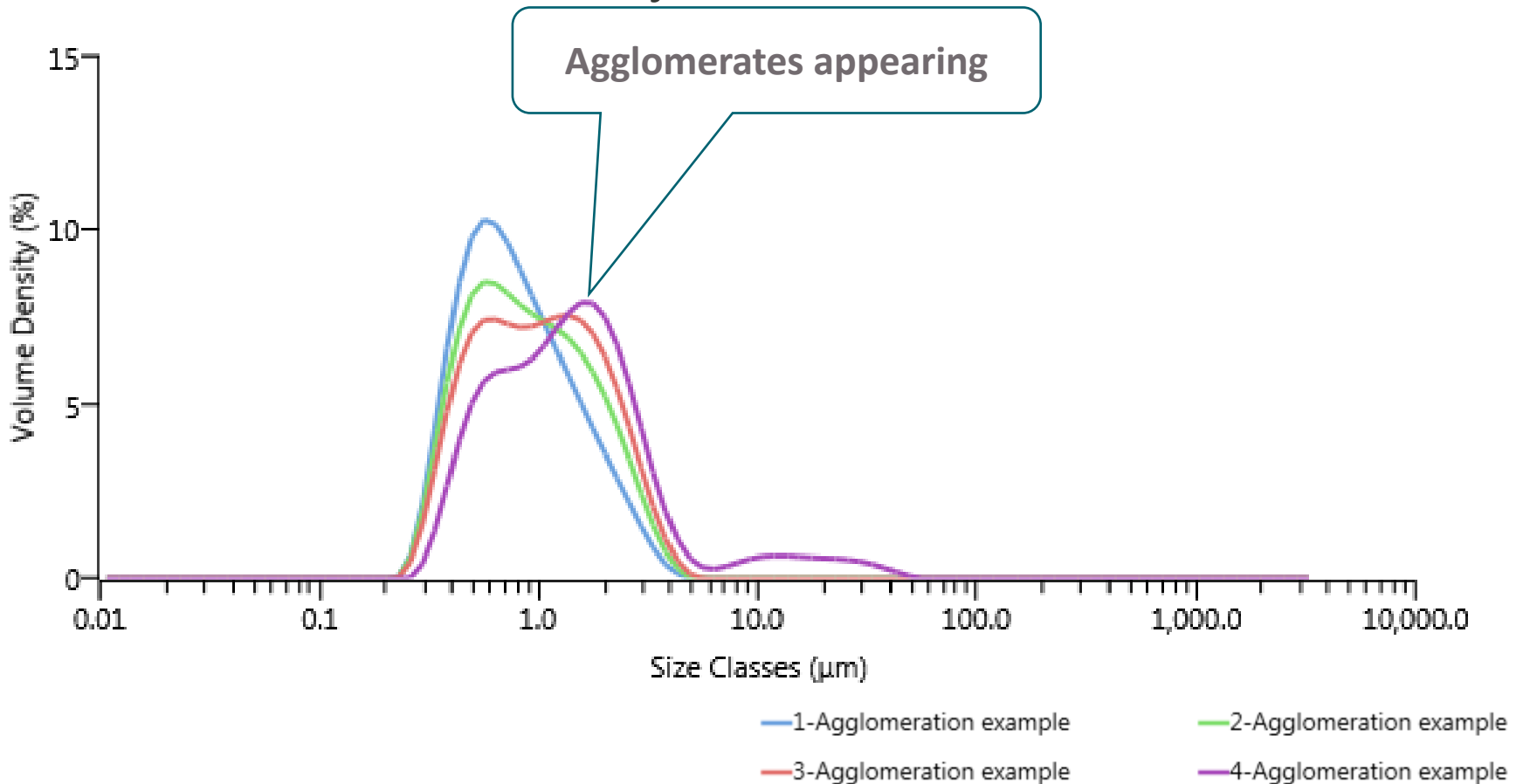
Identifying sample agglomeration: Trend view

- When the sample is agglomerating
 - The obscuration will decrease gradually
 - The Dv90 will increase



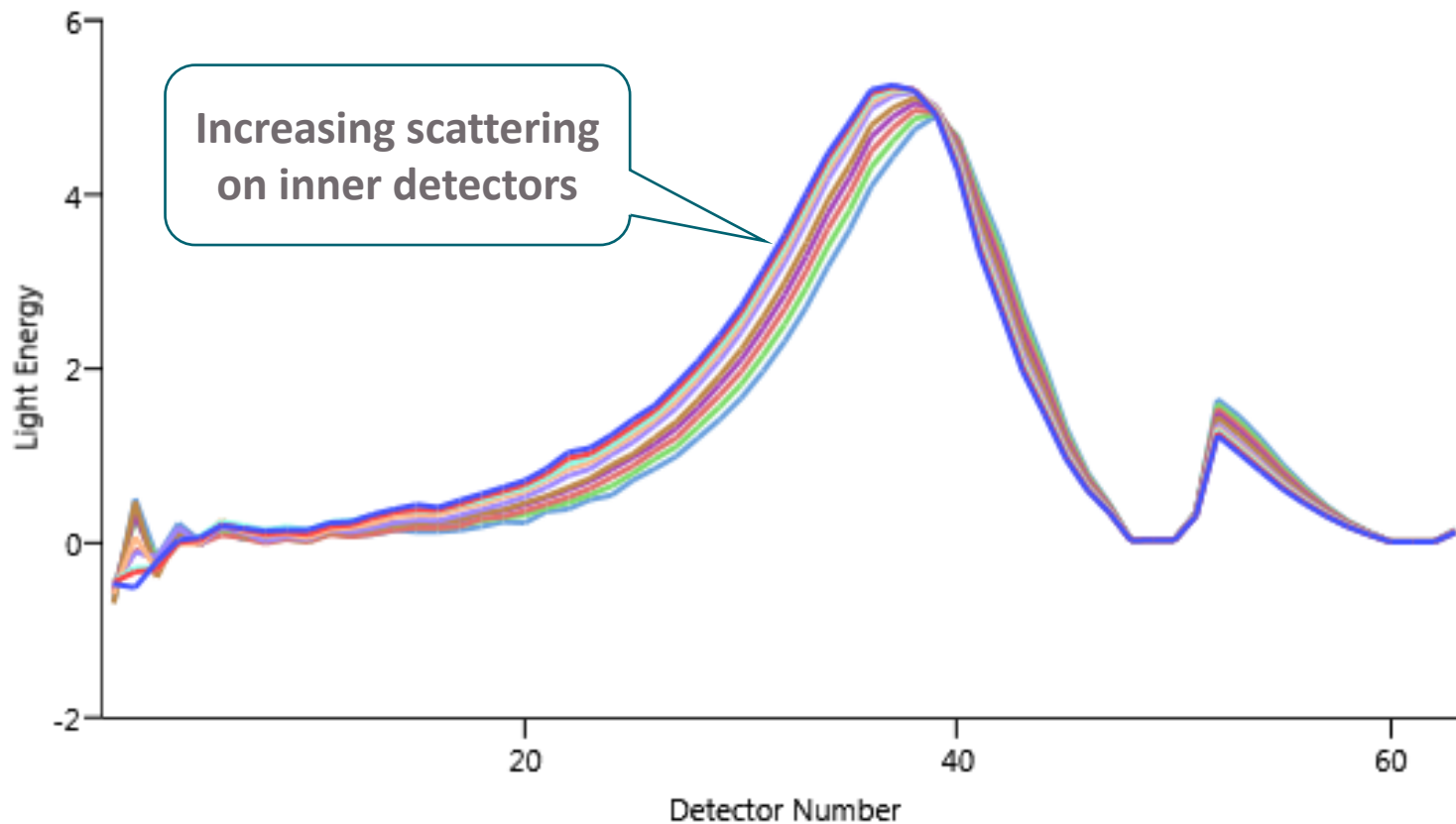
Identifying sample agglomeration: Size distribution

- The volume in the coarse fraction may increase
- Or the coarse fraction may increase in size



Identifying sample agglomeration: Scattering data

- Scattering data will increase and move towards inner detectors.



Remedies for sample agglomeration

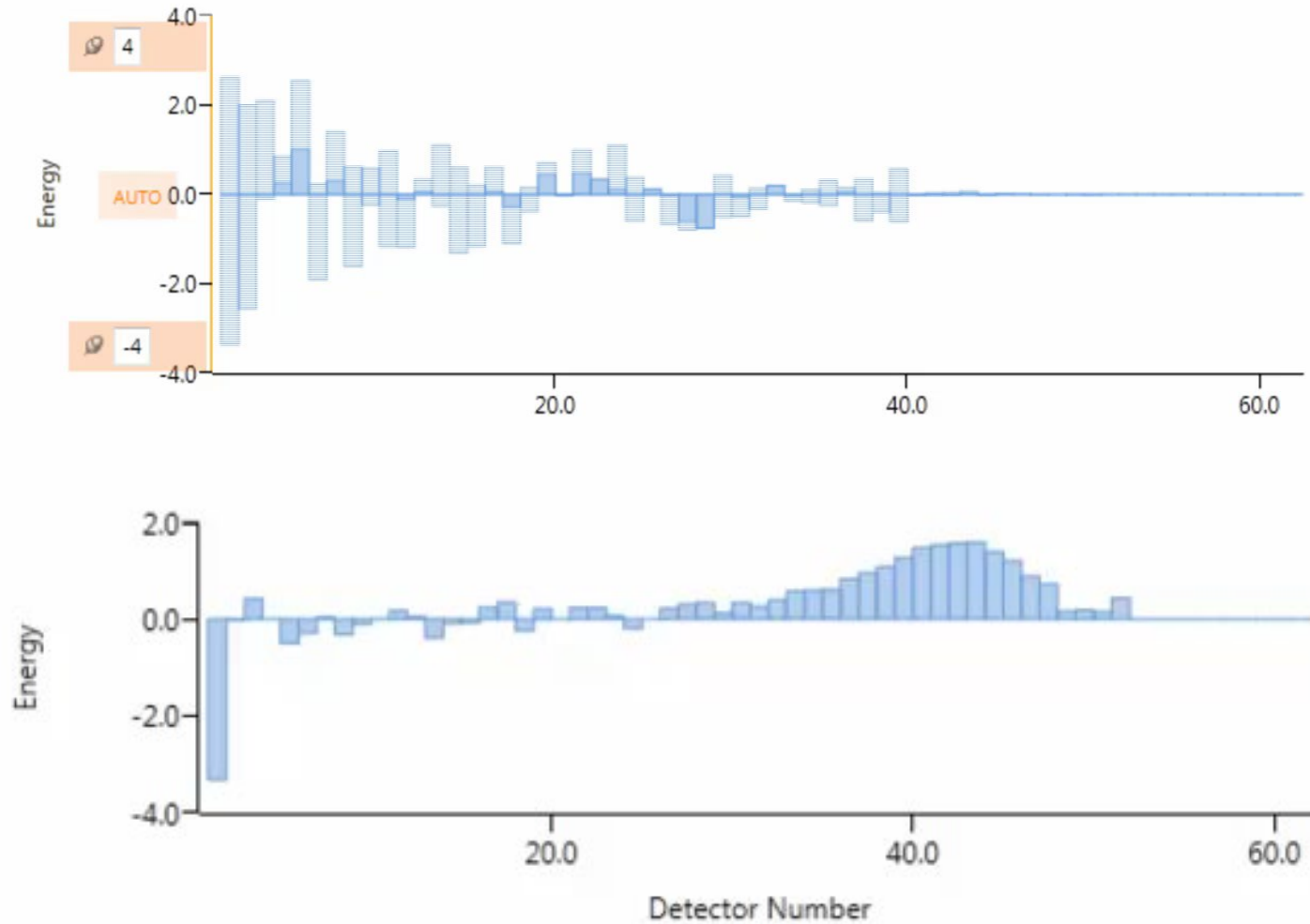
- Surfactants or additives can be used to stabilise dispersions
 - Surfactants decrease interfacial tension
 - List on slide 7
 - Additives increase particle charge
 - Specific adsorption of ions to particle surface
 - E.g. Sodium Hexametaphosphate, Sodium Pyrophosphate, Ammonium Citrate
 - pH can also affect dispersion
- Use low concentrations, a few w/v%
 - as too much can cause agglomeration
 - too much surfactant can also cause foaming



How do measurement conditions affect results

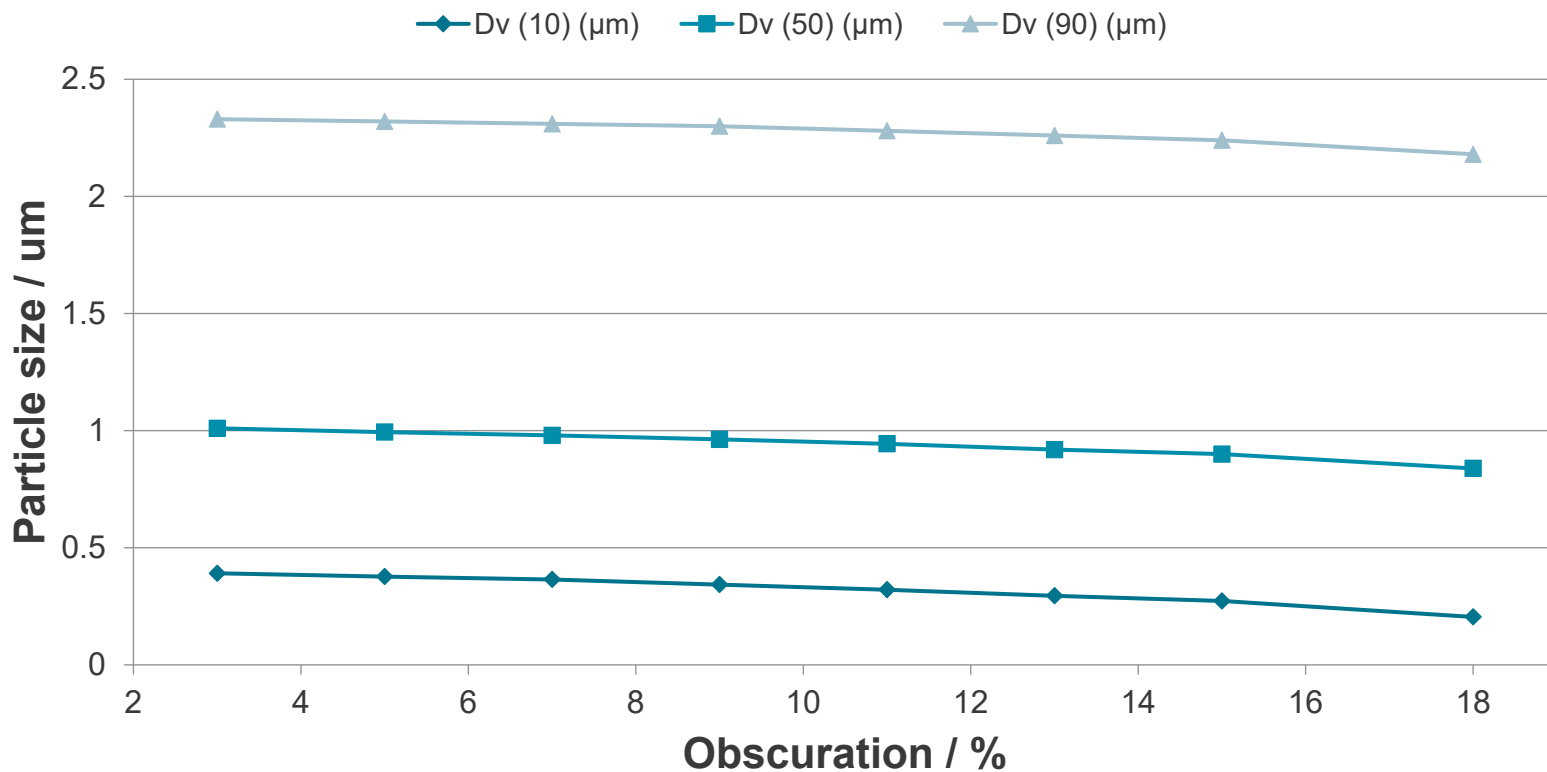
- **Appropriate amount of sample**
 - Good signal to noise ratio
 - Avoid multiple scattering
- **Correct stir speed**
 - Fast enough to prevent sedimentation for large/dense particles
 - Slow enough not to break emulsions
- **Correct measurement duration**
 - Long enough to sample all of the particles in the dispersion unit

Low obscuration limit: signal to noise ratio

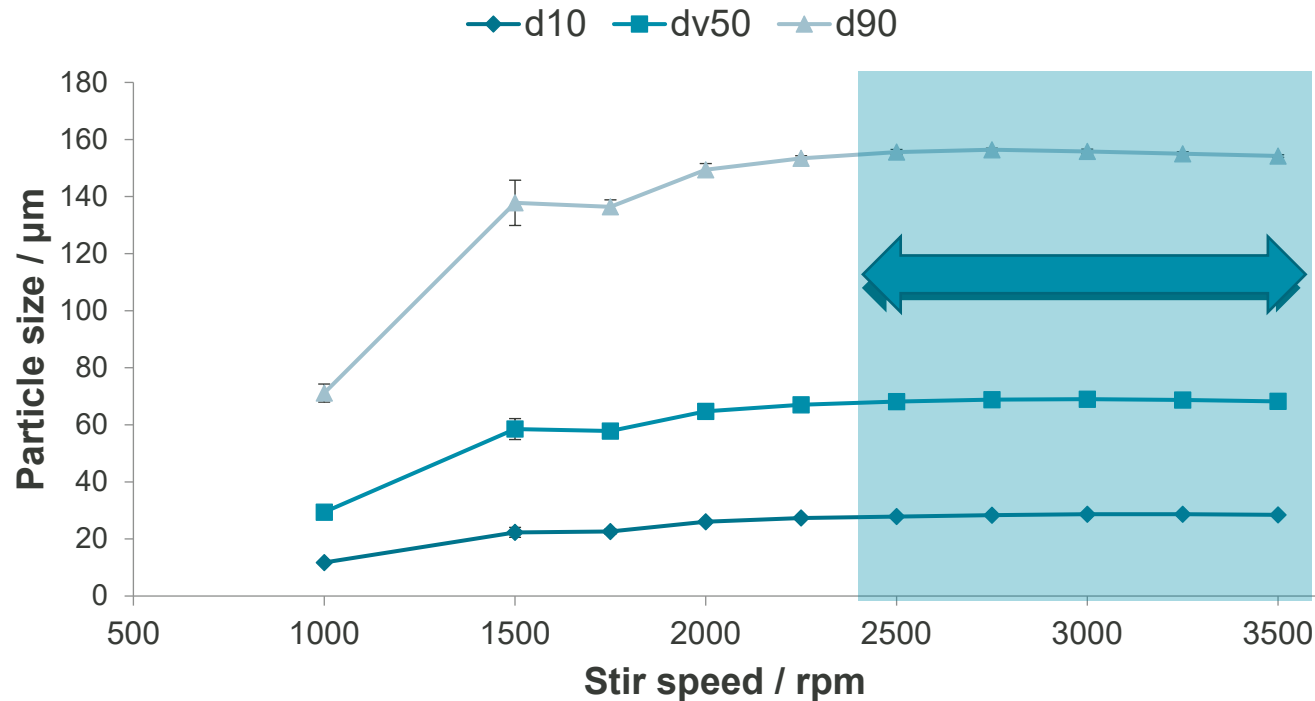


High obscuration limit: multiple scattering

- The upper limit of the obscuration range depends on multiple scattering:
 - Sample should be measured in the range where size is stable with obscuration.

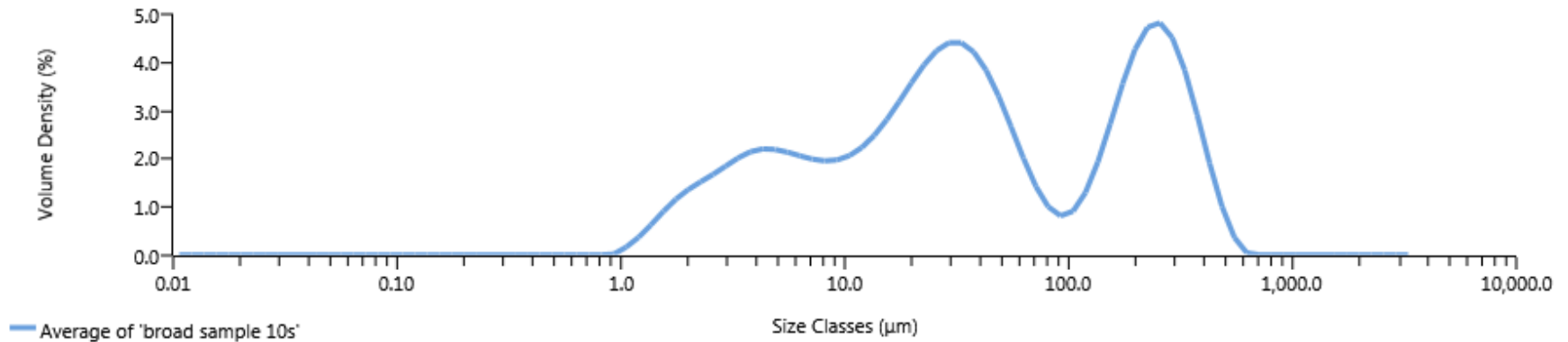


Determine the correct stir speed



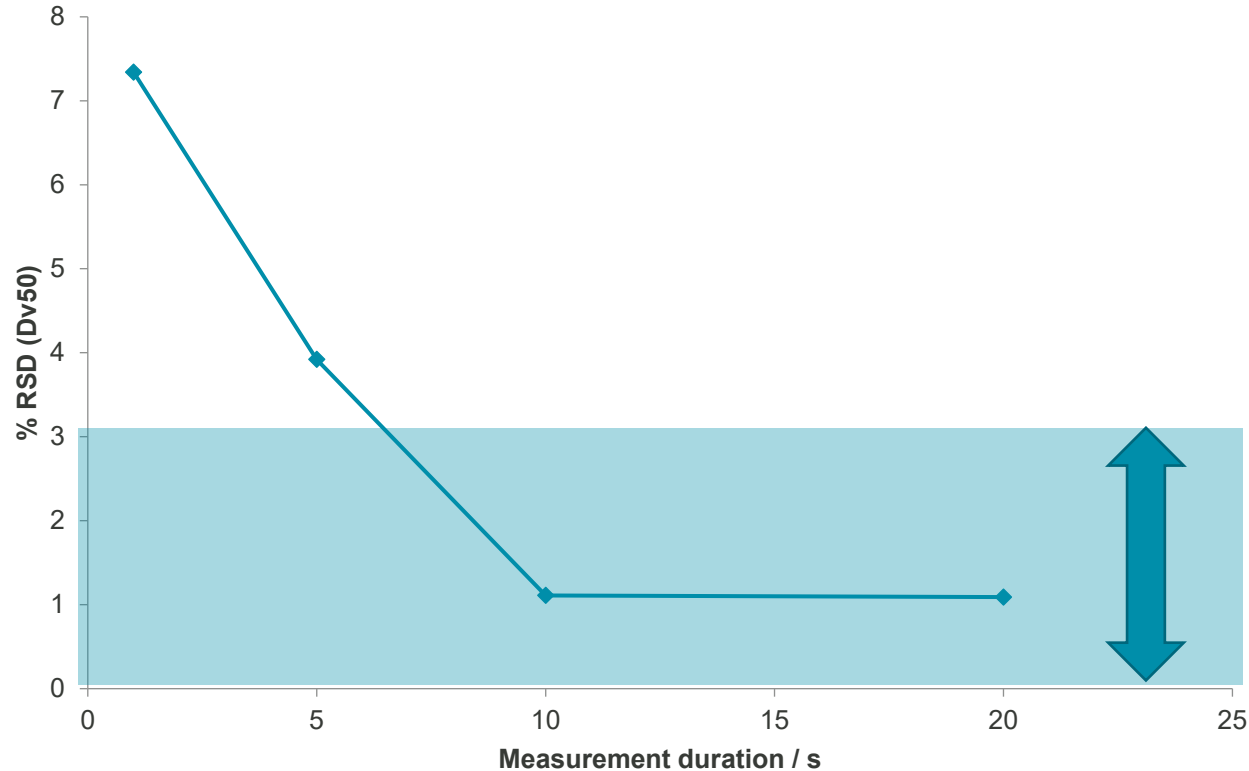
- For coarse or dense materials particle size will increase with stir speed until all particles are suspended
 - A stable particle size is obtained above 2500rpm

Determine the correct measurement duration



- For broad distributions measurement duration must be sufficient to sample all particles in the system.

Affect of measurement duration on variability



- Result variability is reduced as measurement duration is increased
 - Variability is within ISO limits when duration $\geq 10s$

General rules for good measurements

- Measurement and background times
 - Background duration should be at least as long as the measurement duration
 - Make short repeat measurements before, during and after ultrasound to establish dispersion
 - For coarse or polydisperse materials the measurement duration may need to be increased to improve repeatability
- **Ultrasound: generates heat in the dispersant**
 - In organic solvents, allow a pre-measurement delay to let this heat dissipate

The purpose of method development

- A laser diffraction measurement requires

‘a representative sample, dispersed at an adequate concentration in a suitable liquid or gas’

<USP429>

- Method development must define appropriate
 - Sampling
 - Dispersion
 - Measurement conditions



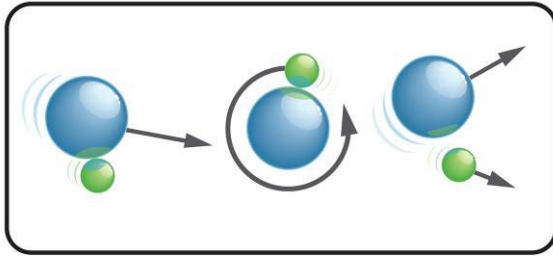
**Malvern
Panalytical**

Mastersizer 3000 Part 4:

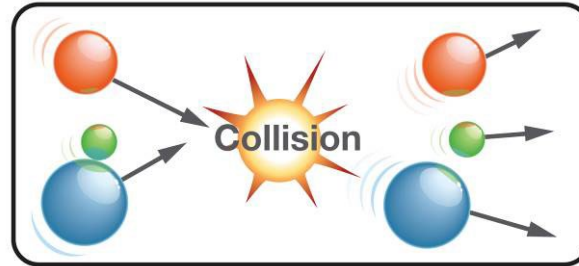
Method development for dry powder dispersion



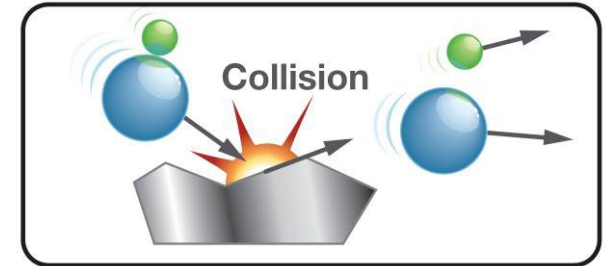
Dry powder dispersion: Mechanisms



Velocity gradients caused by shear stress



Particle-to-particle collisions



Particle-to-wall collisions

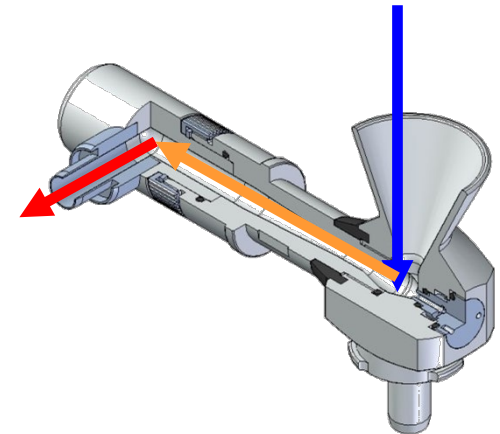
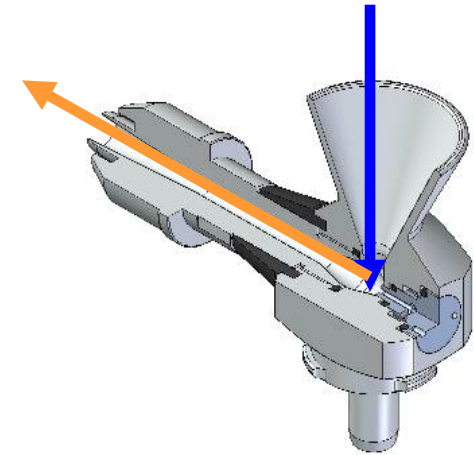
Energy/aggression



- Importance of each mechanism depends on:
 - Disperser geometry
 - Flow rate or pressure drop
 - Material type
- Higher impact energies may improve the dispersion effectiveness
 - Needs to be balanced against the risk of particle break-up

Dry powder dispersion: Disperser design

- Standard disperser
 - Straight through design
 - No direct wall impaction
 - Suitable for most types of sample
- High energy disperser
 - Elbow design
 - Direct impaction surface
 - Suitable for robust aggregated samples



Setting up the feed rate: sample tray

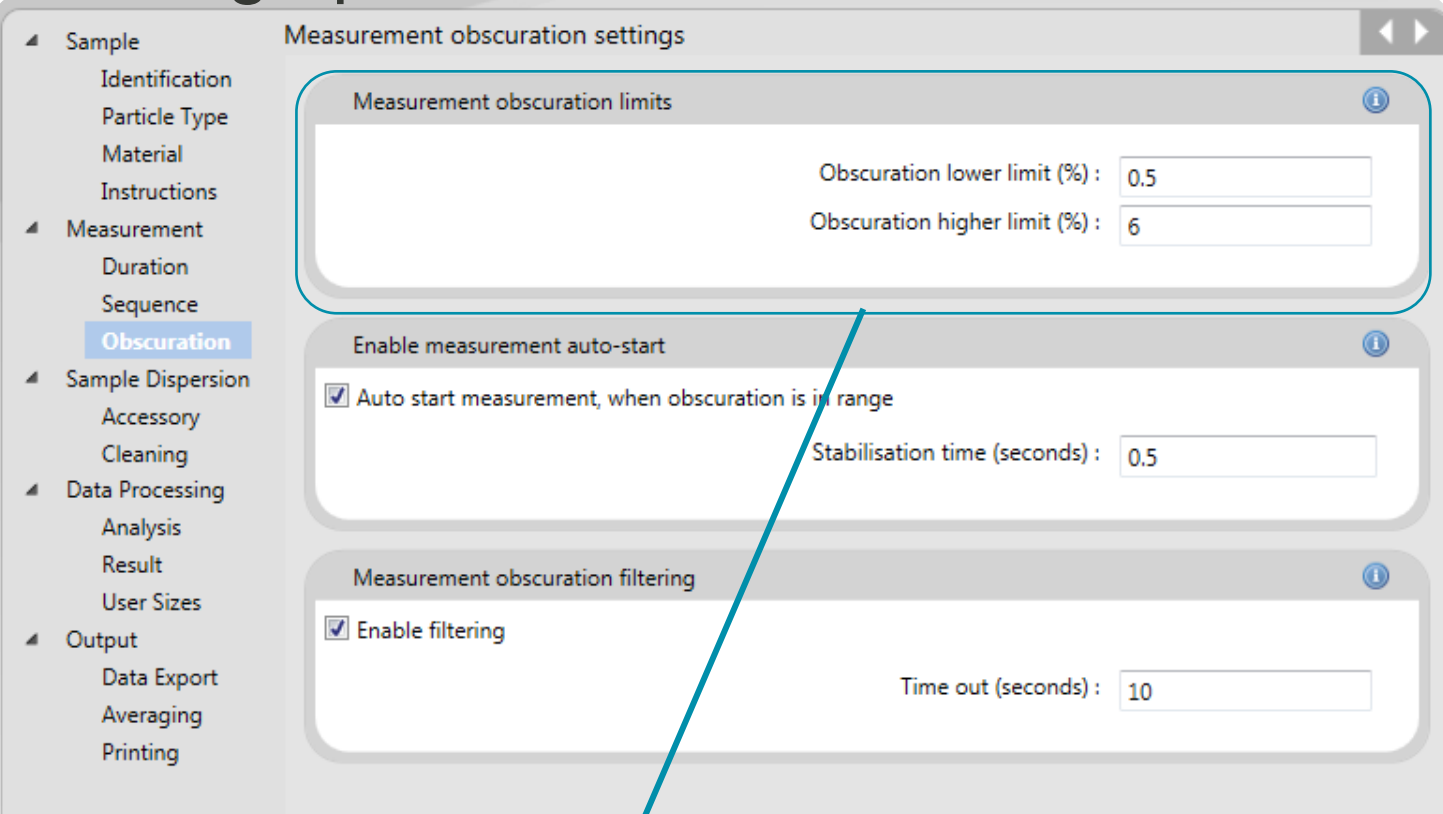
- General-purpose sample tray
 - Designed for bulk powders
 - Hopper – designed to regulate the flow flowing powders
 - Hopper height can be adjusted to control powder flow



Step 1: Setting up the feed rate

- Use a manual measurement to test the feed rate
 - Set up a long measurement
 - This gives you time to adjust the feed rate and see the response
 - Use high pressure, 4bar
 - High obscurations are likely to be observed at higher pressures
- Set up obscuration range
 - Coarse particles: ~0.5% to ~6%
 - Fine particles: ~0.5% to ~3%
- Turn on obscuration filtering
- Increase the feed rate until the obscuration is in range
 - And within range for the majority of the measurement

Step 1: Setting up the feed rate



Measurement obscuration settings

Measurement obscuration limits

Obscuration lower limit (%) : 0.5

Obscuration higher limit (%) : 6

Enable measurement auto-start

Auto start measurement, when obscuration is in range

Stabilisation time (seconds) : 0.5

Measurement obscuration filtering

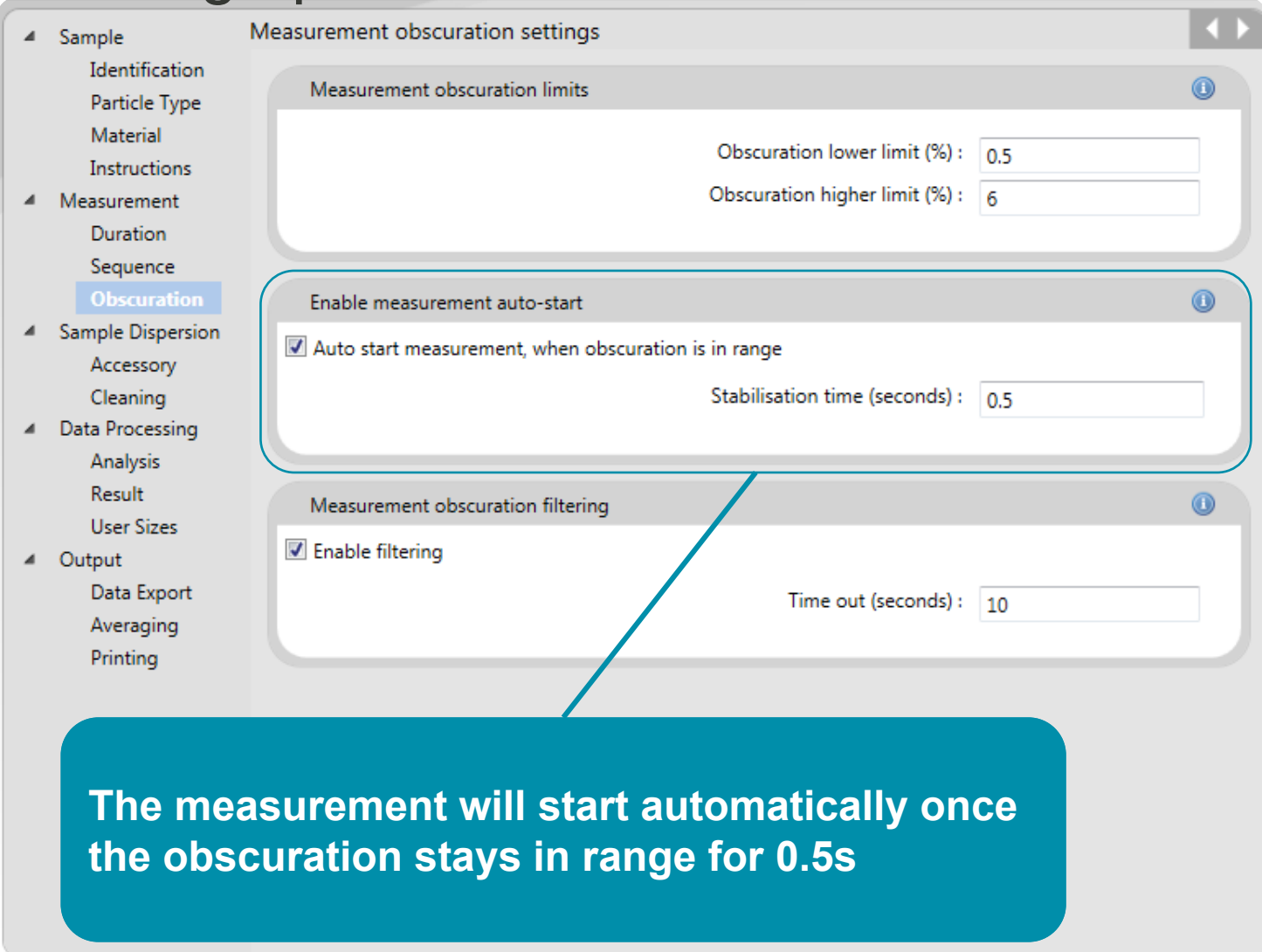
Enable filtering

Time out (seconds) : 10

Set up the obscuration range

- Coarse particles: ~0.5% to ~6%
- Fine particles: ~0.5% to ~3%

Step 1: Setting up the feed rate



Measurement obscuration settings

Measurement obscuration limits

Obscuration lower limit (%) : 0.5

Obscuration higher limit (%) : 6

Enable measurement auto-start

Auto start measurement, when obscuration is in range

Stabilisation time (seconds) : 0.5

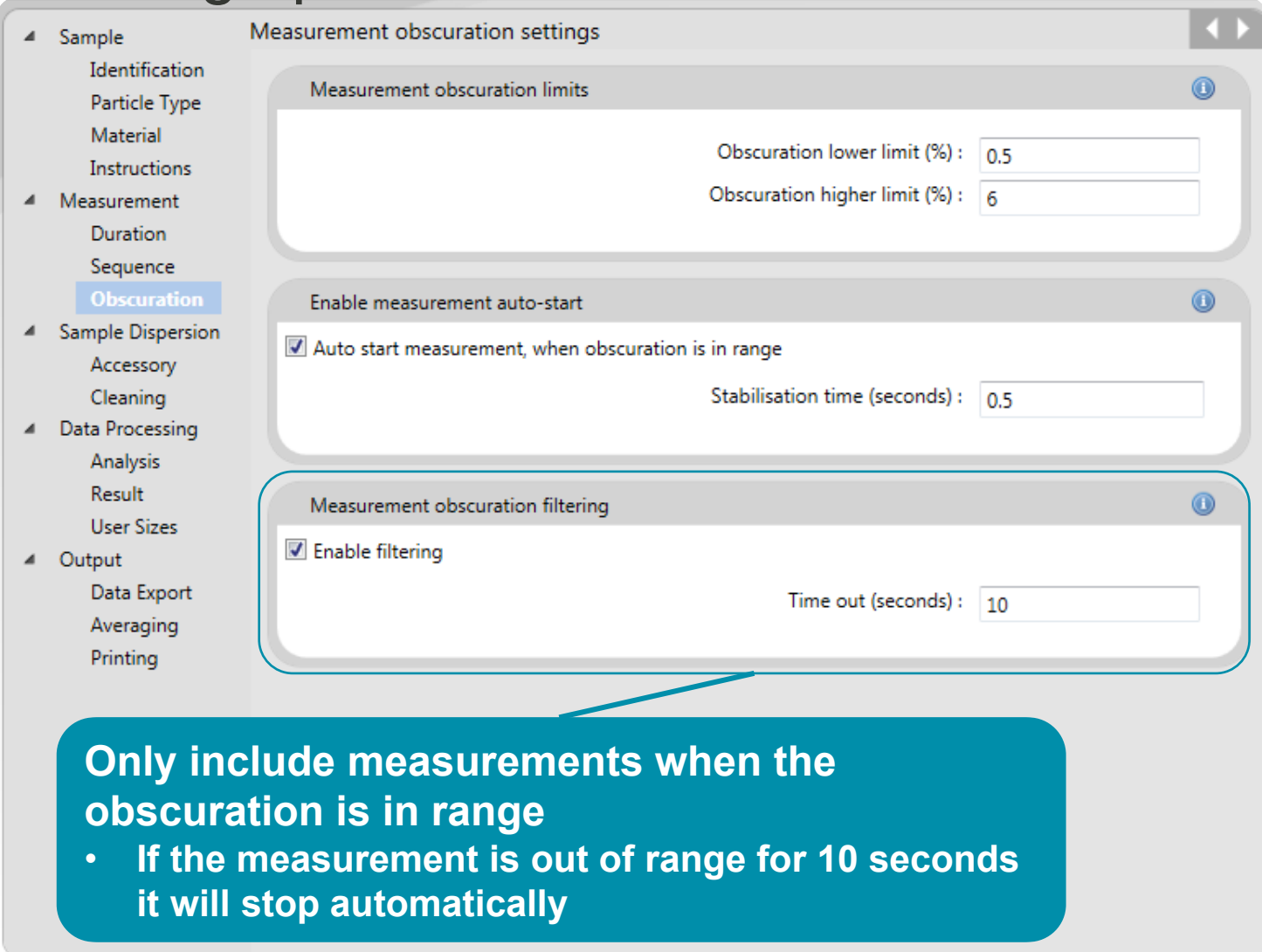
Measurement obscuration filtering

Enable filtering

Time out (seconds) : 10

The measurement will start automatically once the obscuration stays in range for 0.5s

Step 1: Setting up the feed rate



The screenshot shows the 'Measurement obscuration settings' window. The left sidebar contains a tree view with categories: Sample (Identification, Particle Type, Material, Instructions), Measurement (Duration, Sequence, **Obscuration**), Sample Dispersion (Accessory, Cleaning), Data Processing (Analysis, Result, User Sizes), and Output (Data Export, Averaging, Printing). The main area is titled 'Measurement obscuration settings' and contains three sections:

- Measurement obscuration limits:** Obscuration lower limit (%) : 0.5, Obscuration higher limit (%) : 6
- Enable measurement auto-start:** Auto start measurement, when obscuration is in range, Stabilisation time (seconds) : 0.5
- Measurement obscuration filtering:** Enable filtering, Time out (seconds) : 10

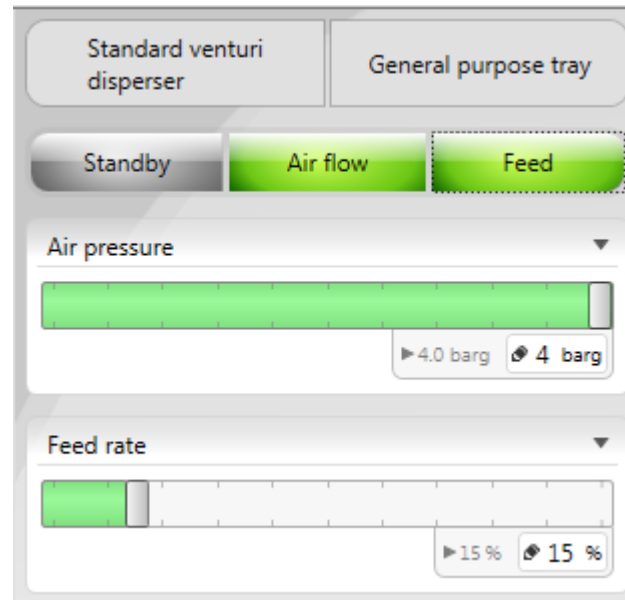
A blue callout box at the bottom contains the following text:

Only include measurements when the obscuration is in range

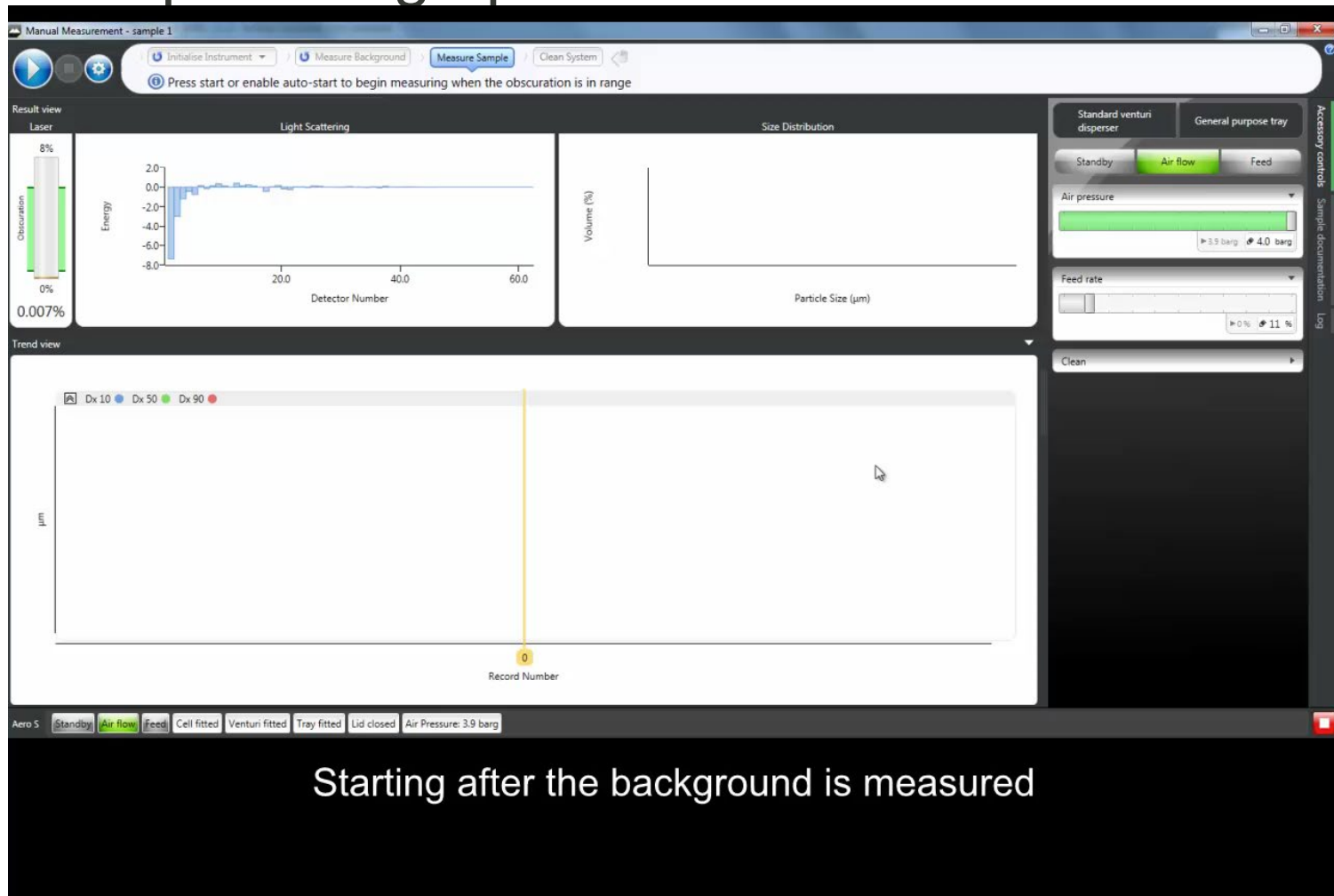
- If the measurement is out of range for 10 seconds it will stop automatically

Step 1: Setting up the feed rate

- Increase the feed rate until the obscuration is in range
 - And stays within range for the majority of the measurement



Video clip: setting up the feed rate



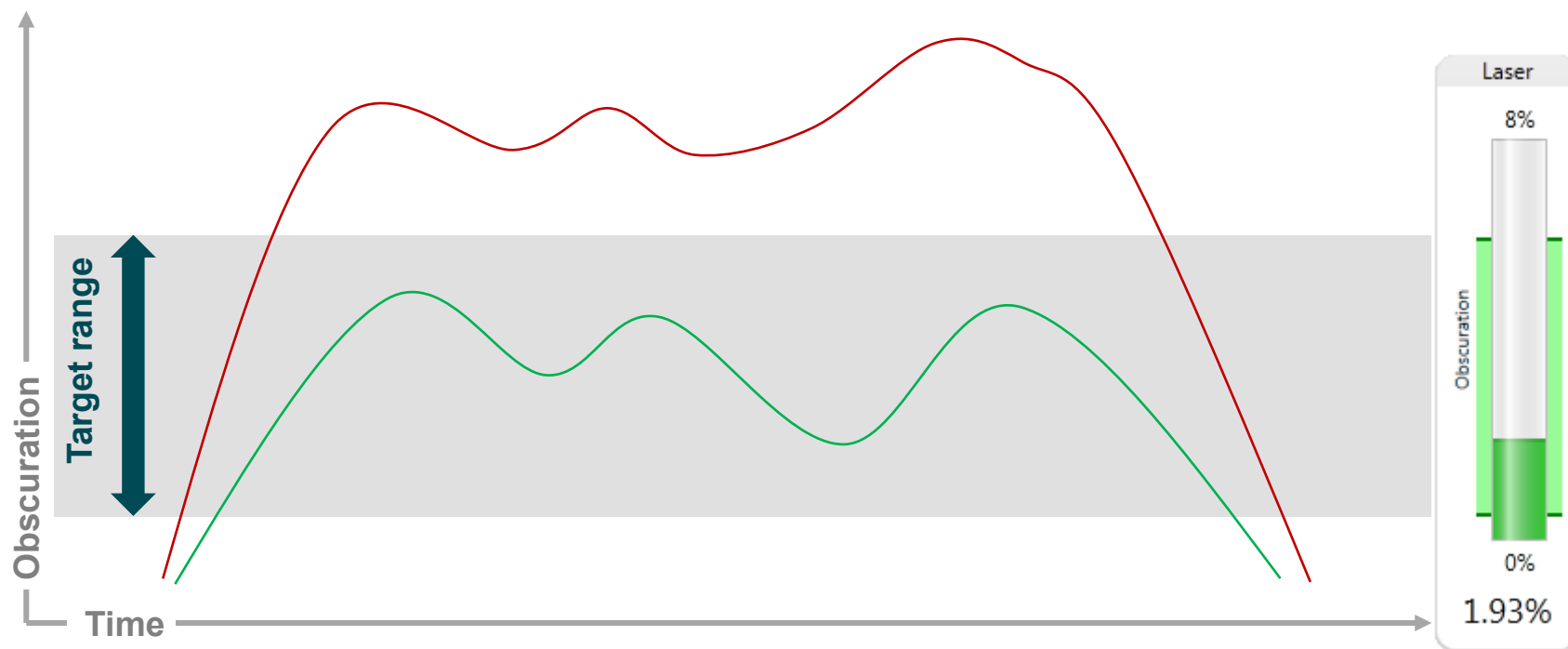
The screenshot displays the Malvern Panalytical software interface for manual measurement. The window title is "Manual Measurement - sample 1". The top navigation bar includes buttons for "Initialise Instrument", "Measure Background", "Measure Sample", and "Clean System". A status message reads: "Press start or enable auto-start to begin measuring when the obscuration is in range".

The interface is divided into several sections:

- Result view:**
 - Laser:** Shows an obscuration level of 8% (indicated by a green bar) and a target of 0.007%.
 - Light Scattering:** A plot of Energy vs. Detector Number (0 to 60.0). The energy values are mostly negative, ranging from approximately -8.0 to 2.0.
 - Size Distribution:** A plot of Volume (%) vs. Particle Size (μm).
- Trend view:** A large plot area with a y-axis labeled μm and an x-axis labeled "Record Number". A single data point is visible at record number 0.
- Control Panel (Right):**
 - Buttons for "Standby", "Air flow" (highlighted in green), and "Feed".
 - Air pressure control: Set to 3.9 barg, with a target of 4.0 barg.
 - Feed rate control: Set to 0%, with a target of 11%.
 - A "Clean" button.
- Status Bar (Bottom):** Shows system status: "Aero S", "Standby", "Air flow" (highlighted), "Feed", "Cell fitted", "Venturi fitted", "Tray fitted", "Lid closed", and "Air Pressure: 3.9 barg".


Starting after the background is measured

Feed rate is selected to keep obscuration in range



Rec 25 - dry sample 8


Obscuration is okay for this particle size.
The fit is good

Overall data quality: Pass 



Rec 24 - dry sample 8

Obscuration is okay for this particle size.
The fit is good
Poor obscuration control, check feed'

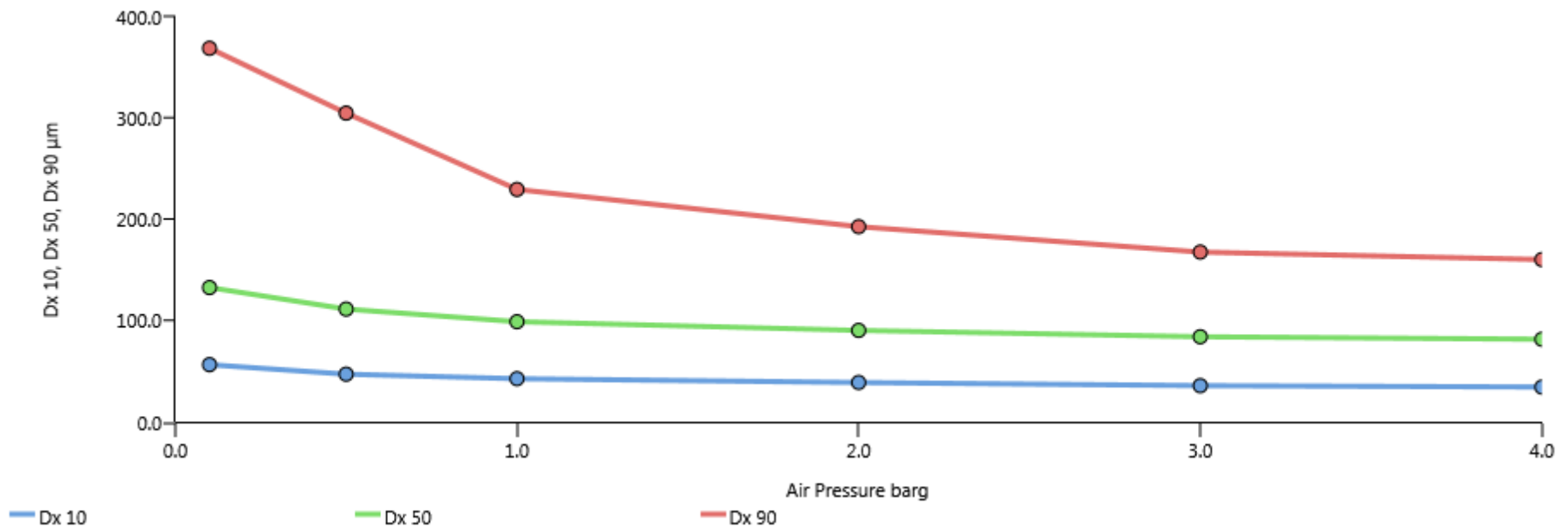
Overall data quality: Fail 

Dry powder dispersion: ISO guidance

- Degree of dispersion is controlled by primary air pressure
 - Monitor change in size distribution with pressure
 - Carry out pressure titration – **Step 2**
- Check that particle comminution (milling or particle break up) has not occurred
 - Compare dry results to a well dispersed wet measurement – **Step 3**
 - Choose the pressure which agrees with the wet results
 - Shows dispersion and not particle breakage

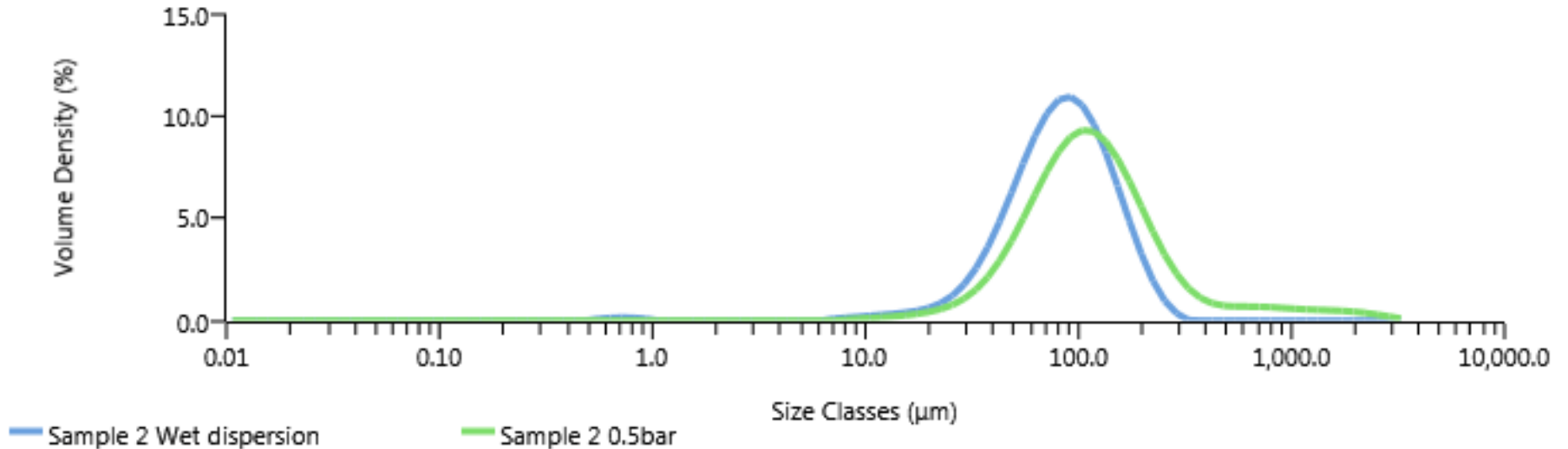
Step 2: Measure a pressure titration

- Make measurements at 4, 3, 2, 1, 0.5 and 0.1 bar.
- Make repeat measurements at each pressure to check for sample segregation



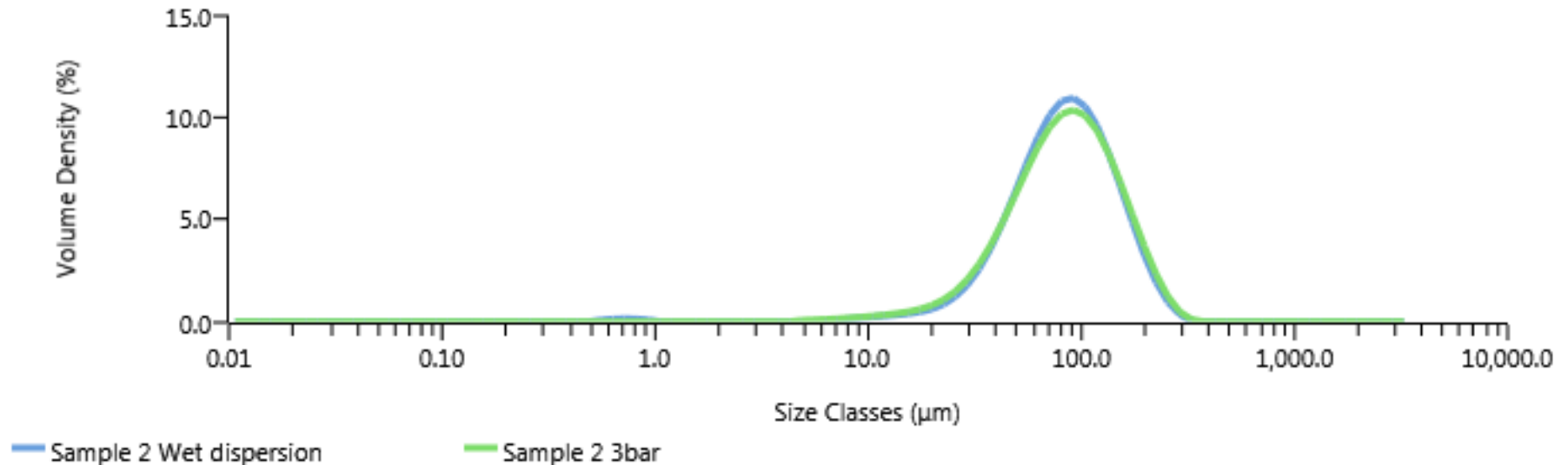
Step 3: Compare dry results to wet

- Low pressure dry result shows larger result
 - Indicates sample is not fully dispersed



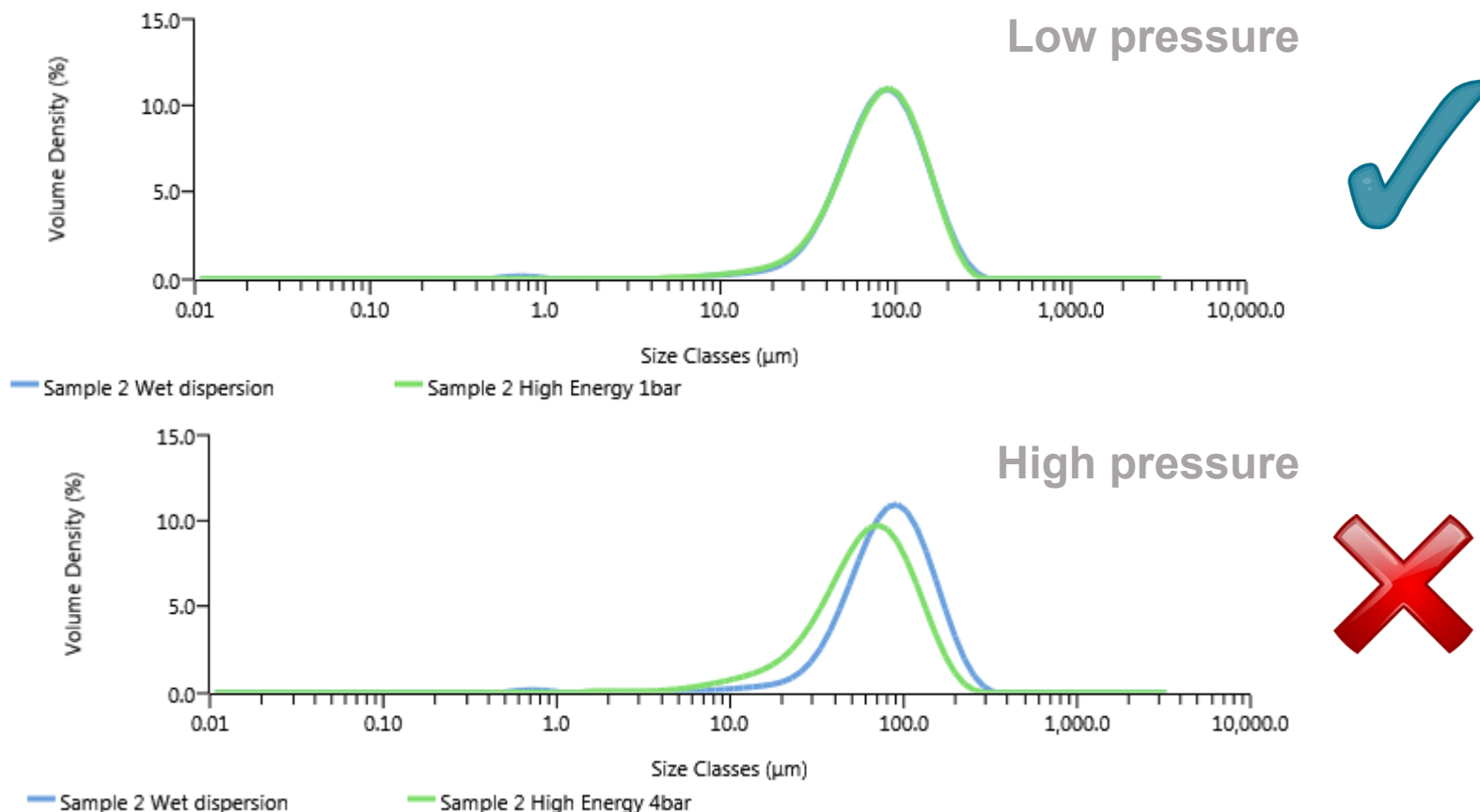
Step 3: Compare dry results to wet

- High pressure shows good agreement
 - Suggests the material is dispersed

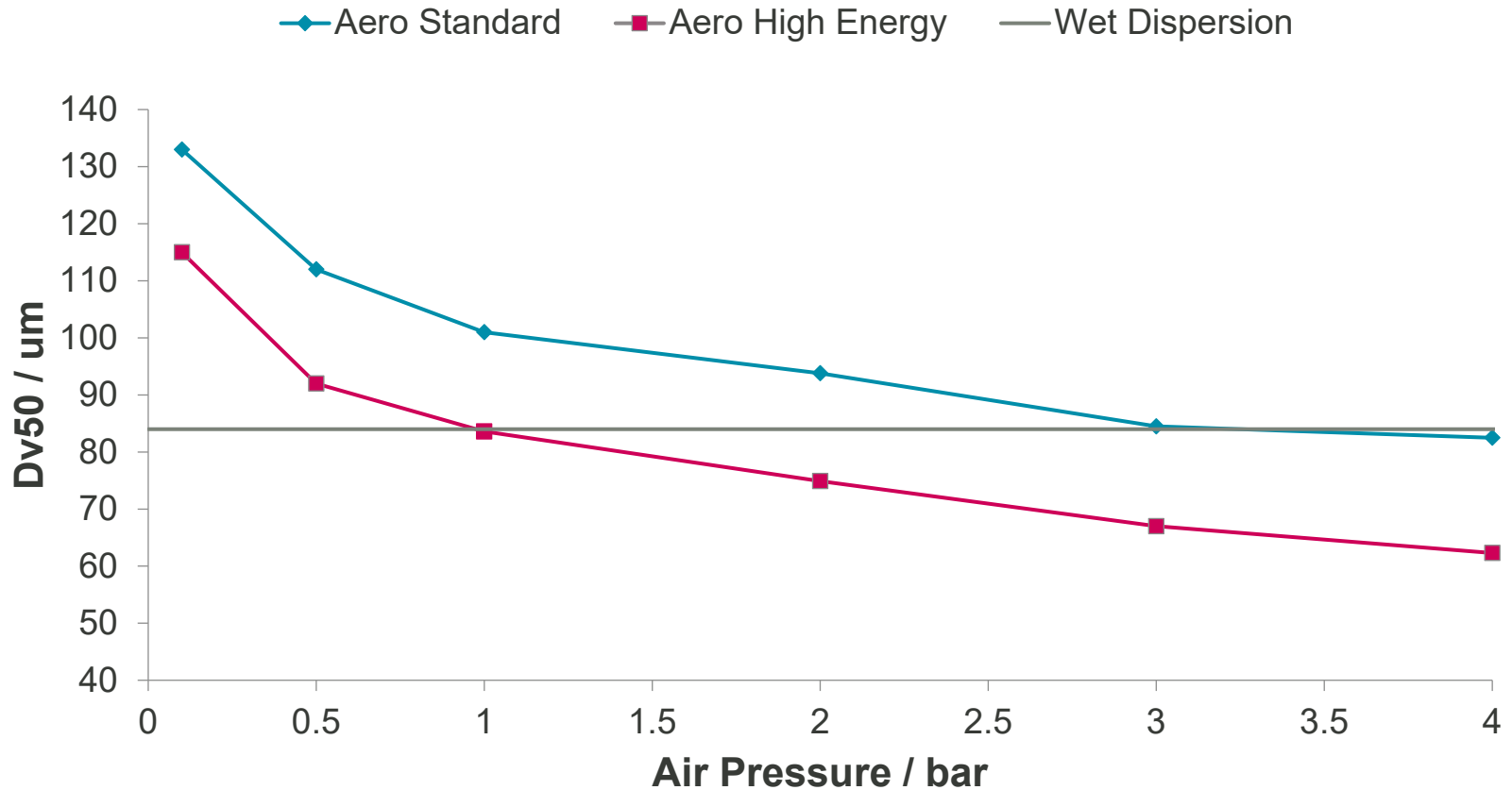


Step 4 : High energy venturi

- For robust, highly agglomerated materials the high energy venturi may be required.

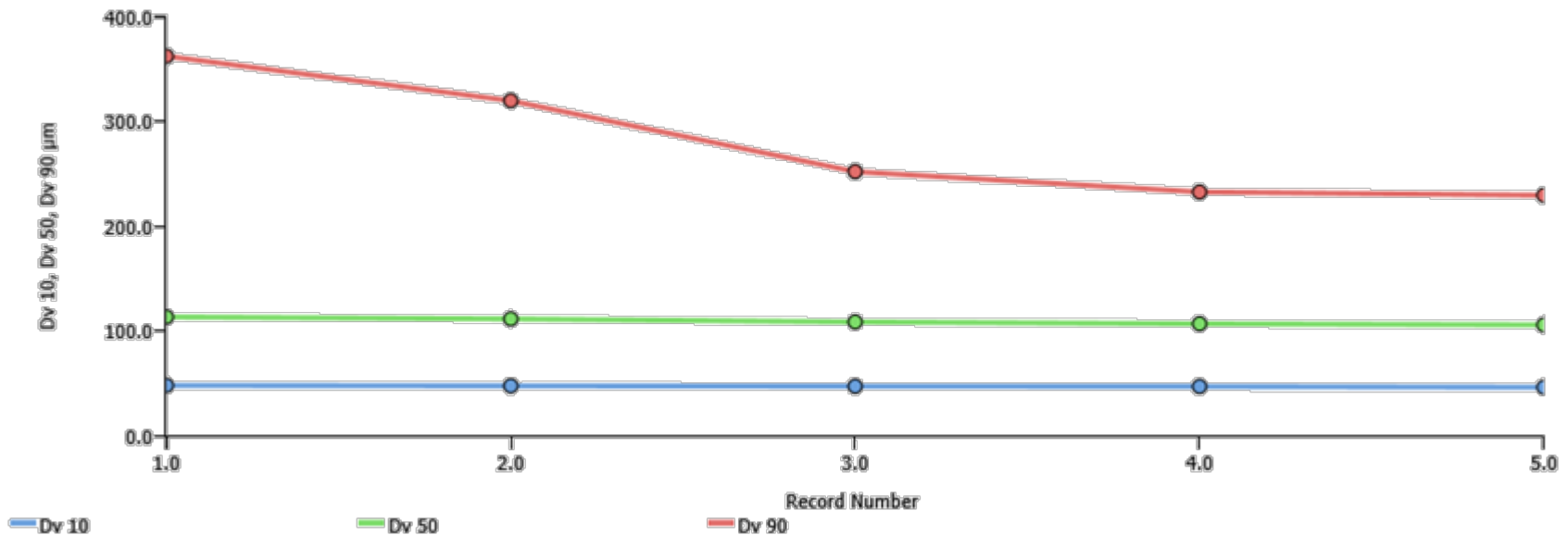


Comparing standard and high energy venturitis



Segregation in dry measurements

- Segregation can occur with free-flowing powders with wide particle size distributions
 - Characterized by a decrease in size over repeat measurements
 - Make several quick repeat measurements at each pressure
 - This can be done as part of the pressure titration



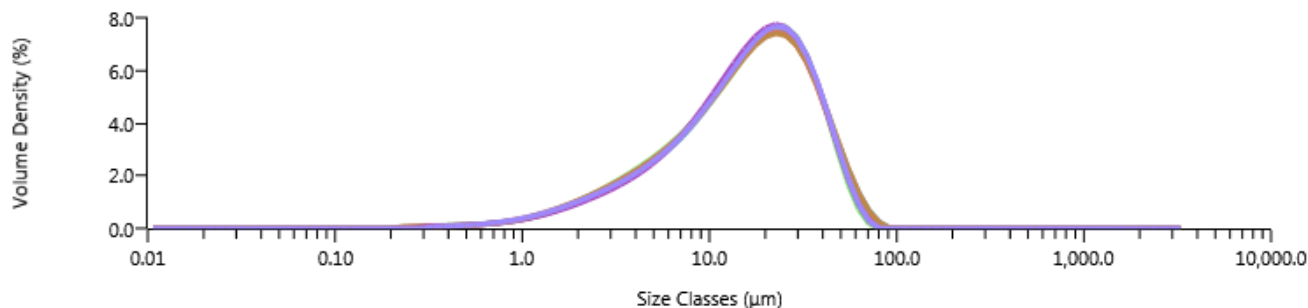
Segregation in dry measurements

- Segregation can occur with free-flowing powders with wide particle size distributions
 - Characterized by a decrease in size over repeat measurements
 - Make several quick repeat measurements at each pressure
 - This can be done as part of the pressure titration
- Always measure the whole sample, either by:
 - Making enough short measurements to use the whole sample and then create an average
 - Make one long measurement long enough to use up all of the sample.

Definitions of repeatability, reproducibility etc.

- Repeatability
 - Same sample run time and time again
 - Critical Factors:
 - Instrument
 - Dispersion i.e. any changes occurring during a measurement
- Reproducibility
 - Another sample, operator, instrument, technique, day, etc.....
 - Critical Factors:
 - Instrument
 - Dispersion
 - Sampling methodology
- Robustness
 - Capacity of method to remain unaffected by small, deliberate, variations in method parameters

Reproducibility: Sampling



| | Record Number | Sample Name | Dx 10 (µm) | Dx 50 (µm) | Dx 90 (µm) |
|------------------|---------------|------------------|--------------|--------------|--------------|
| | 1 | Baby Powder 50mg | 3.62 | 16.6 | 40.4 |
| | 2 | Baby Powder 40mg | 3.47 | 16.2 | 38.1 |
| | 3 | Baby Powder 30mg | 3.51 | 16.2 | 39.4 |
| | 4 | Baby Powder 20mg | 3.79 | 16.5 | 38.8 |
| | 5 | Baby Powder 10mg | 3.57 | 16.5 | 40.4 |
| | 6 | Baby Powder 5mg | 3.61 | 16.5 | 38.9 |
| Mean | | | 3.59 | 16.4 | 39.3 |
| 1xStd Dev | | | 0.112 | 0.180 | 0.917 |
| 1xRSD (%) | | | 3.12 | 1.10 | 2.33 |

- Reproducible results can be achieved for very small sample masses.

General rules for good dry measurements

- Ensure that the obscuration filtering is on
 - Eg 0.5 to 6%
 - Set the feed rate to keep the obscuration in range
- Always measure the whole sample
 - In several short measurements, which are averaged
 - Or one long measurement
- Measure several sub samples from the same batch to assess reproducibility

The purpose of method development

- A laser diffraction measurement requires

‘a representative sample, dispersed at an adequate concentration in a suitable liquid or gas’

<USP429>

- Method development must define appropriate
 - Sampling
 - Dispersion
 - Measurement conditions