

Characterisation for powdered- polymer Additive Manufacturing

Dr Candice Majewski (she/her)

The University of Sheffield

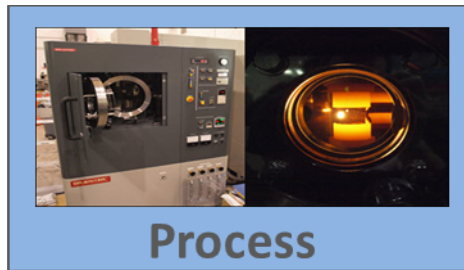
c.majewski@sheffield.ac.uk

MAPP Hub

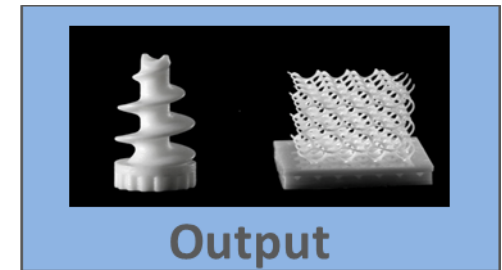
- **Manufacture using Advanced Powder Processes**
 - £20 million research hub, (£10 million funding from EPSRC, £7 million industrial support, over £3 million from collaborating universities)
 - Overall aim to enable Advanced Powder Processes to live up to their potential...
 - See <https://mapp.ac.uk/> for details



Designed for process



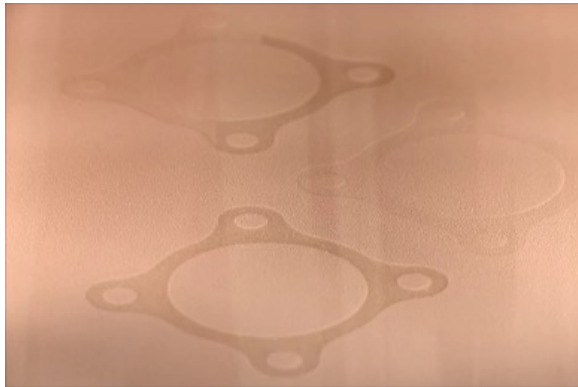
Monitored
Dynamic control via machine
learning



Designed
Quality built in

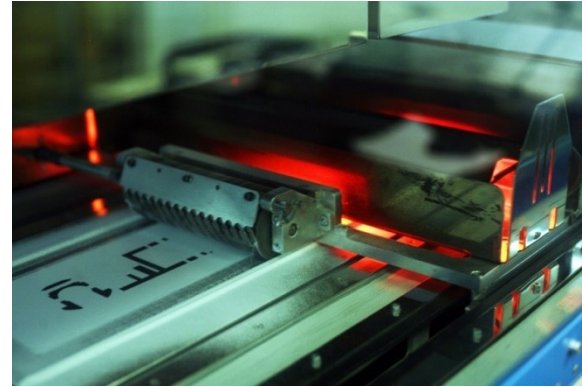
Advanced Polymer Sintering laboratory

- Major focus on powdered polymer Additive Manufacturing
 - Laser Sintering, High Speed Sintering & others
 - Process and material developments
 - Understanding of interaction between process and material



Laser Sintering

Parts built by selectively scanning and sintering cross-sections of powdered material

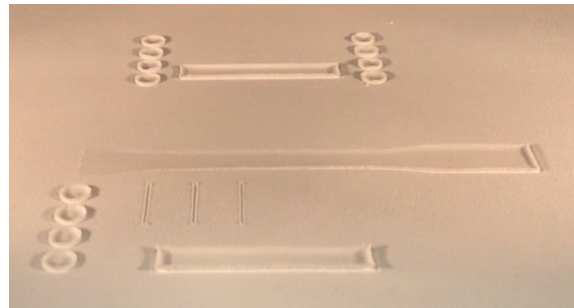
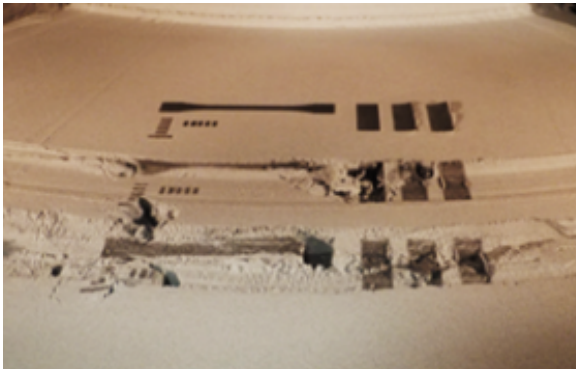


High Speed Sintering

Required cross-section ink-jet printed with a Radiation Absorbing Material, then sintered using an infra-red lamp

Interaction between material and process

- We want to know how our powders and processes interact, and what effect do these interactions have on the resultant parts?
 - Need to process a wider range of materials, more efficiently, and more repeatably...
 - We can't do this until we understand exactly what's going on!



Where do characterisation techniques come in?*

*Spoiler alert... Pretty much everywhere!

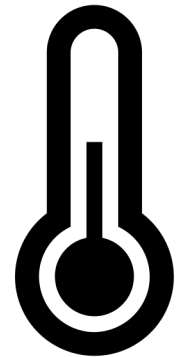
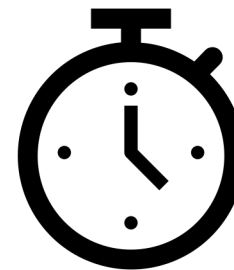
Powder deposition

- A crucial stage... without a smooth, well-packed powder base to start with, we can end up with parts which have:
 - Low density
 - Poor accuracy/surface finish
 - Over-exposure of certain areas of parts (non-homogenous parts)
- We might want to characterise:
 - Particle size and size distribution
 - Particle shape
 - Powder flow
 - Moisture content
 - Static
 - Etc



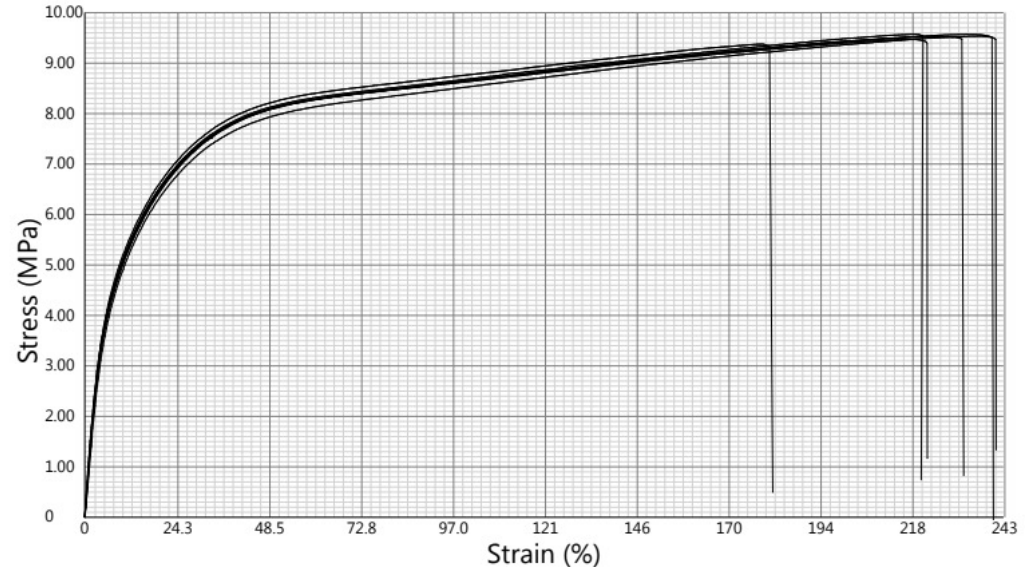
Parameter setting

- How can we determine the ‘best’ parameters to process a given material?
 - How much energy does it take to melt the material?
 - How quickly will particles fuse together?
 - What pre-heat temperatures should we use?
 - At what point will we start to damage or degrade our material?
 - What sort of cooling regime do we need?



Analysing part structure and quality

- What's going on in our parts?* For example:
 - Mechanical properties
 - Microstructure
 - Porosity
 - Flame retardance
 - Fatigue life
 - Etc., etc.



* (Very dependent on intended application)

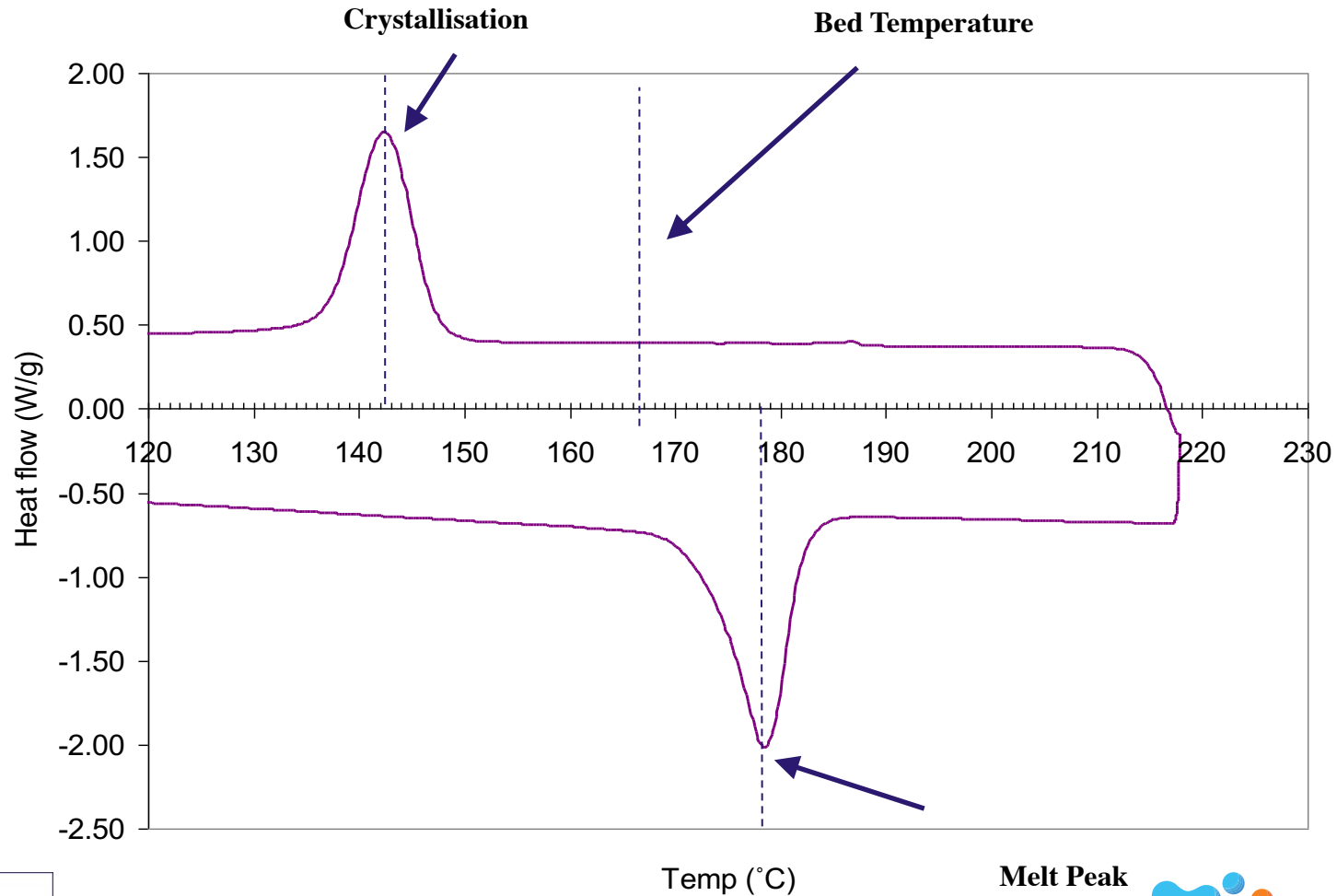
A couple of specific examples

(Please ask about others if you're interested)

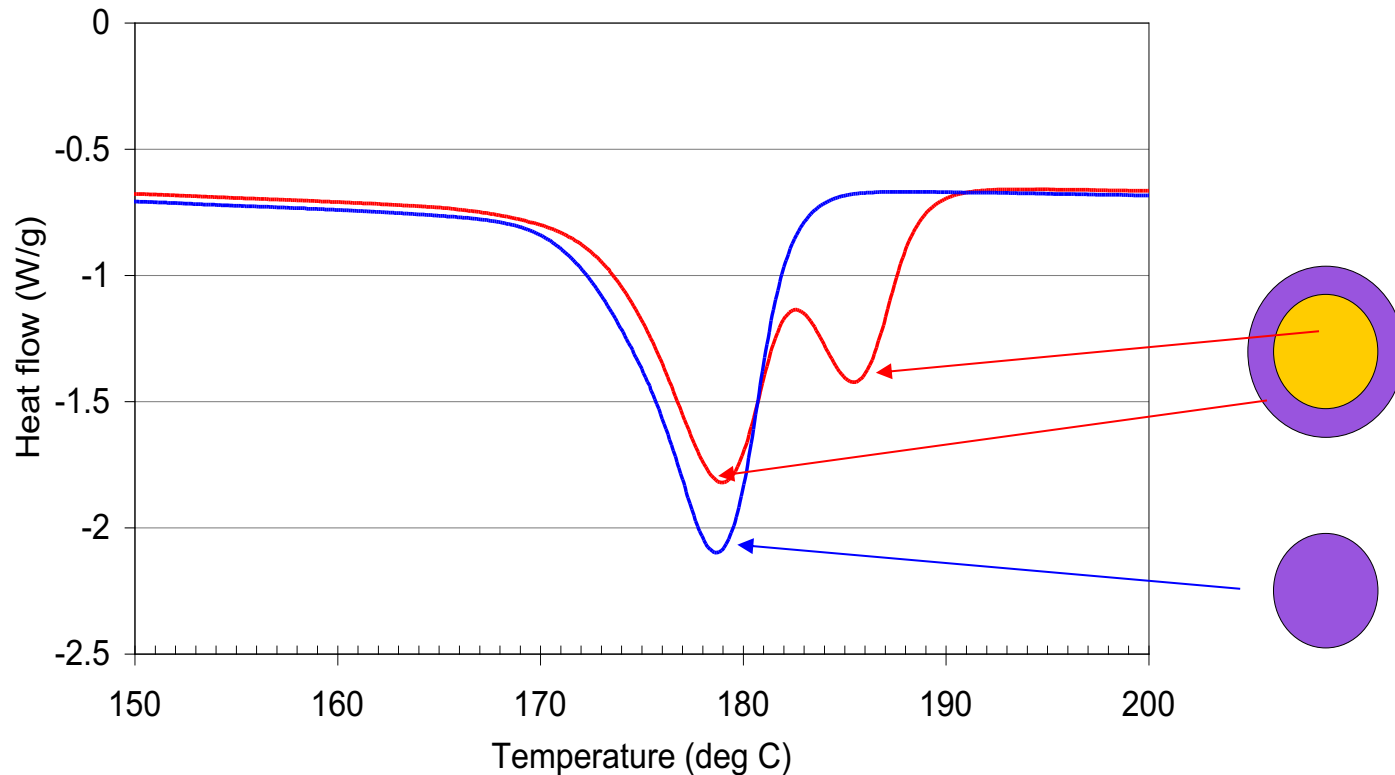
Differential Scanning Calorimetry

- Thermal analysis technique
 - Measure difference in heat flow between a sample and a known reference, with respect to time and temperature
 - For example, if we heat this reference sample at a certain rate, how much energy do we need to put into our test sample to match this?
- Useful for...

Differential Scanning Calorimetry – choosing sintering parameters

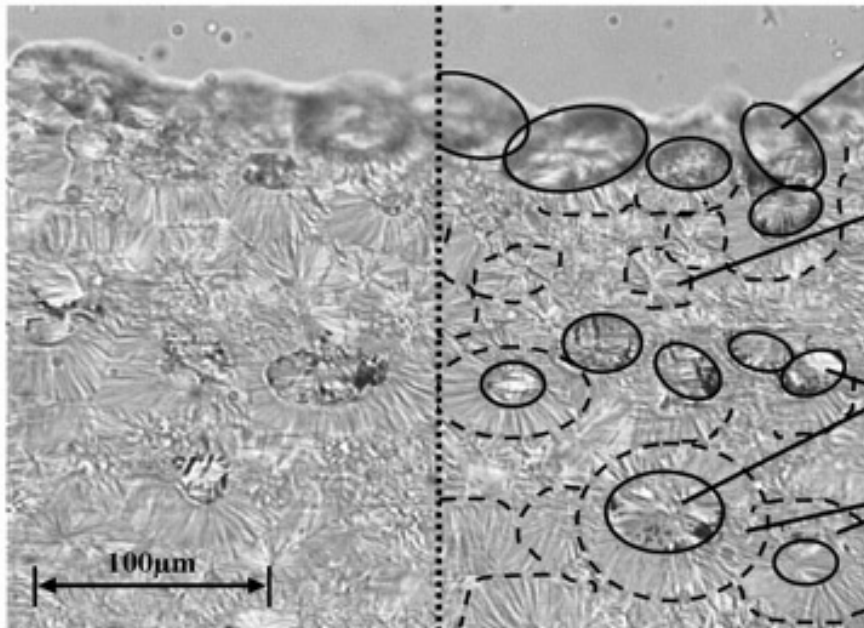


Differential Scanning Calorimetry – microstructure of parts



Differential Scanning Calorimetry – microstructure of parts

- What this shows us...
 - Parts are often comprised of regions with varying proportions of melting.



Unmolten particle fused to edge

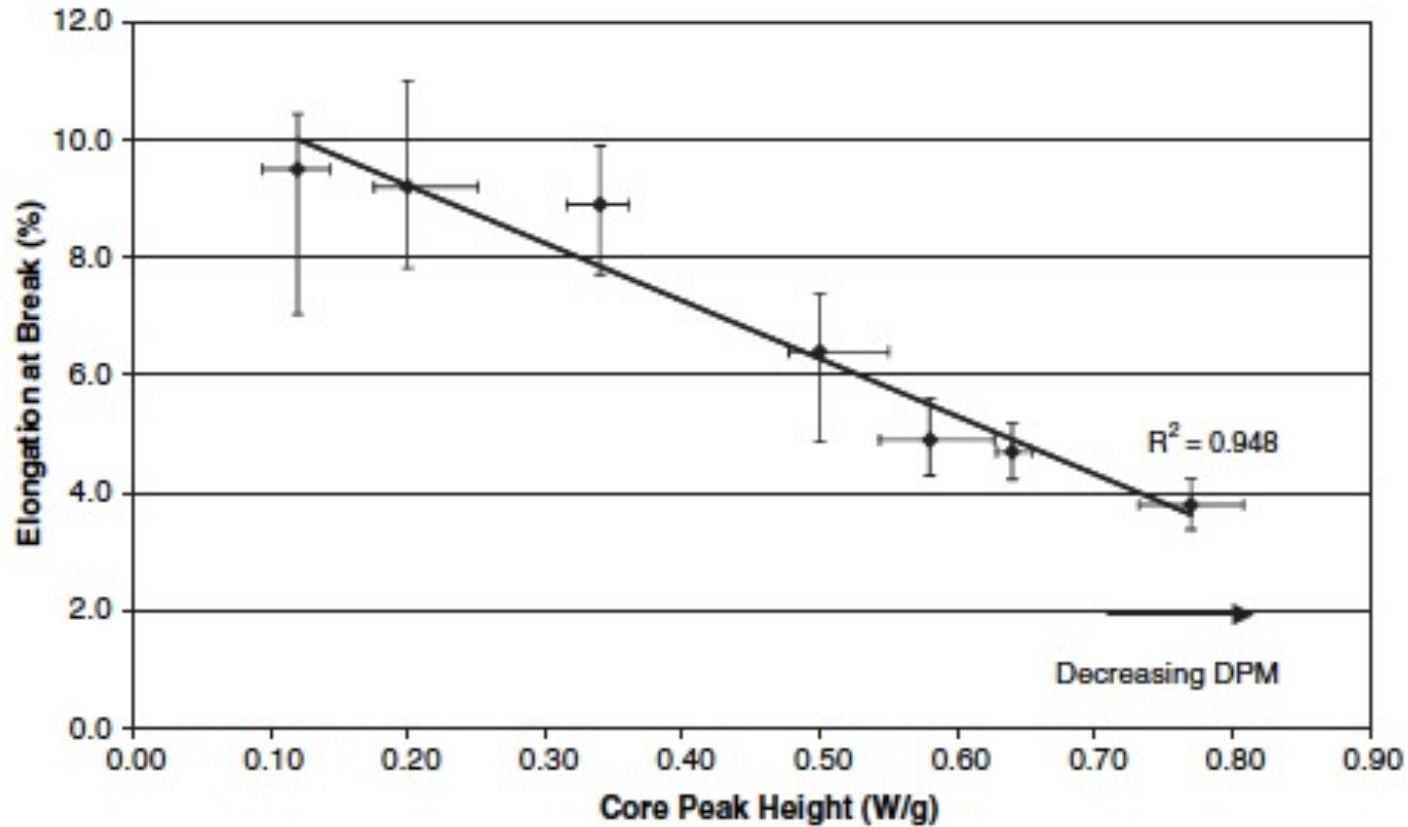
Spherulite from fully melted & crystallised particle

Unmolten particle
core

Spherulite from melted & crystallised
region

Zarringharam, H., Hopkinson, N., Kamperman, N.F., de Vlieger, J.J., 2006, Effects of processing on microstructure and properties of SLS Nylon 12, Materials Science and Engineering A, Vol. 435-436, pp 172-180, ISSN 0921-5093

Differential Scanning Calorimetry – microstructure of parts



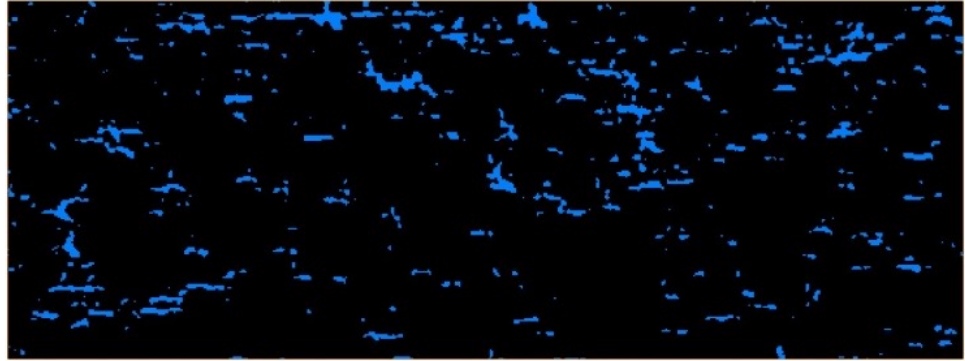
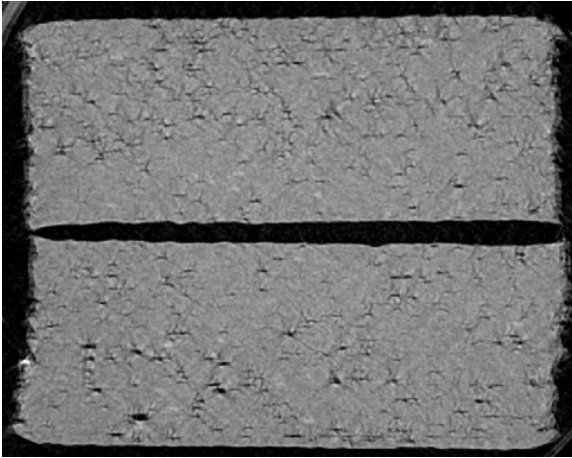
Majewski C, Zarringhalam H & Hopkinson N (2008) Effect of the degree of particle melt on mechanical properties in selective laser-sintered Nylon-12 parts.

X-ray computer micro-tomography (Micro-CT)

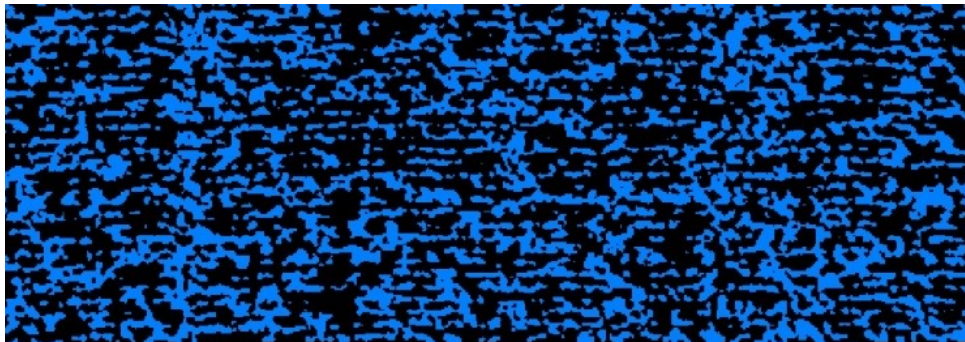
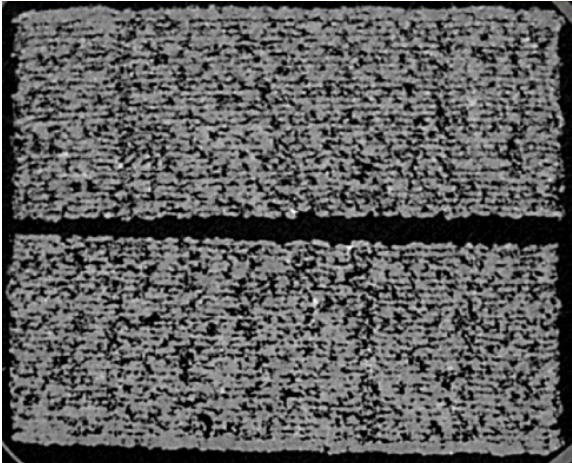
- A non-destructive 3D imaging technique
 - Can help in analysing the internal and external geometry of complex components

- Useful for...

Micro-CT – understanding porosity

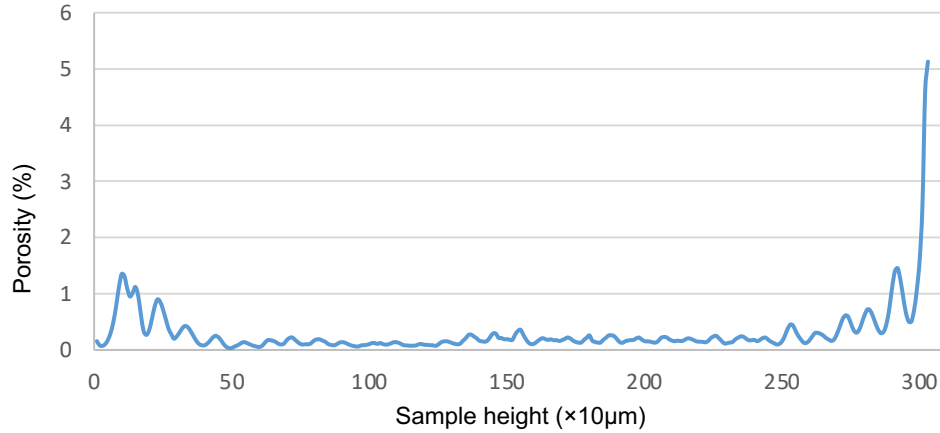


High energy - **Porosity: 6.79% (vol)**

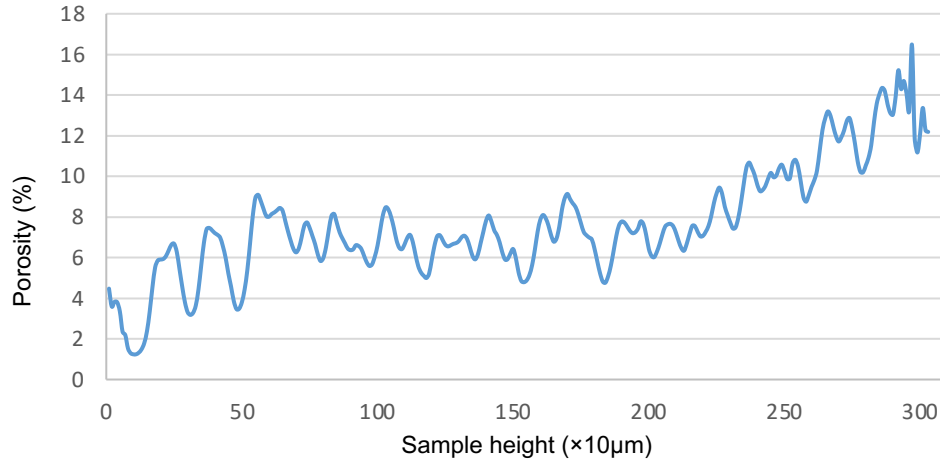


Low energy - **Porosity: 32.08% (vol)**

Micro-CT – understanding porosity

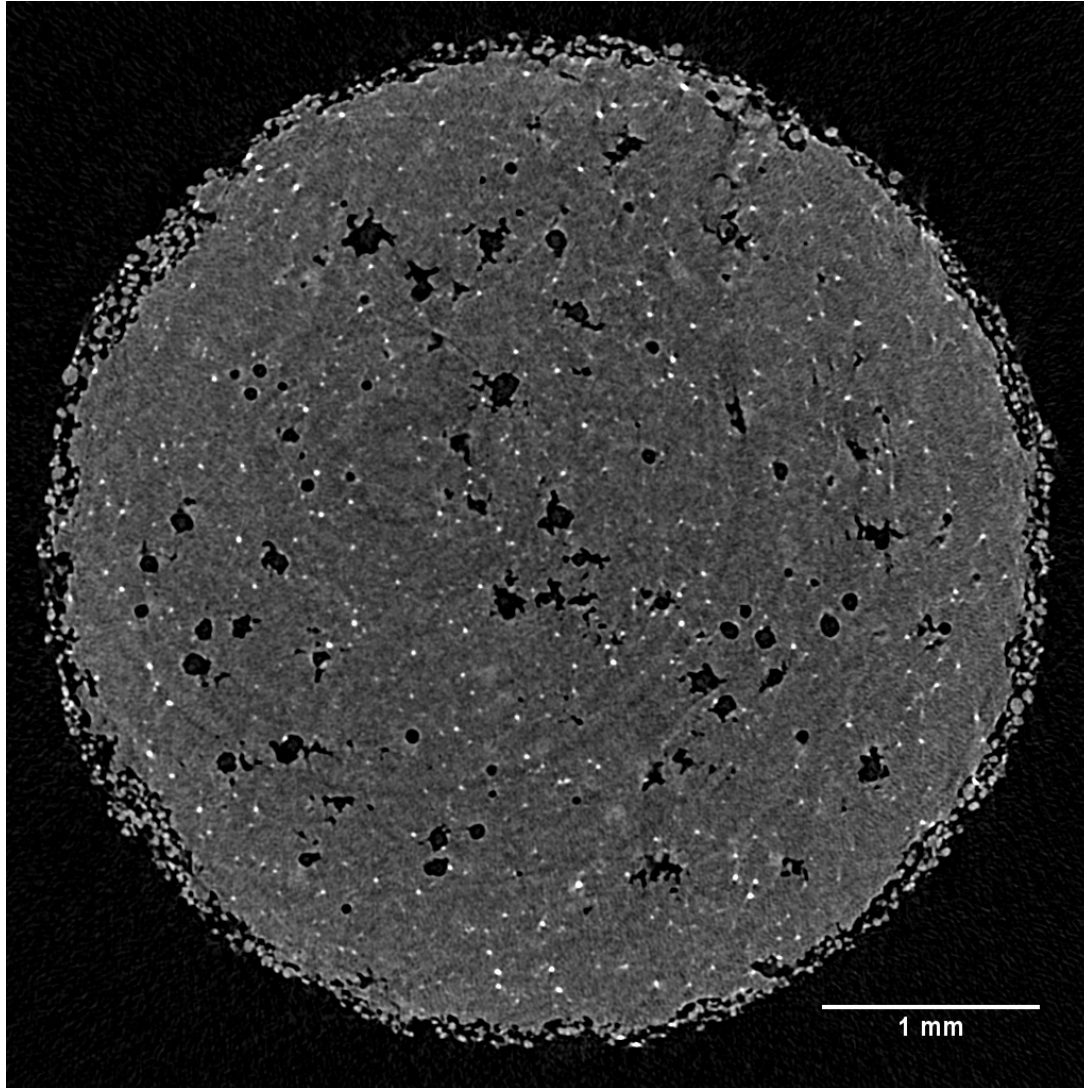


Higher energy - Porosity: 0.58% (vol)

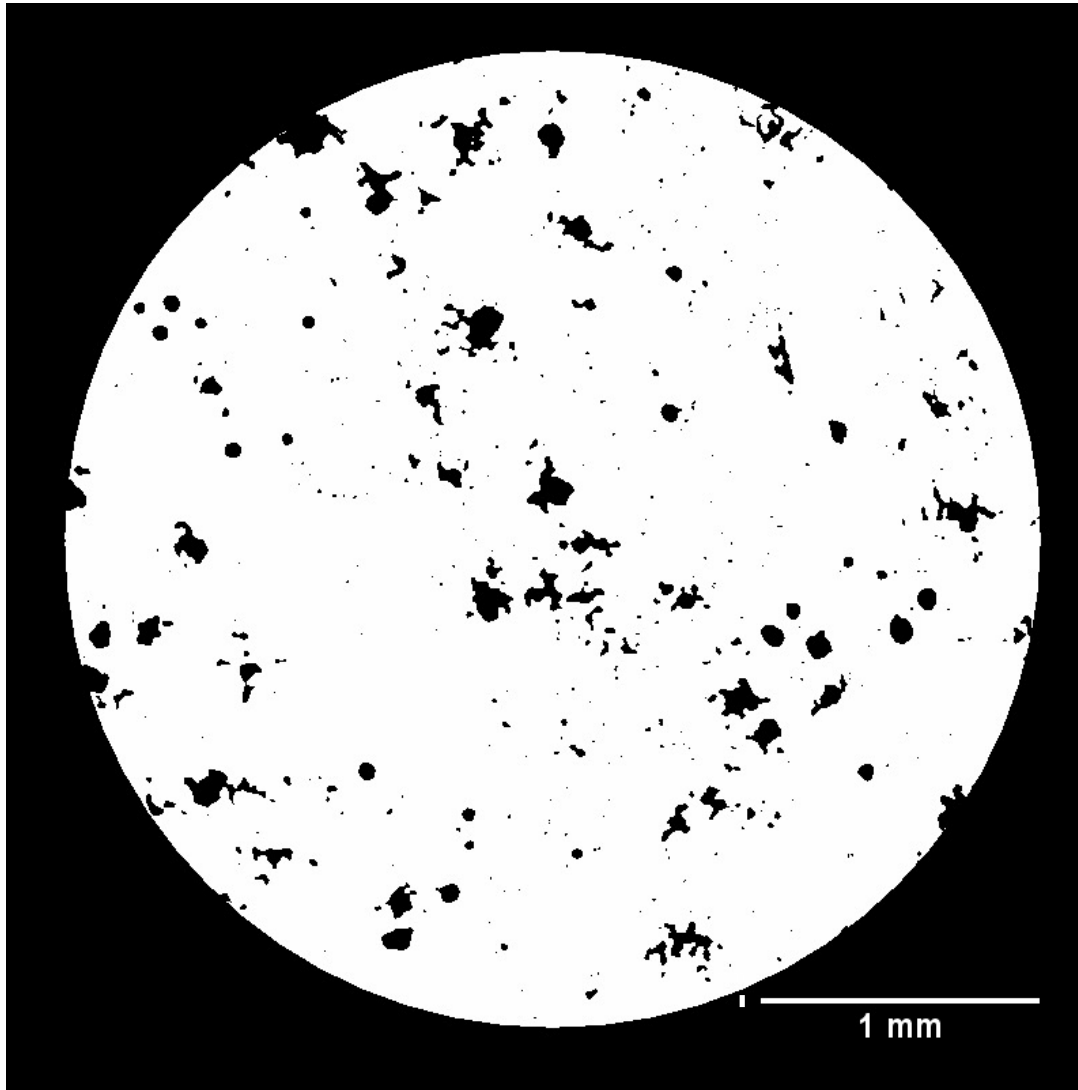


Lower energy - Porosity: 7.36% (vol)

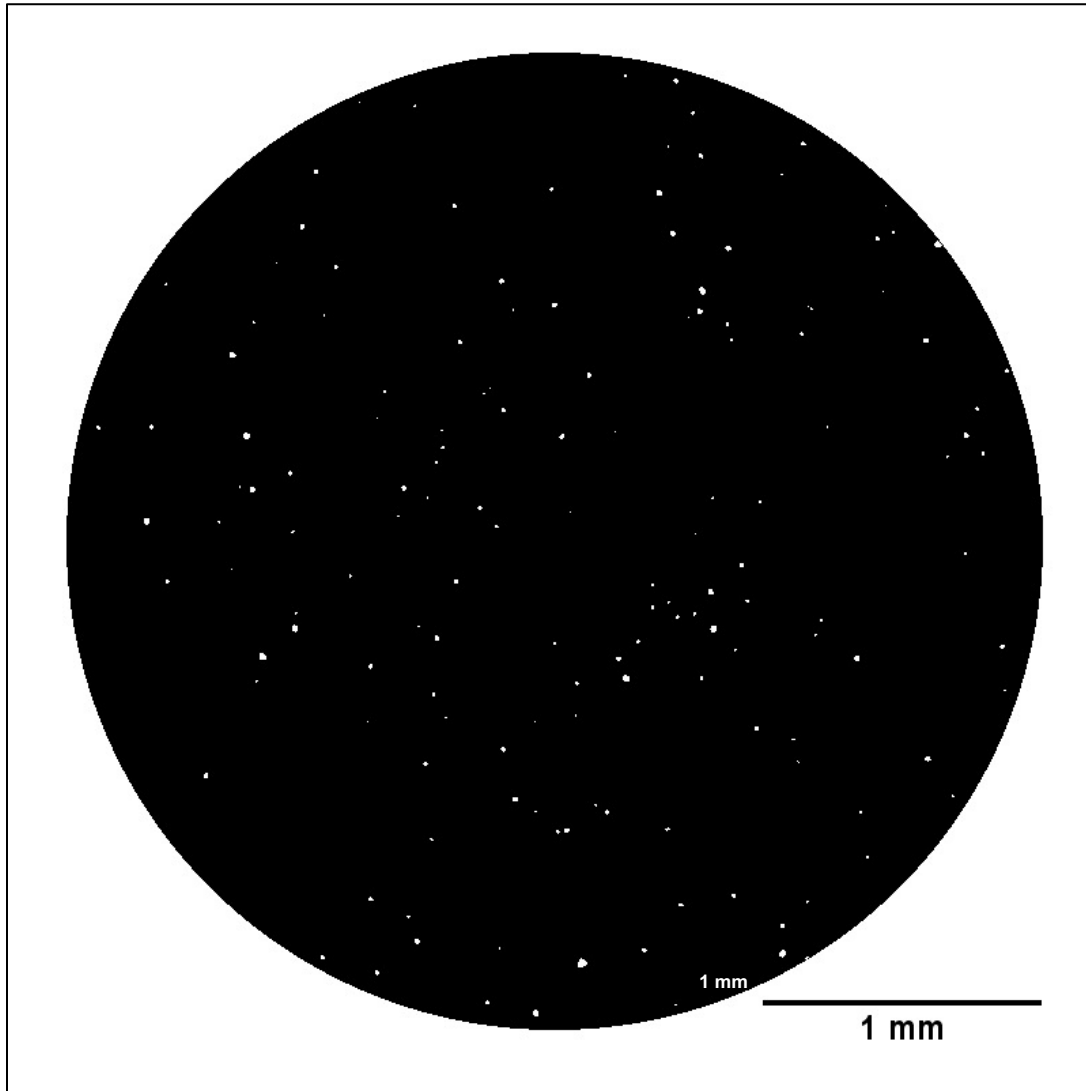
Micro-CT – homogeneity



Micro-CT – homogeneity



Micro-CT – homogeneity



Any Questions?

Dr Candice Majewski (she/her)
The University of Sheffield
c.majewski@sheffield.ac.uk