

# SOFTWARE UPDATE NOTIFICATION (SUN) – HYDRO SIGHT SELF-DEMO SOFTWARE:PSS0009-02



**PARTICLE SIZE**



**PARTICLE SHAPE**

## Introduction

This document details the release of software PSS0009-02: the Hydro Sight Self-Demonstration software package v1.00. Installation instructions and PC requirements are provided.

## Installation

It is assumed that you have authority to install or update software within your facility. It is also assumed that you have Administrator rights for the system upon which the software is installed, as this is a requirement of the installation process. If you do not have this authority please consult with your I.T. support department before proceeding.

## Recommended System Requirements

The minimum requirements for running this software are highlighted in table 1 below. Although the software can run using Windows 8™ (64-bit), it has been tested under Windows 7™ (32-bit and 64-bit). Windows 7 is therefore the preferred operating system.

*Note:* Some 3rd party software or OS patches may prevent the Hydro Sight software from running correctly. It is not possible to test for compatibility with all windows programs.

## Supported Languages

The Hydro Sight software suite is available as an English language application.

Table 1: Recommended system requirements for the Hydro Sight software.

Feature	Specification
Processor Type	Intel Core i7 3770 Processor,
Memory	4GB
Hard Disk Storage	1Tb HDD,
Additional Storage Media	DVD +/-R/RW drive
Display	1 x 22" Widescreen Flat Panel Monitor for software
Connectivity	None
Operating System	Windows 7™ (32 bit and 64 bit), Windows 8™ (64-bit)

## Installation Instructions

The software is supplied as a self-executing installer. Once downloaded, the installer must be run to set up the software.

It is required that the sample videos packaged with the software are copied to the PC for use in the demonstration. Copy the three pre-recorded sample dispersion videos to the folder called 'Sample Videos' located in here C:\Users\Public\Videos, if necessary create a folder of this name.

## Uninstall Procedure

The software can be uninstalled using the standard Add/Remove feature in the Windows 'Control panel'.

## Known Issues

The following software bugs have been discovered within the software, and will be investigated as part of a future release. Please follow the suggested work-around for

Issue	Work around	Comment
Can only select sample video on application startup	Restart application to select new sample video	

### Malvern Instruments Ltd

Enigma Business Park • Grovewood Road  
Malvern • Worcestershire • UK • WR14 1XZ

Tel: +44 (0)1684 892456  
Fax: +44 (0)1684 892789

### Malvern Instruments Worldwide

Sales and service centers in over 50 countries  
for details visit [www.malvern.com/contact](http://www.malvern.com/contact)

© Malvern Instruments Ltd 2014

more information at [www.malvern.com](http://www.malvern.com)