

OMNISEC SOFTWARE: v10.40 (PSS0051-06) SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the installation and known issues with OMNISEC v10.40. This release is a major release to introduce new features, primarily Windows 10 support and deletion of data. In addition some issues are fixed. Installation instructions and PC requirements are provided.

Installation

It is assumed that you have authority to install or update software within your company's SOPs. If you do not have this authority, please consult with your I.T. support department before proceeding.

It is assumed that you have Administrator rights for the computer. This is required by the installation process.

Minimum System Requirements

Intel Core i7 processor, 4 GB RAM, 500 GB free hard disk space, 1600 x 900 screen resolution running in 16 bit color mode, CD-ROM drive, 1 free USB port, Windows 7 or 10 64bit Professional operating system.

Recommended System Requirements

Intel Core i7 processor (Quad Core HT, 3.1 GHz Turbo, with HD Graphics), 8 GB RAM, 500 GB free hard disk space, 1920 x 1080 full HD screen resolution running in 16 bit color mode, CD-ROM drive, 1 free USB port, Windows 7 or 10 64 bit Professional operating system.

Supported operating systems

OMNISEC software v10.40 is compatible with Windows 7 or 10. Only 64 bit Professional versions are supported.

Windows XP, Windows Vista, and Windows 8 are not supported.

Supported Languages

English

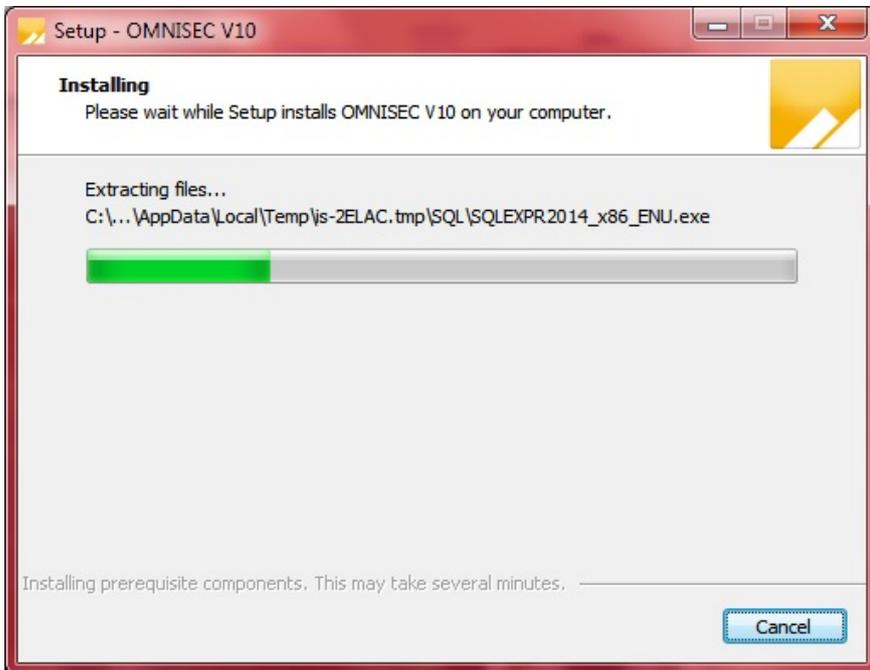
Installation Instructions

Windows 7 or 10 64 bit Professional

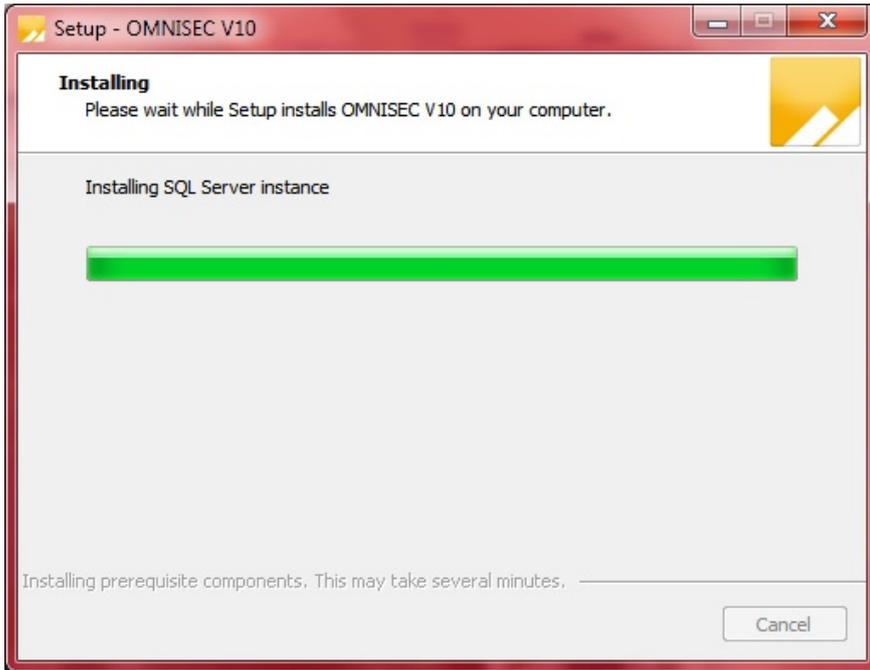
In most cases, OMNISEC should install automatically when the CD is inserted in the drive. If it doesn't, the user can double-click on the installer file, OMNISEC1040.exe, to begin the installation.



When the Setup window appears, select Next to process. You can select cancel if you would like to stop the setup and continue another time.



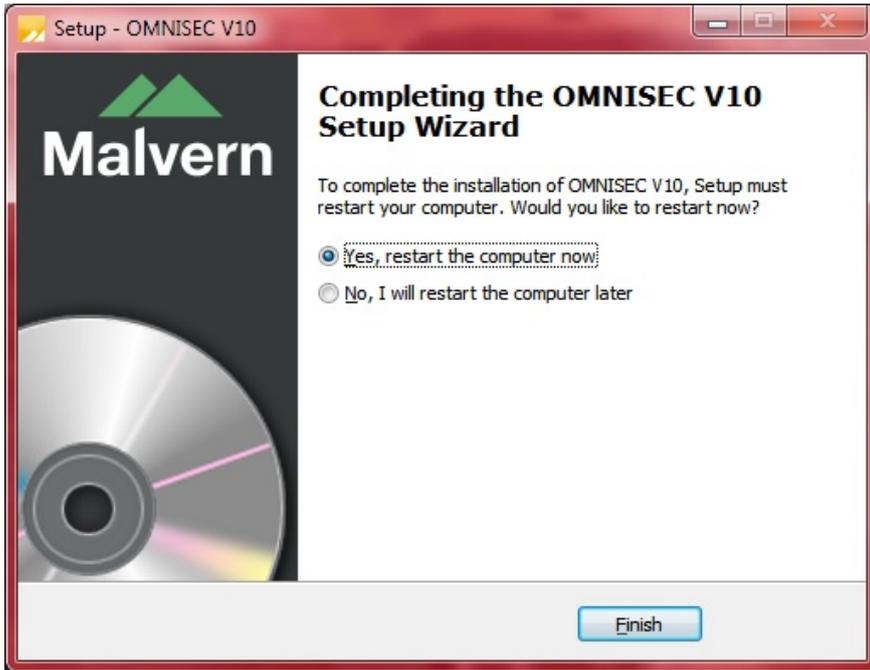
Once you have followed all of the instructions and clicked "Install", the software will begin the installation process.



If this is the first time OMNISEC v10 is being installed, it will install a 2014 SQL Server instance. This process may take 10 to 20 minutes to complete.



During the installation, a Windows Security window may appear asking if you would like to install a device. Select "Install" to install the device.



The last section in the Setup window will inform you that the computer needs to be restarted to complete the setup. Once the computer has been fully restarted, OMNISEC v10 will be ready to use.

Connecting the OMNISEC to a PC

With the software installed, the instrument should be connected via the USB port, and the system switched on. If present, turn on OMNISEC RESOLVE first and wait 30 seconds for it to complete its startup. Next, turn on OMNISEC REVEAL and wait 1 minute to allow it to complete its startup and connect to OMNISEC RESOLVE. When the two instruments are connected, the light bars on both instruments should be green.

If this is the first time connecting the instrument to the PC, there will be a message that appears saying “Installing device driver software”. The names used to identify the drivers will be “CMULx25”, “CHROME TAU A” and “CHROME TAU B”. Allow the installation to proceed and do not disconnect the instrument from the PC during this process. Once the drivers have been installed, Windows will indicate that they are ready to use.

Updating the OMNISEC REVEAL firmware

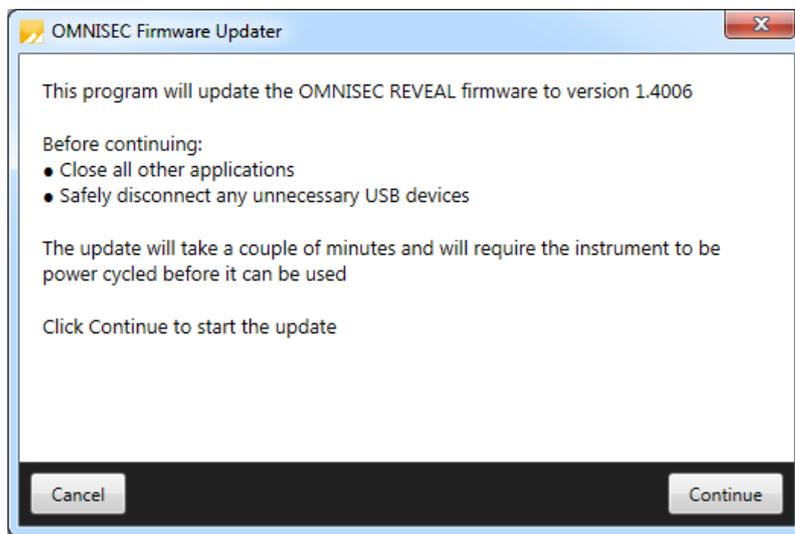
The first time you launch OMNISEC V10 with an OMNISEC REVEAL connected, you may be prompted to install a new version of the instrument firmware.



Updating the firmware improves system reliability and compatibility with the version of OMNISEC V10 you have installed on your computer. The update process takes up to 10 minutes and the instrument will need to be power cycled after the update is complete.

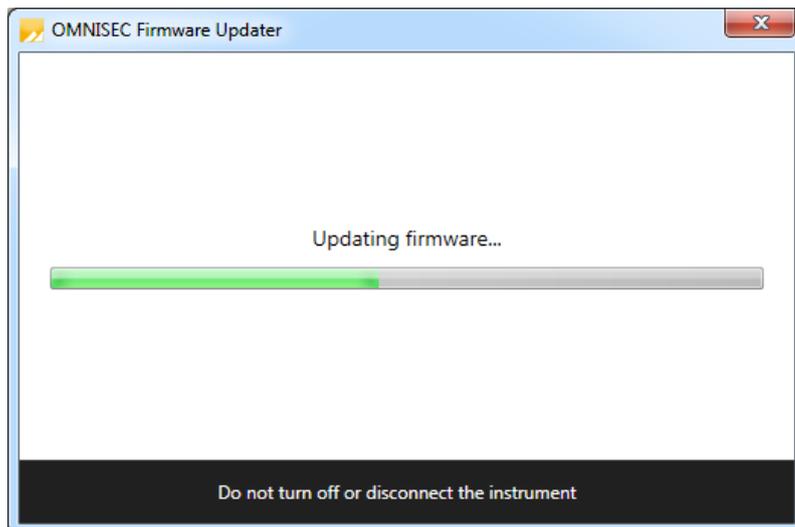
Click **OK** to close the OMNISEC application. After a few seconds the OMNISEC Firmware Updater program should automatically launch.

Note: If the OMNISEC Firmware Updater fails to open, or if the updater says it could not detect the instrument, disconnect the instrument, relaunch OMNISEC V10 and reconnect the instrument. You should receive the update prompt to try again.



When the OMNISEC Firmware Updater launches, you should carefully read the information presented to you before continuing.

Click **Continue** to start the update process.



The update process takes several minutes. It is recommended that you do not do anything else on your computer during the update.

After the firmware update has completed, you will be prompted to power cycle the instrument. Once the power up cycle has completed and you see the green lights on the front of RESOLVE and REVEAL click **Exit** and relaunch OMNISEC V10.

Note: If the update process fails, you don't see two green lights on the instrument or you are still being prompted to update the firmware, try power cycling the instrument and restarting OMNISEC V10 again.

Uninstall Procedure

The software can be uninstalled using the 'Programs and Features' section of the Windows 7 'Control Panel' or the 'Apps' section of Windows 10 'Settings'.

New Features

The new features in this release are listed in table 1.

Table 1: New features implemented in OMNISEC 10.40

Reference	Feature
76222	Windows 10 support
76203	Acquired sequences can now be deleted from the Data Selector window

Fixed issues

The main issues fixed in this release of the software are listed in table 2.

Table 2: Bug fixes implemented in OMNISEC 10.40

Reference	Issue	Comment
67103	Record View cannot be navigated with the keyboard up and down keys past an injection with analysis history	Fixed
72938	The Cancel button in the Calculation Method window does not function when creating a new calculation method	Fixed
73548	The Save button is enabled even when no changes have been made to an existing analysis	The Save button will no longer be enabled when no changes have occurred to an existing analysis
75980	When copying a Results by Peak report table, the Sample Name is not included	Fixed
77006	When attempting to add a new sample to a sequence, if invalid data has been entered, the cell focus does not reset to the cell with the invalid data	The cell focus will now automatically change to the first cell with invalid data
77867	The Export button in the Data Selector window does not become enabled after selecting data from the Advanced data selector window	Fixed

78071	After zooming in the PDA Data view, the 3D Spectrograph colour scale does not adjust to the visible range of data	The colour scale of the 3D Spectrograph now adjusts to the range of data currently visible
78636	When multiple analyses are selected in the Record View, making an adjustment to the baselines or limits of the 'active' analysis in the Raw Data view causes an extra item in the legend of the Distribution Plot to appear	Fixed
78673	Custom Sample Types are not saved to the Sample Types database when importing data	Any Sample Types used in imported data which do not exist in the Sample Types database are now automatically imported
78839	When expanding the analysis history of an injection in the Record View, the latest analysis for that injection gets selected	The latest analysis for an injection with analysis history is no longer loaded when expanding the analysis history, maintaining the current selection
79003	Partially selected sequences are not deselected when clicking the Deselect button in the Data Selector window	Fixed
79283	Record audit trails are not included in archives when exporting data	Audit trails for sequences, samples, injections and analyses are now included in archives when exporting data, and get imported when importing those archives
79286	Copying data from the Overlay Data chart has an incorrect CSV header	Fixed
80096	The legend in the Overlay Data chart displays incorrectly after making an adjustment to the active analysis	Fixed
80641	Importing an archive might get stuck and never complete	Fixed
81117	Copied charts are clipped when using a HiDPI display	Fixed
82349	On REVEAL-only systems, MALS data can drift out of sync across injections	Fixed
84017	Data export is not audited	Export operations are now recorded with a Record Exported event in the System Audit Trail and record audit trails of the data being exported. These events include the file path of the original export location
86325	Methods cannot be created on chromatograms with no light scattering detector	Fixed
86721	Errors analysing data when using a non-standard detector configuration	Fixed

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.

Table 3: Known issues in OMNISEC v10.40 software

Reference	Issue	Workaround
64935	Band broadening version 2 is set as default	When creating a method, in the Band broadening tab, use the drop-down menu to select which band broadening version to use
65981	The channel buttons in the Ribbon do not enable/disable appropriately for the selected view	Click on a different view, then click on the original view again
66441	The legend in the distribution and overlay charts may not be in the same order as in the Record View	No workaround
66452	Sequence & Sample name can be set to empty in the Edit Run Conditions window, causing errors in analyses	Do not set a blank sequence or sample name
66520 67578 67510	Creating and editing Column Sets does not work as expected	Delete and recreate column sets from scratch instead of trying to edit them
66897	MW values are rounded to 4 significant figures in the Calibration Standards database	No workaround
69856	Application crashes when trying to print or print preview a report where a table is longer than the page	Break long report tables into separate smaller tables
72004	Extrapolation may be incorrect with negative DP within limits	No workaround
72077	No ability to manually zoom out on the UV Spectrum Map chart	Use the Zoom Back and Full Zoom Out functions in the Ribbon Bar
72124	Crash in the Report Selector when clicking the Left arrow button when no report is selected in the right-hand pane	Make sure the report you want to remove is properly selected before clicking the Left arrow button
72239	The Delete button may be disabled for a user created Sample Type / Calibration Standard / Solvent	Rename the record, save and try again
72240	When deleting a Sample Type / Calibration Standard / Solvent, clicking the X button on the message box asking for confirmation will delete the record anyway	Do not click the X button on message boxes, always choose from the available buttons below the message
72681	Audit Trail event details cannot be copied	No workaround
72688	Current selection in the Data Selector window is lost when toggling the Exclude Baseline Monitors or Exclude Quick Runs options	No workaround

72707	Certain warning messages may be too long to fit in the status bar and become truncated, with the Reset button being inaccessible	If possible, make the window wider to see the whole text and the Reset button
72746	Changing the pump pressure units in Options may not automatically update the pump pressure displays in the reports	After changing the pump pressure units in Options, restart the software to ensure that all pressure displays use the new units
72751	The selected items in the Data Selector window does not stay in sync with the data currently open in the Record View	No workaround
72788	The View Audit Trail option is disabled when right clicking on a sequence in the Recently acquired sequences or Recently analysed sequences lists in the Data Selector window	Open any sequence that you want to view the audit trail for first. Right click on any injection from the sequence in the Record View and click View Sequence Audit Trail
72866	Find Limits fails where the RI signal is inverted and there is a single negative peak with no positive peaks	Set limits manually
72939	When editing an existing calculation method, the Cancel button in the Calculation Method editor window reverts any changes but does not close the window	Close the window by clicking the X button in the window title bar
73334	The Save button does not disable when relevant permissions expire and may cause a crash when clicked	Do not use time limits on permissions in the MAC
74039	No error is displayed when changing 21 CFR feature options if the user does not have sufficient privileges to change the feature setup	When setting up 21 CFR feature options, run the application as an Administrator
74280	The Concentration field in the Edit Run Conditions window does not allow enough precision for low concentrations	No workaround
74953	When importing archives with baseline monitors, the imported baseline monitors are displayed even though the Exclude Baseline Monitors option is selected	Deselect, then reselect the Exclude Baseline Monitors option to refresh the list
75977	The top of the MALS Fit Order view can be cut off	Resize the view
75978	The MALS fit order slider remains on the Raw data view after switching from a MALS method to a non-MALS method if MALS Model/Fit View window is active	Do not have the MALS Model/Fit View as an active window
75979	Baseline fit order may be incorrect for the number of set baseline points	Delete the baselines and reset them
75981	Legends in the overlay, distribution and derived data plots display nonsensical Analysis numbers to differentiate between multiple selected analysed records	No workaround
76127	When exiting the application after the instrument has been disconnected, the user is still asked whether they want to return the pump to standby flow rate	No workaround

76199	Opening baseline monitors with UV data in the PDA view may cause an Out of Memory crash	Do not open baseline monitors with UV data in the PDA view
76290	If a user clicks Cancel when asked if they want to save their changes to a sequence after attempting to exit the application, they are still asked if they want to return the pump to standby flowrate	No workaround
76528	Fraction collector triggers on using Baseline monitor	Disconnect fraction collector trigger lead from fraction collector or switch fraction collector off prior to using Baseline monitor function.
77868	Unrun samples are missing from partially run sequences when exported as OMNISEC 10.20 format archives	Export as 10.30 format archives where possible
78994	'Results by sample and peak' report table column headers get truncated	No workaround
79089	After a daylight savings time transition, the time displayed in report printouts still uses the old time zone	After the beginning or end of daylight savings, restart the software to ensure that the new time offset is used
79243	The application may crash when copying data from the Derived Data chart	No workaround
80018	The existing selection in the Data Selector window may be lost when selecting data from the Advanced data selector window	No workaround
80150	Warning messages may reset without the user clicking the Reset button	No workaround
80199	Sequences continue to run when the pump has been stopped due to a safety event	No workaround
80541	An injection run can be started when the pump is disabled due to a safety event	Ensure all safety events have been reset before starting an injection run
80724 80784	Copying data from the Derived Data chart produces incorrect CSV	No workaround
80766	IV polydispersity fit is calculated against MW polydispersity instead of IV polydispersity	No workaround
80783	Incorrect calculation method may be loaded when reloading calibrated data	No workaround
80844	Crash when switching between view modes in the Run Conditions Summary Report	No workaround
80910	Error when using the UV Wavelength slider in the Acquisition view before the data acquisition starts	Do not use the UV Wavelength slider until data starts being acquired

81100 88256	Application freezes when the viscometer is failing to zero	No workaround
82160	Application will fail to launch if certain configuration data gets corrupted	Reinstall the application. If the problem continues, contact your customer support representative
82353	Baselines may not appear on the Raw Data chart even though the points are set	Switch to a different record then switch back. If the problem remains, restart the software.
82355	The target flow rate is lost when clicking the Standby Flowrate button	No workaround
82358	MALS specific sample information, like laser power level, wavelength and noise, may not display correctly	No workaround
82359	Application may fail to launch if the instrument is currently in safety mode	Disconnect the instrument from the PC, launch the software, then reconnect the instrument
82361	Changing the active light scattering detector in an existing method creates an invalid method before it can be recalibrated	No workaround
83308	When exiting the application while editing a report with unsaved changes, the user may not be asked if they want to save their changes	Make sure any modifications to reports are saved or discarded before exiting the application
84020	The temperature controller of the MALS detector may unexpectedly turn off	No workaround
84099	Analysis error when setting close integration limits	No workaround
84680 86721	"Key Not Found" error when opening data for analysis	No workaround
84820	Crash when opening a report created by a different user	No workaround
84821	Non-alphanumeric characters are removed when setting a sequence name	Avoid using non-alphanumeric characters in sequence names
86583	After deleting the last run sequence, the samples for that sequence are still visible in the Open Sequence window from the Sequence view	No workaround
86783	The Save Report As function allows the user to save with the name of another existing report	No workaround
86785 88543 88565	Using the Search function with the Add function in any Database window may cause the application to crash or cause the window to become inoperable	If the window becomes inoperable, try closing the window, then reopen it. Do not use the search feature in these windows. Try sorting the data instead by clicking on the column headers

87328	Crash when quickly changing between different Themes in the Options window	Do not change the theme when any data is open or being acquired
87330	Crash when another Malvern Panalytical instrument is connected to the same PC as an OMNISEC	Do not connect any other instruments to the PC when an OMNISEC is connected
87729	"Cannot get cut point for the peak" error when calculating or calibrating after abandoning previous changes	No workaround
39602 88519	After extended periods of working with large datasets, the software may run out of memory	Save analyses regularly and restart the software if the problem occurs
88520 88616	Multiple data series may get tied to one axis in the Overlay Data chart	No workaround
88615	Retention volume may fail to calculate on the first sample of a sequence when using a method with manually set detector offsets	No workaround
89039	When changing the Sample Type for a sample in the Sequence editor, the highlighted Sample Type in the Database window is not the currently set Sample Type	No workaround
H 39617	After some crashes, the software may not restart	Disconnect the OMNISEC system from the PC and reconnect it. Alternatively, re-install the software. No data will be lost by doing this
H 39480	It is possible for the Zoom buttons in the ribbon bar to be inappropriately greyed out	Use the right click menu in the chart you wish to zoom out of
H 39446	The Copy button does not work in the Print Preview window	Copy the image directly from the report pages or the charts
H 39364	The Help button in the Report Selector does not function	Open the help file from the Start Page or another area
H 39315	Internet Explorer may block the help file from appearing depending on security settings	Allow Internet Explorer to display blocked content
H 39314	The search feature in the columns database only works for column sets, not columns	Manually search the columns list
	When sample type, dn/dc or concentration are updated, it updates this value for all injections in that sample, however, only the currently selected injection is recalculated	Select other injections for that sample and recalculate and save them to bring all the results up to date. This requirement is noted in the 'Edit Run Conditions' dialog

MALVERN PANALYTICAL

Groewood Road, Malvern,
Worcestershire, WR14 1XZ, UK

Lelyweg 1, 7602 EA Almelo,
Netherlands
P.O.Box 13, 7600 AA Almelo,
Netherlands

Tel: +44 1684 892456
Fax: +44 1684 892789

Tel: +33 546 534 444
Fax: +33 54 534 598

info@malvernpanalytical.com
www.malvernpanalytical.com

Disclaimer: Although diligent care has been used to ensure that the information in this material is accurate, nothing herein can be construed to imply any representation or warranty as to the accuracy, correctness or completeness of this information and we shall not be liable for errors contained herein or for damages in connection with the use of this material. Malvern Panalytical reserves the right to change the content in this material at any time without notice. Copyright: © 2018 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.