

SOFTWARE UPDATE NOTIFICATION (SUN) – OMNISEC SOFTWARE v10.31: PSS0051-05



MOLECULAR WEIGHT



MOLECULAR SIZE



MOLECULAR STRUCTURE

Introduction

This document details the installation and known issues with OMNISEC v10.31. This release is a maintenance release to fix a number of critical issues and improve performance.

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 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 64 bit Professional operating system.

Supported operating systems

OMNISEC software v10.31 is compatible with Windows 7. Only 64 bit Professional versions are supported.

Windows XP, Windows Vista, and Windows 8 and Windows 10 are not supported.

Supported Languages

English

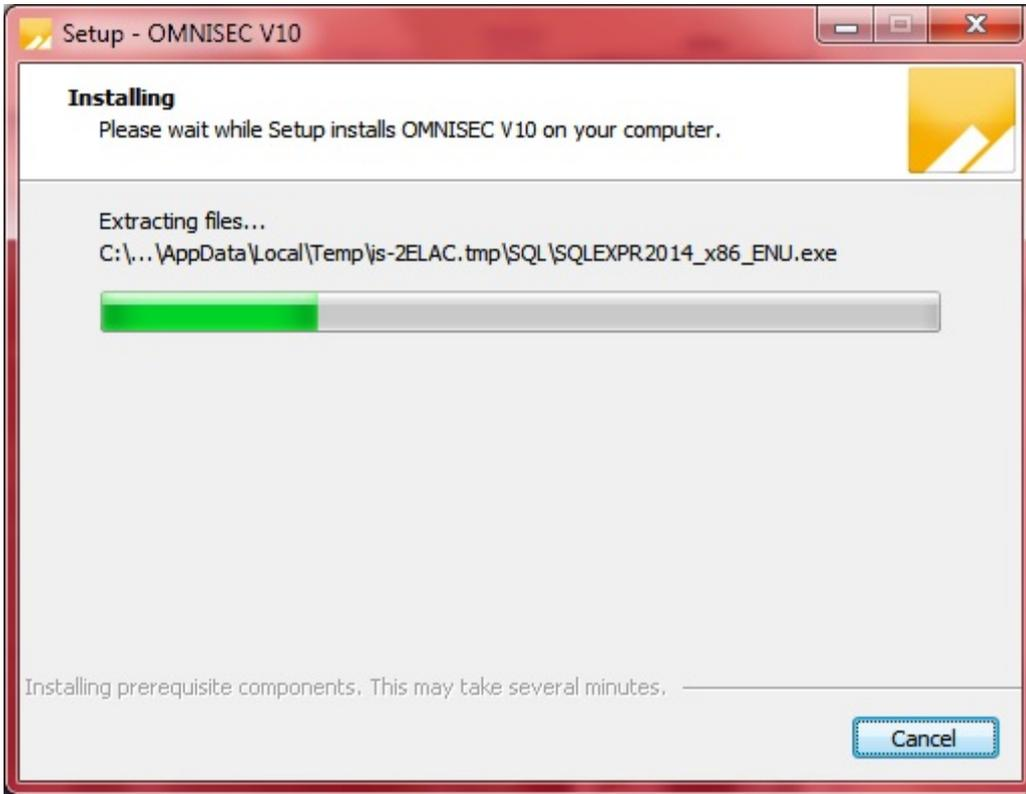
Installation Instructions

Windows 7 64 bit Professional

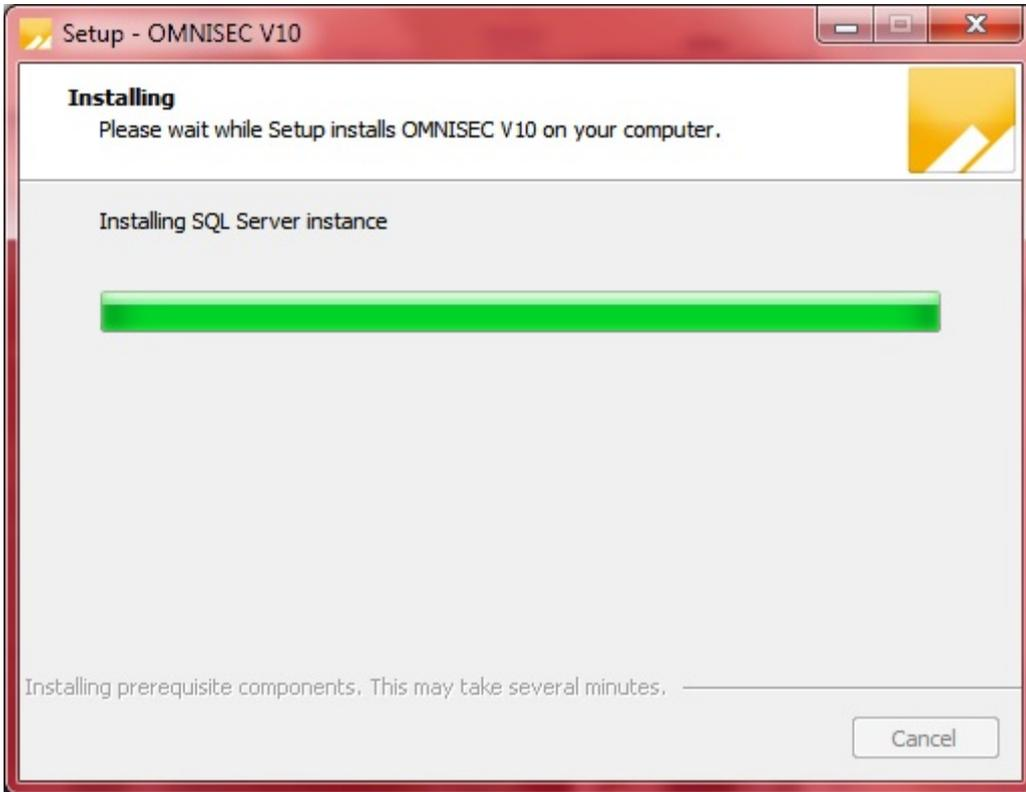
In most cases, OMNISEC should install automatically when the CD is inserted in the drive. If it doesn't, the user can double-click on the installer file, OMNISEC1031.exe, to begin the installation.



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.



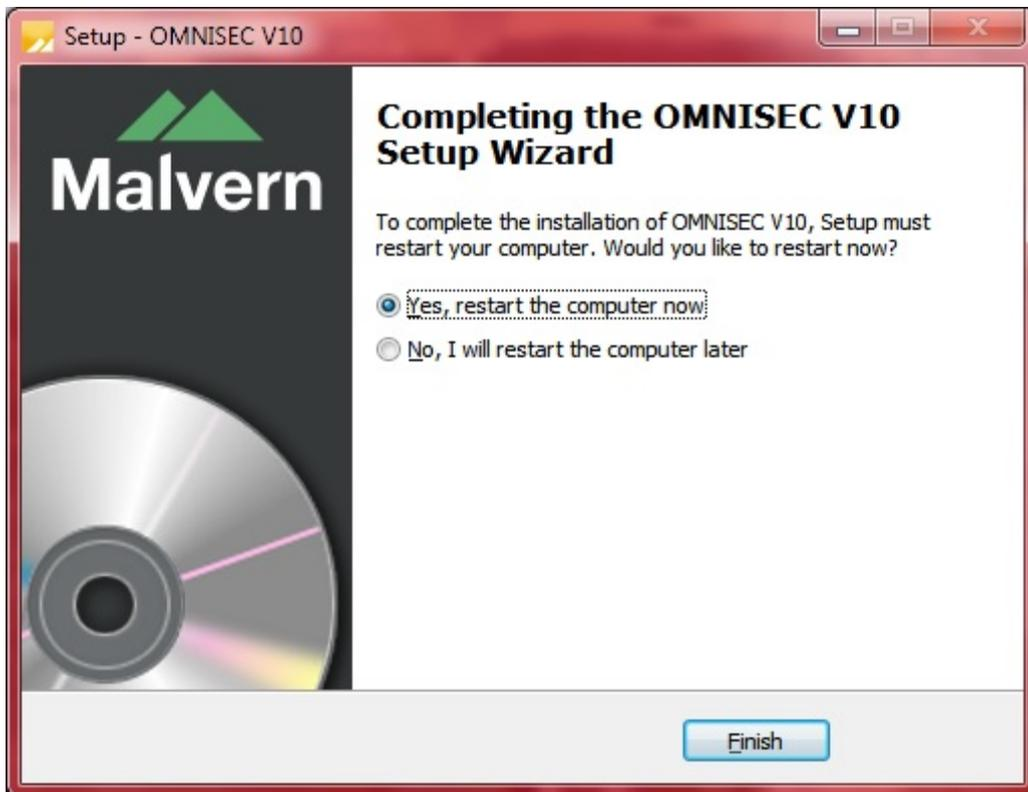
Once you have followed all of the instructions and clicked "Install", the software will begin the installation process.



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.

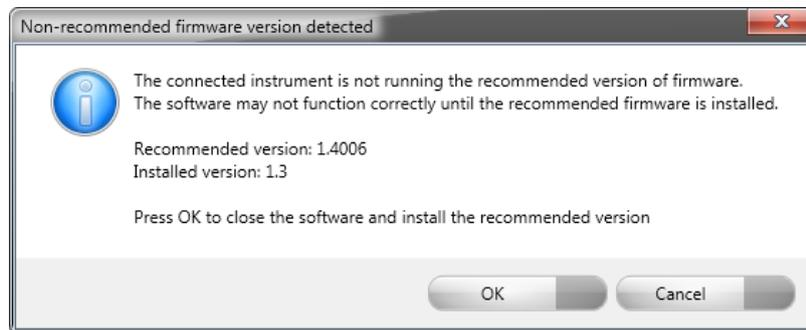
Connecting the OMNISEC to a PC

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

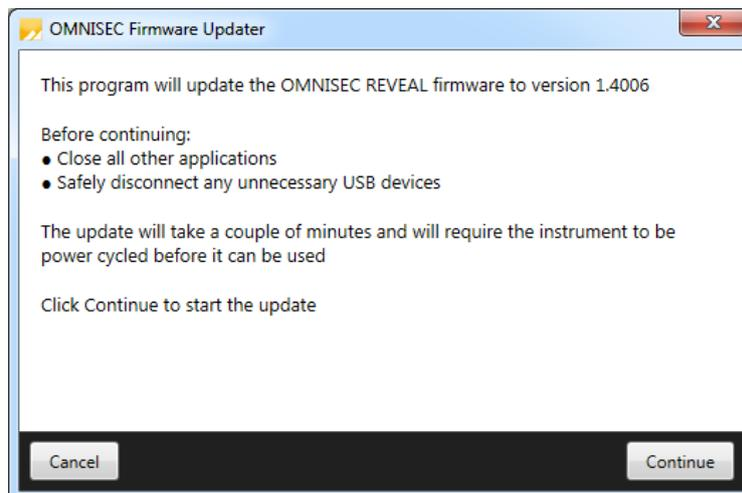
The first time you launch OMNISEC 10.31 with an OMNISEC REVEAL connected, you may be prompted to install a new version of the instrument firmware.



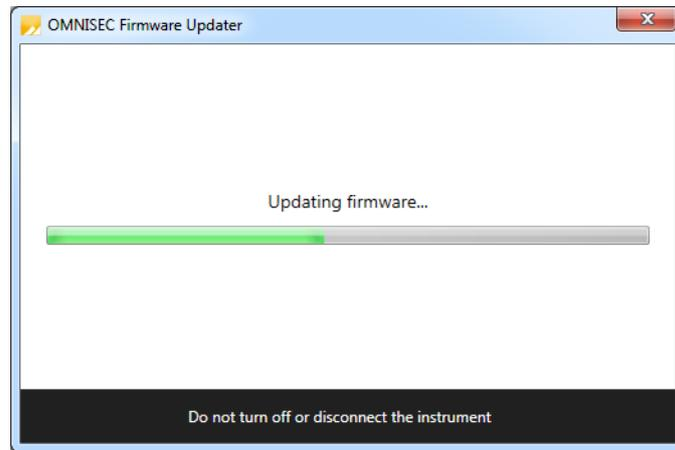
Updating the firmware improves system reliability and compatibility with the version of OMNISEC V10 you have installed on your computer. The update process takes up to 10 minutes and the instrument will need to be power cycled after the update is complete.

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'.

New Features

This release does not introduce any new features.

Fixed issues

The main issues fixed in this release of the software are listed in table 1.

Table 1: Bug fixes implemented in OMNISEC 10.31

Reference	Issue	Comment
66731	Sequences with a large number of samples are very slow to open	Performance when opening data for analysis has been improved
75008	Signature details may disappear in a report when the Abandon Changes button is clicked on an already signed analysis	Fixed
76857 77280 77281 77282	The application may become generally slow and unresponsive as more data is acquired	General performance improvements
77279	The Calculation Method browser window can take a long time to load	Performance improved
77784	Cannot import archives exported from OMNISEC software versions earlier than 10.20	Fixed
78135	Incorrect time zone displayed in report printouts	Fixed
78440	When opening acquired injections/samples from a sequence which is still running, the selection in the Data Selector window is lost when the next injection in the sequence completes	The selection is now maintained and newly acquired injections and samples from the running sequence are automatically selected
78784	When selecting an item in the Data Selector window, the related items in the sequence/sample/injection hierarchy may not automatically select/deselect as expected	Fixed
78802	On REVEAL-only systems, the live acquisition data is not plotted at the correct rate as defined in the Sequence editor	Fixed
79090	Changing the computer's time zone settings may cause false data integrity breaches	Fixed
79218	Crash when copying overlaid data from the Distribution Plot	Fixed
80533	Band broadening uses the wrong limits for calculating maxima	Fixed
80899	Unable to save analyses on data acquired from REVEAL units without a RALS/LALS detector or viscometer	Fixed
81199	Application freezes for several seconds when selecting an analysis if the Derived Data or Distribution Plot is enlarged	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 2: Known issues in OMNISEC v10.31 software.

Reference	Issue	Workaround
66520 67510 67578	Creating and editing Column Sets does not work as expected	Delete and recreate column sets from scratch instead of trying to edit them
82358	MALS specific sample information, like laser power level, wavelength and noise, may not display correctly	No workaround
82355	The target flow rate is lost when clicking the Standby Flowrate button	No workaround
80641	Importing an archive might get stuck and never complete	After waiting a reasonable amount of time where the progress bar has not changed, forcefully close the OMNISEC application using Windows Task Manager, then restart the software and try again.
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
78070	Report table columns may not be wide enough to display the full names of records	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.
76199	Running long measurements with the UV detector enabled may cause Out of Memory crashes	Restart the software before and after running an acquisition or baseline monitor with the UV detector enabled
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
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
75981	Legends in the overlay, distribution and derived data plots display nonsensical Analysis numbers to differentiate between multiple selected analysed records	No workaround
75980	When copying a Results by Peak report table, the Sample Name is not included	No workaround
75979	Baseline fit order may be incorrect for the number of set baseline points	Delete the baselines and reset them
75978	The MALS fit order slider remains on the Raw data view after switching from a MALS method to a non-MALS method	No workaround

75977	The top of the MALS Fit Order view can be cut off	Resize the view
74280	The Concentration field in the Edit Run Conditions window does not allow enough precision for low concentrations	No workaround
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.
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
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
72938	When creating a new calculation method, the Cancel button in the Calculation Method editor window is disabled	Close the window by clicking the X button in the window title bar to discard a new method
72866	Find Limits fails where the RI signal is inverted and there is a single negative peak with no positive peaks	Set limits manually
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
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
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
72707	The Reset button next to warning messages in the Status Bar may become clipped or hidden	Resize the window to make it wide enough to display the Reset button
72681	Audit Trail event details cannot be copied	No workaround
72242	The Malvern icon is missing next to pre-defined Sample Types	No workaround
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
72239	The Delete button may be disabled for a user created Sample Type / Calibration Standard / Solvent	Rename the record, save and try again
72004	Extrapolation may be incorrect with negative DP within limits	No workaround
69862	Report printing exhibits incorrect page setup behaviour	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
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

66441	The legend in the distribution and overlay charts may not be in the same order as in the Record View	No workaround
66262	When MALS is enabled, the software may refuse to start if the MALS unit is not connected to the PC or is not turned on.	If it is not possible to reconnect the MALS unit, disable the Network Interface Adapter used to connect to the MALS in the Windows Network Connections Control Panel. Open the software and then disable MALS in the Options window.
65981	The channel buttons in the Ribbon do not enable/disable appropriately for the selected view	No workaround
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.
39602	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
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.
39446	The Copy button does not work in the Print Preview window	Copy the image directly from the report pages or the charts.
39364	The Help button in the Report Selector does not function	Open the help file from the Start Page or another area.
39315	Internet Explorer may block the help file from appearing depending on security settings	Allow Internet Explorer to display blocked content.
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 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

© Malvern Instruments Ltd 2017

more information at www.malvern.com