

SOFTWARE UPDATE NOTIFICATION (SUN) – VISCOTEK SOFTWARE: CPS0038 – CPS0042



MOLECULAR WEIGHT



MOLECULAR SIZE



MOLECULAR STRUCTURE

Introduction

This document details the changes made to the OmniSEC software package in version 5.11. This release is a feature release which contains a small number of significant changes and improvements over version 5.10 as well as a selection of bug fixes.

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 Dual Core or hyper threaded processor, 2GB RAM, 150MB free hard disk space, 1024 x 768 screen resolution running in 16 bit color mode, CD-ROM drive, up to 3 serial ports and 1 free USB port depending on hardware configuration, Windows XP Pro (SP3 or later) operating system.

Recommended System Requirements

Intel Core 2 Duo, 2GB RAM, 160GByte hard disk drive, 1024 x 768 screen resolution running in 32 bit color mode, CD-ROM drive, up to 3 serial ports and 1 free USB port depending on hardware configuration, Windows 7 operating system.

Supported operating systems

OmniSEC 5.11 is compatible with Windows XP Pro (SP3) and Windows 7. Both the 32 bit and 64 bit versions of Windows 7 are supported.

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

Supported Languages

English

Installation Instructions

1. Windows XP

It is recommended that older versions of OmniSEC be uninstalled before version 5.11 is installed. This should be done following the uninstall procedure below. Double-click on the OmniSEC installer and follow the on-screen instructions.

When the Zetasizer μ V is connected, the USB driver will be installed and the following message will appear twice:



Figure1: Windows XP driver installation error message

This warning can safely be ignored as the software installation has been fully tested on Windows XP. If this warning is encountered the button, 'Continue Anyway' should be selected.

2. Windows 7

It is recommended that older versions of OmniSEC be uninstalled before version 5.11 is installed. This should be done following the uninstall procedure below. Double-click on the OmniSEC installer and follow the on-screen instructions.

Double-click on the OmniSEC installer and follow the on-screen instructions. When the Zetasizer μ V is connected, the USB driver will be installed and the following message may appear twice:



Figure2: Windows 7 driver installation error message

This warning can safely be ignored as the software installation has been fully tested on Windows 7. If this warning is encountered, 'Install this driver software anyway' should be selected.

3. Connecting the Zetasizer to a PC

With the software installed the instrument should be connected via the USB port, and the system switched on.

The Found New Hardware Wizard should open to install the drivers for the Zetasizer μ V. The name used to identify the hardware will be "Malvern Zetasizer".

This should be left with the default selection of 'Install the software automatically' and the 'Next' button should be selected so that file transfer begins. Once file transfer has completed the 'Finish' button should be selected to complete the installation.

You may see the same security warnings as described in the previous section. The warning can safely be ignored as the software installation has been tested on Windows XP and Windows 7. If the warning is encountered the 'Continue Anyway' or 'Install this driver anyway' option should be selected.

Windows will indicate that the new hardware is installed and ready to use.

Uninstall Procedure

The software can be uninstalled using the standard 'Add/Remove Programs' feature in the Windows XP 'Control panel' or in the 'Programs and Features' section of the Windows 7 'Control Panel'.

New Features

The new features added to the software are shown in table 1:

Table 1: New features added to OmniSEC 5.11

Reference[s]	Feature	Comment
PBI 39906	Network saving of data.	OmniSEC software can be set to save data to a network file location. OmniSEC 5.11 has been updated so that should the network connection be lost at the time the data is saved, that data will be saved locally, rather than lost.

Fixed issues

The main issues fixed in OmniSEC 5.11 are listed in table 2.

Table 2: bug fixes implemented in OmniSEC 5.11.

Reference(s)	Issue	Comment
PBI 29578	Return of autosampler to the HOME position.	The autosampler needle will now return to the HOME position after every injection. This reduces the occurrence of minor power fluctuations which could interfere with pump flow rate stability.
Bug 39579	Tools, Options crash.	Fixed a bug where opening the tools, options menu while there was no data open would cause the software to crash.

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 software.

Reference(s)	Issue	Workaround
Bug 37192	Result certification does not work. This feature was disabled in version 4.7.	Do not upgrade the software beyond version 4.6 if this feature is required.
Bug 36913	If results including a MALS detector are processed for conventional calibration, a baseline must still be set on the MALS detector channel.	Set a baseline on the MALS channel
Bug 36904	The calculated concentration when running a copolymer analysis will be reported as the input concentration unless the input concentration is set to zero, in which case, the calculated value is reported correctly	Set the input concentration to zero for copolymer sample analysis
Bug 37004	In the report designer, Rg(z) is incorrectly reported as zero.	There is currently no workaround for this bug
Bug 36918	In the report designer, if two detector channel line styles are modified at the same time, only one of them will be refreshed in the image	Modify one channel at a time
N/A	If the GPCmax is not connected to COM1 or one of the ports on the 6-way multi-COM port and the GPCmax toolpad is opened, the software can crash to windows with no error message.	THE GPCmax must be connected to COM1 or one of the ports on the 6-way multi-COM serial-to-USB converter.
N/A	False positives on the vapor sensor can occur if the threshold in the SEC-MALS settings is set too low.	The threshold should be set to a high enough level that false positives do not occur. The default value is 3000
N/A	When running the Zetasizer μ V with a computer on windows XP and certain processors, if the PC is shutdown, it will not shutdown but re-start.	The Zetasizer μ V should be turned off before the PC is shutdown.
N/A	MALS data from v5.00 re-calculated with 5.01 or 5.02 or 5.1 or 5.11 will be incorrect.	Any MALS data collected will need to be re-calibrated and re-calculated to obtain the correct answer.

N/A	Question marks replace some characters in the Model/Fit view if the system is running Chinese language settings	There is currently no workaround for this bug
------------	---	---

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 2013

more information at www.malvern.com