

SOFTWARE UPDATE NOTIFICATION (SUN)

– MicroCal Origin Add-On Dissociation Model Update

SOFTWARE: MCL6001



MOLECULAR STRUCTURE

Introduction

This document details the update of MicroCal Analysis software Dissociation model. 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) or Windows 7™ (64-bit), it has been tested under Windows 7™ (32-bit). Windows 7 is therefore the preferred operating system.

Table 1: Recommended system requirements for the MicroCal Analysis software.

Processor Type Intel Core2 Processor,
Memory 4GB
Hard Disk Storage 300 GB HDD,
Additional Storage Media DVD +/-R/RW drive
Display Flat Panel Monitor for software
Connectivity None
Operating System Windows 7™ (32 bit)

Supported Languages

The MicroCal Analysis software is available as an English language application.

Installation Instructions

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

Uninstall Procedure

The software cannot be uninstalled using the standard Add/Remove feature in the Windows 'Control panel'. It will be removed as part of the Origin Add-on Disk.

Fixed issues

The main issues fixed in the patch.

Reference(s)	Issue	Comment
00043	<p>There has been some debate as to the appropriateness of the dimer dissociation model employed by MicroCal in the data analysis software provided with the instruments. The discussion focused on the correct approach to account for the displaced volume upon each injection applied in the dissociation analysis. We have monitored the debate and revisited the expressions in MicroCal dimer dissociation model. In doing so, we have found an error in our approach and decided to use the expressions described by Alan Cooper et al (ref). The dissociation model supplied with all new instruments will be supplied with this new model and the manuals revised accordingly.</p> <p>The dimer dissociation model has not been widely applied in the literature but we acknowledge that this may impact some of the results in that published data. We would recommend that our customers review the literature data to ensure its correctness and relevant impact on future work. In most cases the discrepancies in the data will be within typical error limits for these types of experiments but there may be instances where this is not the case and as such, we recommend using the new model to check for this.</p> <p>Reference</p> <p>Deborah McPhail and Alan Cooper. J. Chem. Soc., Faraday Trans., 1997, 93(13), 2283-2289.</p>	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 for

Issue	Work around	Comment
Windows 7 Compatibility settings	<p><i>Run as Admin</i> needs to be applied to <i>Origin70.exe</i> and <i>MicroCalAnalysisLauncher.exe</i></p> <p>The <i>UAC (Universal Access Control)</i> feature of Windows should be set to <i>Never Notify</i> or project files (.opj) cannot be launched although the files can be opened directly from within Origin.</p>	Compatibility

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