

SOFTWARE UPDATE NOTIFICATION (SUN)

Viscotek DSV SOFTWARE: PSW0109



SOLUTION VISCOSITY

Introduction

This document details the changes made to the ARV software package for the Viscotek DSV in version 2.07. This release is a maintenance release and so includes only minor changes from version 2.06.

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.

Recommended System Requirements

PC running Windows XP 32 bit, Window 7 32 bit, Windows 7 64 bit.

Supported Languages

- English

Installation Instructions

Windows XP and Windows 7

It is recommended that ARV version 2.06 be uninstalled before version 2.07 is installed.

Run the installation package by starting the executable file.

Connecting the DSV to the PC

With the software installed, and DT3016 board installed, connect the Y501 as described in the Viscotek Y501 Viscometer Instrument Manual (MAN0423-2.0) and click on ARV icon to launch program.

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

No new features have been implemented in this version of the software:

Feature Modifications

The one change that has been made in this release is:

Reference(s)	Issue	Comment
	<p>The priority function in the sample queue prioritized sample dissolution but not sample measurement.</p> <p>This has been changed to prioritize sample dissolution and sample measurement ahead of other fully dissolved samples awaiting measurement.</p>	

Known Issues

There are no known issues in this version of the software.

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