

# PEAQ-DSC SOFTWARE: v1.60 (PSW0005-14) SOFTWARE UPDATE NOTIFICATION

## Introduction

This document details the release of PEAQ-DSC Control Software v1.60. This is a maintenance release to fix known issues and improve performance.

### Note:



Please check for any updates to this Software Update Notification at <https://www.malvernpanalytical.com/en/support/product-support/microcal-range/microcal-dsc-range/microcal-peaq-dsc>

## 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.

If a previous version of the software is detected, it will automatically be updated by this installer.

## Minimum System Requirements

Intel Core i5 Processor, 8 GB RAM, 300 GB hard disk drive, 1366 x 768 screen resolution.

## Supported operating systems

Windows 10.

## Supported Languages

The PEAQ-DSC Software suite is available as an English language application.

## Installation Instructions

The software is supplied as a self-executing installer from malvernpanalytical.com. Once downloaded, the installer must be run to set up the software from your local drive. If required, .NET 4.6 will automatically be installed.

If running in "Offline mode", a license key will be required to run the software. A license key can be exported from the instrument (PC controller must be connected to online instrument) or by contacting MicroCal. Instructions can be found in "Exporting a license file" section from the Software overview in the PEAQ-DSC User Manual.

## Uninstall Procedure

The software can be uninstalled using the standard Add/Remove feature in the Windows Control Panel.

## Known issues

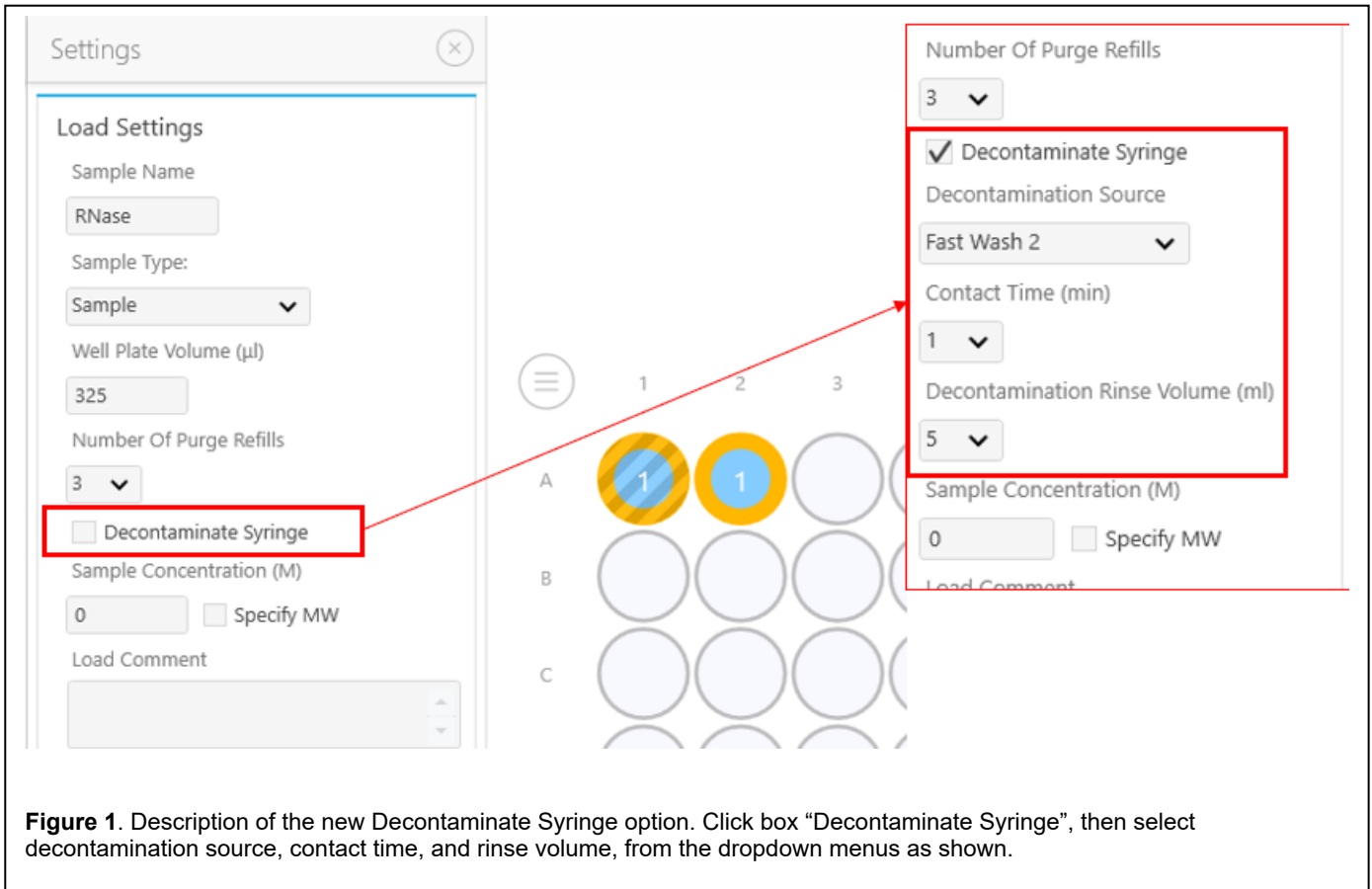
The known issues in this release of software are listed below.

Issue	Comment
Regional Compatibility	<p>To run the instrument, regional settings must be set to EN-US in the Windows 'Control panel'. NOTE: the system will be configured from the factory with EN-US settings.</p> <p>The software may be run offline using the user's preferred regional settings. Offline mode is the default installed configuration</p>
Out of Memory from large measurement file.	<p>Errors such as 'out of memory' or 'invalid format' might appear when a user attempts to load a too large .dmes file.</p> <p>While critical (.dmes) file size will vary for each PC due to different hardware, it is good practices to keep a .dmes file no larger than ~150 MB (~200 records) to ensure the file can be properly saved and loaded by the software. The user should be mindful of this when configuring sequences and targeting existing measurement files.</p>
Windows Taskbar	<p>When the Windows taskbar is set to auto-hide, maximizing the PEAQ-DSC software will prevent access to the taskbar. A user can disable the auto-hide of the taskbar.</p>
Disabled Windows Themes	<p>When Windows Themes service is disabled, the PEAQ-DSC software may render incorrectly. Users should keep Windows Themes service enabled.</p>
Incorrect clean message.	<p>Depending on how the system is stopped, the message instructing to clean manually, and saying "The system was scheduled to clean this...", may not be accurate. (In some cases, it was cleaned). If you are unsure, clean the instrument from the Maintenance workspace.</p>
Read-Only Measurement	<p>In rare instances, a measurement may be left as read-only after the completion of a sequence. The read-only flagged can be toggled through the measurement's file properties dialog accessed through Windows Explorer.</p>
Out of Memory from large sample raw data file.	<p>A very large raw data file (.dscx) may show an 'out of memory' or 'invalid format' warning when attempting to save. The critical (.dscx) file size will vary for each PC due to different hardware but will be approximately 25 hours of total scanning.</p> <p>There are safeguards in place to prevent this, but given the variability of the limit, the user should be mindful of this when configuring samples.</p>

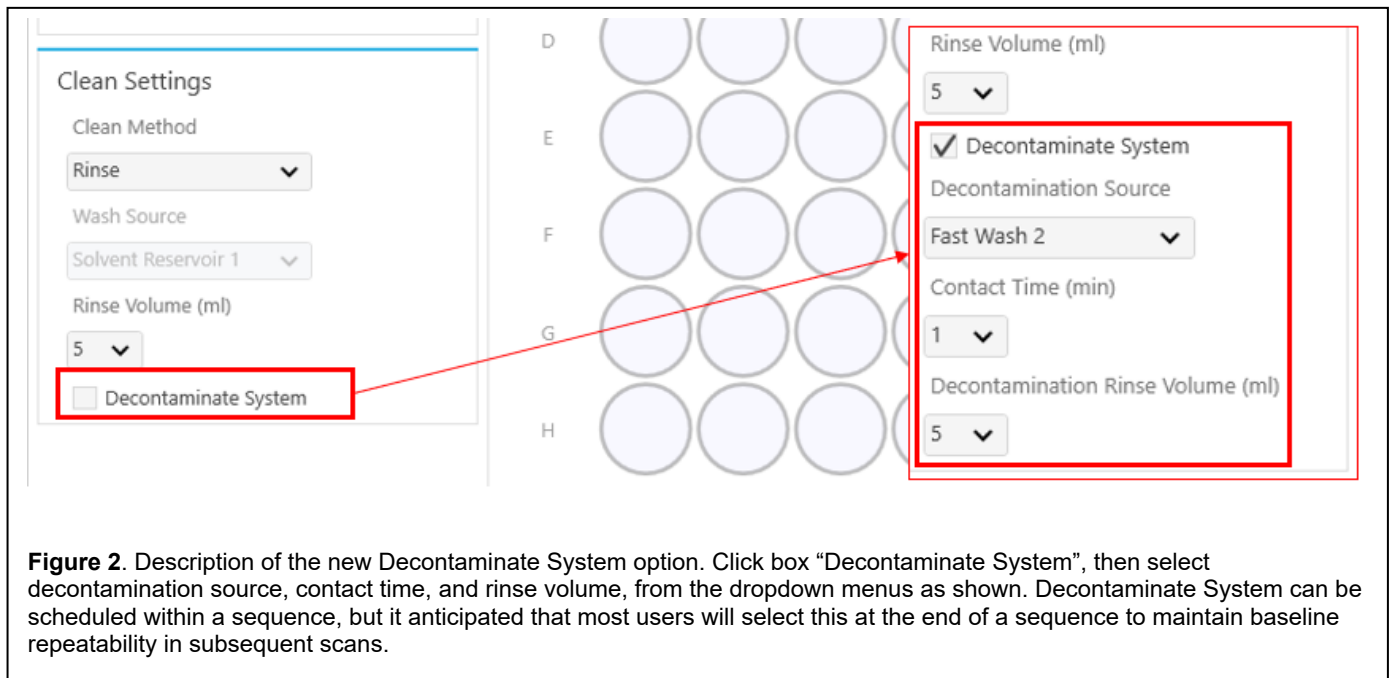
## Main issues

The main issues addressed in this release of software are listed below.

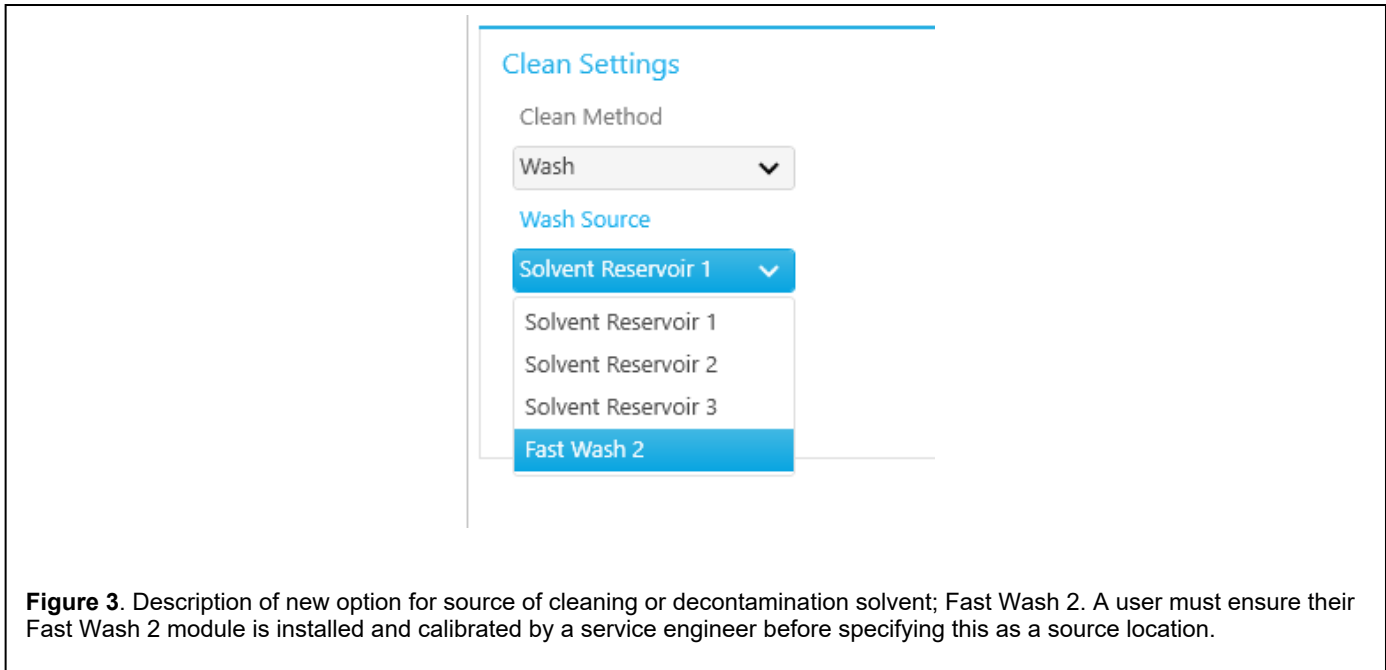
Issue	Comment
Decontamination option	<p>Using the existing autosampler, a user may choose to decontaminate their <i>syringe</i>, after loading the sample, and/or decontaminate the <i>entire system</i> at the end of a sequence. Choosing either options will permit the user to choose a location from which to load decontaminate solution and a contact time over which to soak either the syringe or the system.</p> <p><b>Syringe decontamination:</b> Here, the autosampler will finish loading the sample, initiate the scan, and begin decontaminating the syringe. Because of the concurrency of the scanning and decontaminating operations, it is suggested the decontamination's duration (i.e. contact time) be shorter than the samples' scans. This constraint is validated during SOP definition. See Figure 1, below.</p> <p><b>System decontamination:</b> Here, the autosampler will finish the specified cleaning method and begin decontaminating the system (syringe and cells). Because the system decontamination will likely delay the next sample load, it is suggested any system decontamination is scheduled at the end of a sequence. In the event a system decontamination is scheduled mid-sequence, any baseline repeatability will be disrupted. If applicable, this warning will be communicated to the user during SOP definition. See Figure 2, below.</p> <p><b>Fast Wash 2:</b> A user must ensure their Fast Wash 2 module is installed and calibrated by a service engineer before specifying this as a source location.</p>
Wash from Fast Wash 2	<p>Traditional wash procedures, and the newly added decontamination option, can utilize Fast Wash 2, if the module is configured. If not, please contact your Malvern Panalytical specialist. See Figure 3, below.</p> <p><b>Fast Wash 2:</b> A user must ensure their Fast Wash 2 module is installed and calibrated by a service engineer before specifying this as a source location.</p>
Failed to save data	<p>A bug was fixed wherein folder permissions were being checked but were not specific to the current user. Data may have failed to be saved without warning.</p>
Permission file	<p>A bug was fixed wherein the Malvern Access Configurator's permissions.xml file did not display the current software version. This did not affect function but was a cause for confusion.</p>
Exception while editing running sequence	<p>A bug was fixed wherein a user may interrupt or lose sight of the currently running sequence's scheduled list. This may happen if repeatability rescans or detergent scans have already been completed, the user has pending changes to the sequence, and the user chooses to select a different sequence (or create a new sequence). If the pending changes are discarded after being prompted, the bug described may occur.</p>
Sample duration validation warning	<p>A bug was fixed wherein sample duration was validated against a 25 hour limit (see known memory issue), but the exception notification indicated a recommended maximum duration of 15 hours. These values were made consistent at 15 hours.</p>



**Figure 1.** Description of the new Decontaminate Syringe option. Click box “Decontaminate Syringe”, then select decontamination source, contact time, and rinse volume, from the dropdown menus as shown.



**Figure 2.** Description of the new Decontaminate System option. Click box “Decontaminate System”, then select decontamination source, contact time, and rinse volume, from the dropdown menus as shown. Decontaminate System can be scheduled within a sequence, but it anticipated that most users will select this at the end of a sequence to maintain baseline repeatability in subsequent scans.



## MALVERN PANALYTICAL

Malvern Panalytical Ltd.  
Groewood Road, Malvern,  
Worcestershire, WR14 1XZ,  
United Kingdom

Malvern Panalytical B.V.  
Lelyweg 1, 7602 EA Almelo,  
Netherlands

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

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

[info@malvernpanalytical.com](mailto:info@malvernpanalytical.com)  
[www.malvernpanalytical.com](http://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: © 2021 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.

MP DisCop MP DisCop MP DisCop MP DisCop MP DisCop MP DisCop