

PEAQ-ITC Automated Control SOFTWARE: v1.21 (PSW0002-3) SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the installation and known issues with PEAQ-ITC Automated Control v1.21. This is a maintenance release to fix known issues, improve performance and establish Windows 10 compatibility.

Minimum System Requirements

Intel Core i5 Processor, 8 GB RAM, 250 GB hard disk drive, 1366 x 768 screen resolution.

Supported operating systems

Windows 7 and Windows 10 are supported.

Supported Languages

English

Installation Instructions

The software suite comes on an USB drive or may be downloaded from MalvernPanalytical.com. You must run the setup program, Setup MicroCal PEAQ-ITC Automated Control Software, from your local drive and only from an account with administrative privileges. If a previous version is detected, it will be automatically removed by the installer.

The driver installers will automatically be launched from this installer.

Uninstall Procedure

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

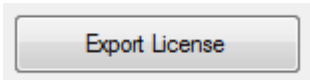
New Features

Single Application

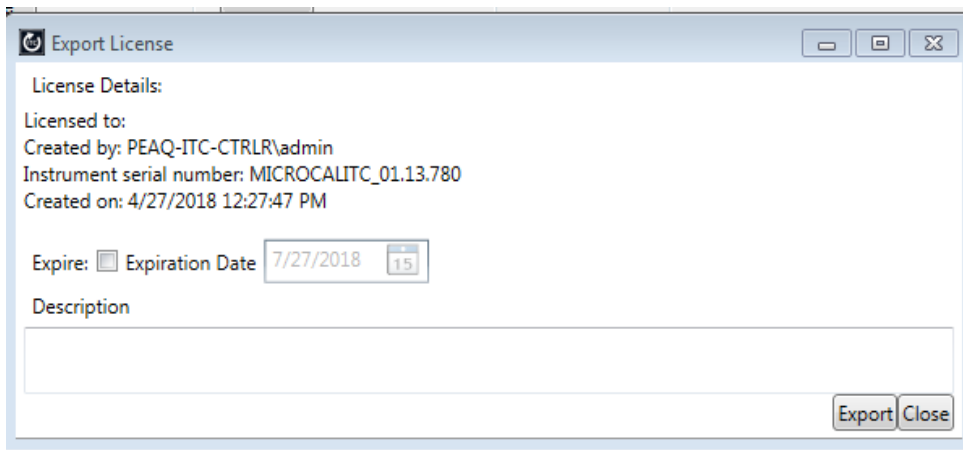
- This software now runs as a single application. The PEAQ-ITC Automated Control Software will now be the single source of instrument interaction.
- To ensure you open the most recent version of the software, please remove any old shortcuts to the MicroCal Automated Operation Software.
- Please note that some operations are now found in the System tab of the new version of the software (i.e., pipette operations, temperature control). The old software can still be shown/hidden if desired. Updating run parameters during a titration still require the old software.

Licensing

- The latest version (and v1.20) of the PEAQ-ITC Analysis software requires a license. While a PEAQ-ITC instrument is online, a license can be generated from the System Tab.

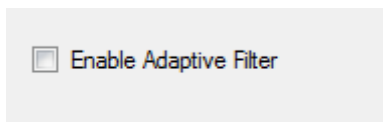


- Clicking the “Export License” button will show the following dialog for creating a license:

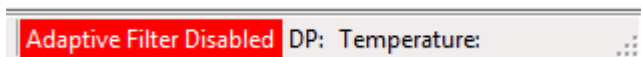


Adaptive Filter Disabling

- The Adaptive Filter was created to automate portions of data analysis and create smoother baselines, resulting in more data-rich sampling during peak events and more accurate peak data than legacy filter parameters.
- Some users have requested the ability to use legacy filter parameters on PEAQ-ITC instruments to systematically compare datasets.
- It is now possible to navigate to the System tab and check the “Adaptive Filter” box to disable the adaptive filter.



- The software will need to be restarted once the filter setting is changed.
- When the Adaptive Filter is disabled, users are notified by a red warning label in the instrument status bar.



- When the filter is disabled, the user will be permitted to specify a Filter Period in the experimental parameters.

Method Parameters

of Injections

Cell Temperature (°C)

Reference Power (µcal/s)

Initial Delay (s)

Stir Speed (rpm)

Filter Period (s)

Feedback Mode/Gain

None
 Low
 High

Known issues

The known issues in this release of the software are listed in Table 1.

Work Around	Issue	Comment
	Tube Rack Lid	It's possible to close the tube door when the lid is out of place. The software correctly will not start in this state; however, the warning dialog message quickly disappears before the user is able to read it.
	Skip Pipette Clean and Load Validation	The validation does not check if the pipette volumes in the experiments can be accommodated by the single pipette load.
Resave the method without special characters.	SIM Method	The SIM method contains special characters and will not validate.

Table 1: The known issues in PEAQ-ITC Automated Control Software v1.21

Fixed issues

The main issues addressed in this release of the software are listed in Table 2.

Reference	Issue	Comment
1909	Operation on Windows 10, 64 bit	The software and drivers are compatible with Windows 10.
3948	License for PEAQ-ITC Analysis	The PEAQ-ITC Analysis software will now require a license. There is a button on the System tab to export a license. This button will only be enabled if the instrument is online.
4570	Adaptive Filter Option	It is now possible to collect data with the adaptive filter disabled. When the adaptive filter is disabled, the instrument filtering will revert to the iTC200/VPITC style filter period. A warning label will be displayed when adaptive filter is disabled. The analysis software will have a visual designation of files that were run with Adaptive Filter disabled.
4921	Manual Log Entry	The user can manually enter any statement and write it to the AutoPEAQITC.log.
5199	Export license button	A bug was fixed where the button might remain disabled.
5196	Service Tool	A bug was fixed where the Create Support Files tool did not extract the serial comm logs.
4778	PEAQ-ITC Library	A convenient PEAQ-ITC library is created for accessing data, logs, and settings. A library is simply a shortcut that is visible in Windows explorer. Libraries may need to be made visible by right-clicking in Windows Explorer and selecting 'Show all libraries'.

Table 2: Main Issues addressed in PEAQ-ITC Automated Control Software v1.20 and v1.21

MALVERN PANALYTICAL

Groewood Road, Malvern,
Worcestershire, WR14 1XZ, UK
(MP addbody)

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

Lelyweg 1, 7602 EA Almelo,
Netherlands
P.O.Box 13, 7600 AA Almelo,
Netherlands

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

info@malvernpanalytical.com
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: © 2018 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 MP DisCop