

# PEAQ-ITC Analysis SOFTWARE: v1.21 (PSW0051-3) SOFTWARE UPDATE NOTIFICATION

## Introduction

This document details the installation and known issues with PEAQ-ITC Analysis 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](http://MalvernPanalytical.com). You must run the setup program, Setup MicroCal PEAQ-ITC Analysis Software, from your local drive and only from an account with administrative privileges. If a previous version of the software is detected, it will automatically be removed by the installer.

## Uninstall Procedure

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

## New Features

Licensing

- This version (and v1.20) of the PEAQ-ITC Analysis Software will require a license.
- Please see the [PEAQ-ITC Control Software Update Notification](#) or the [PEAQ-ITC Automated Control Software Update Notification](#), as applicable for your instrument to generate 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.
- Please see the [PEAQ-ITC Control Software Update Notification](#) or the [PEAQ-ITC Automated Control Software Update Notification](#), as applicable for your instrument.

## Fixed issues

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

Table 1: Main issues addressed in PEAQ-ITC Analysis v1.20 and v1.21

Reference	Issue	Comment
184	Leaving Fit workspace	Editing the fitting parameters required the user to click the Fit button before leaving the workspace. This is no longer required. Instead, the user is prompted to update their result upon leaving.
189	Results Table [syr] Column Fix	An issue was fixed in Results Table where [syr] column was inaccurate for data marked as 'Check Data'.
198	Enzyme Fitting Errors Scaling Fix	An issue was fixed where fitting results errors were not scaled properly.
3108	Zoom on Empty Chart Fix	A bug was fixed where an exception would be thrown when zooming on a chart with no curves displayed.
3948	License for PEAQ-ITC Analysis	This version of the PEAQ-ITC Analysis Software will require a license. A license can be exported from either PEAQ-ITC manual or automated control software when 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.
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'.
4884	Enzyme Kinetics Inhibitor toggling fix	An issue was fixed where toggling Kinetics Inhibitor in the Enzyme fitting model (toggled back to original state) would not always yield the same fit as the initial fit. The fitting algorithm was optimized and will now yield consistent results after toggling back to the same state.

## MALVERN PANALYTICAL

Groewood Road, Malvern,  
Worcestershire, WR14 1XZ, UK

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