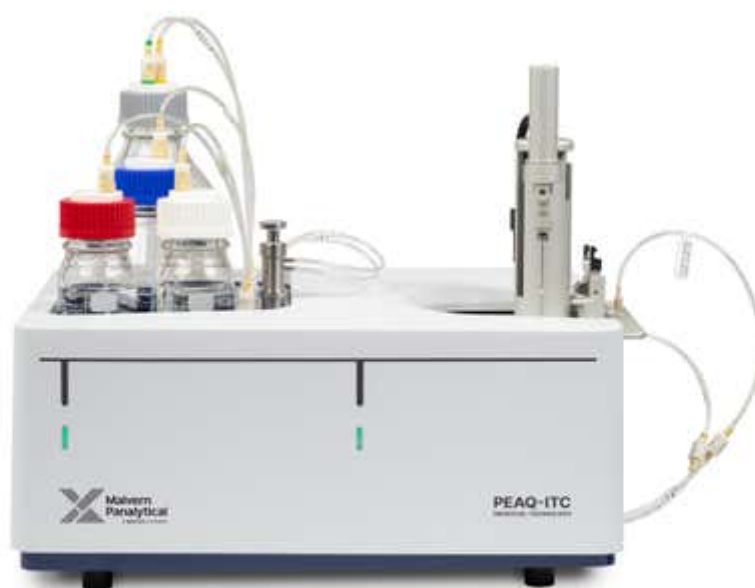


PEAQ-ITC Analysis SOFTWARE: v1.52 (PSW0051-9)

Software Update Notification



Introduction

This document details the release of PEAQ-ITC Analysis software v1.52 (PSW0051-9) for the MicroCal PEAQ-ITC instrument. It covers software issues that have been fixed and introduces new features.

The information within this document is required to perform a risk analysis to determine if the software should be installed. In this risk analysis the benefits of the new features provided and resolved software issues must be weighed against the risk of new issues that may be introduced to vital areas of the software or possible changes to the results of future analysis. Installation instructions are provided.

Note:

Please check for any updates to this Software Update Notification at www.malvernpanalytical.com/en/support/product-support/microcal-range/microcal-itc-range/microcal-peaq-itc

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. The software has been tested under Windows 10 and Windows 11.

Table 1: Recommended system requirements for the PEAQ-ITC Analysis software

Feature	Specification	Feature	Specification
Processor Type	Intel Core i5 Processor	Display	1366 x 768 screen resolution
Memory	8GB RAM	Operating System	Windows 10, Windows 11
Hard Disk Storage	250GB HDD	USB Ports	2 free USB ports

Supported Languages

The PEAQ-ITC Analysis software is only available as an English language application.

Installation Instructions

The software suite may be downloaded from 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.

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'.

Changes

The main updates included in this software release are listed in Table 2. These include new features, improvements, and fixed issues.

Table 2: Changes implemented in PEAQ-ITC Analysis software version 1.52

Reference	Change	Details
343947	.NET 8 upgrade	Updated manual PEAQ-ITC control and analysis software to target .NET 8 and include the updated runtime in the installer package.
331611	Windows 11 support	Updated the PEAQ-ITC manual installer to automatically install the latest (6.74) version of Instacal, enabling Windows 11 support.
228067	Matlab runtime installer improvements	Run the Matlab Runtime installer in the foreground and retain its executable in the program directory, improving error visibility and enabling manual re-execution.
228120	Enzyme MIM baseline fitting	Added a configurable offset method to enzyme MIM baseline fitting, defining the offset as the average initial delay and auto-recalculating results upon change.
228108	Windows analysis service installation	Updated the analysis service to support local Windows installation, allowing clients to submit experiment models for titration fitting and receive results.
228117	Wash module feedback improvements	Enhanced the vacuum-leak error message with additional guidance and disabled sensor-gating when detergent is used to maintain signal reliability.

This page is left intentionally blank



Groewood Road, Malvern,
Worcestershire, WR14 1XZ,
United Kingdom

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

Lelyweg 1,
7602 EA Almelo,
The Netherlands

Tel: +31 546 534 444
Fax: +31 546 534 598

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: © 2025 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.