

# PEAQ-ITC Automated Control SOFTWARE: v1.52 (PSW0002-9)

Software Update Notification



## Introduction

This document details the release of PEAQ-ITC Automated Control software v1.52 (PSW0002-9) for the MicroCal PEAQ-ITC Automated 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-automated](http://www.malvernpanalytical.com/en/support/product-support/microcal-range/microcal-itc-range/microcal-peaq-itc-automated)

## 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 Automated Control software

Feature	Specification	Feature	Specification
<b>Processor Type</b>	Intel Core i5 Processor	<b>Display</b>	1366 x 768 screen resolution
<b>Memory</b>	8GB RAM	<b>Operating System</b>	Windows 10, Windows 11
<b>Hard Disk Storage</b>	250GB HDD		

## Supported Languages

The PEAQ-ITC Automated Control 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 Automated Control 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 Automated Control software version 1.52

Reference	Change	Details
<b>343406</b>	Simultaneous pump operation	Fixed an issue in which the waste pump waited for the syringe pump to finish when a script used the “Do not wait for operation to complete” flag, ensuring both pumps operate concurrently as expected.
<b>344418</b>	.NET 4.6.1 runtime removal	Removed the bundled .NET 4.6.1 runtime from the installer, relying on the controller PC’s existing framework.
<b>346132</b>	Final TA Valve Rinses overflow	Fixed an issue where the Final TA Valve Rinses script overflowed the wash and load station by not emptying it before refilling and by dispensing the titrant arm cannula line into the station; the station is now emptied first, and the cannula line is drained to waste.
<b>348097</b>	Remove purge/refill functionality	Removed unused purge-refill option to increase robustness and 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](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: © 2025 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.