



PEAQ-ITC Analysis SOFTWARE: v1.41 (PSW0051-7) SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the release of software PSW0051-7: v1.41 of the software for the PEAQ-ITC system. It covers software issues fixed and new features introduced. This information 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.

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.

Minimum System Requirements

Intel Core i5 processor, 8 GB RAM, 250 GB free hard disk space, 1366 x 768 screen resolution, 2 free USB ports.

Supported operating systems

Windows 7 and Windows 10 are supported.

Supported Languages

English

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 'Programs and Features' section of the Windows 'Control Panel'.

New Features

This release does not introduce any new features.

Work items

The main items addressed in this release of the software are listed in table 1.

Table 1: Items addressed in PEAQ-ITC Analysis Software v1.41

Reference	Issue	Comment
6316	Valid path check	Paths are being checked more rigorously when being entered as a preference. If the current user does not have write access to the desired path, the entry box will display the error, rather than ultimately saving a file to the 'recovery' location if write access is denied.
6340	Method folder creation	A bug was fixed wherein the public Methods folder was not being created during installation. This prevented it from being the default path on initial startup.
6341	Indicate support files created	A dialog box will now appear once the support files are saved to the desktop.

This page is left intentionally blank

MALVERN PANALYTICAL

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

Lelyweg 1,
7602 EA Almelo,
The Netherlands

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

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: © 2021 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.