

# PEAQ-ITC Analysis SOFTWARE: v1.30 (PSW0051-5) SOFTWARE UPDATE NOTIFICATION

## Introduction

This document details the installation and known issues with PEAQ-ITC Analysis Software v1.30. This is a maintenance release to fix known issues and improve performance.

## 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.

## 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 Software v1.30

Reference	Issue	Comment
6039	Incorrect .ini file	This version maintains compatibility with the new control software that fixes a bug wherein the wrong .ini configuration could be used if the PC is disallowed access to the instrument's onboard configuration.

## MALVERN PANALYTICAL

Groveswood 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  
[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