

PEAQ-ITC Analysis SOFTWARE: v1.40 (PSW0051-6) SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the installation and known issues with PEAQ-ITC Analysis Software v1.40. 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. 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
4779	Export Support Files	This version fixes a bug around exporting support files using the Create Support Files utility. This utility is removed and replaced by a button, named "Create Support Files", in the Service Settings view.
6049	Maximum DP Bin Setting	This version removes to maximum DP value as a binning criterion.

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
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: © 2020 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