

Halo Manager Software: v2.23 (PSW0321-01)

SOFTWARE UPDATE NOTIFICATION

Introduction

This document details the release of software PSW0321: v2.23 of the software for the Halo system. Installation instructions and PC requirements 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.

Recommended System Requirements

The recommended computer system requirements for running this software are highlighted in table 1 below.

Table 1: Recommended system requirements for the Halo Manager software.

Feature	Specification
Processor Type	Intel Core i5 processor
Memory	8 GB RAM
Hard Disk Storage	300 GB free hard disk space,
Display Resolution	1920 x 1080 full HD screen resolution
Connectivity	1 free USB port,
Operating System	Windows 10 64 bit Professional operating system. * Preferred Operating System

Supported operating systems

Halo Manager software v2.23 is compatible with Windows 7 32 and 64 bit, and Windows 10 64 bit.

Supported Languages

- English

Installation Instructions

The software is supplied as a self-executing installer. Once downloaded, the 'Install Halo Manager' link must be run to set up the software. It is not necessary to uninstall a previous version of the software before running the installer.

Uninstall Procedure

The software can be uninstalled using the 'Programs and Features' section in 'Control Panel'.

Fixed issues

Table 1: Bug fixes implemented in Halo Manager v2.23

Reference	Issue	Comment
ASDHAL-1	Using more than 32 characters in a project name could cause a Halo to crash.	The project name field is now limited to 32 characters.
ASDHAL-4	Using Special Characters during project or location creation could cause a white screen crash of the Halo.	The following characters are now disallowed for Project Name or Location Name: < > [] { } ? * : " / \ ' @ ~ `

Known Issues

The following software bugs have been discovered within the software, and will be investigated as part of a future release. Please follow the suggested work-around.

Table 2: Known issues in Halo Manager v2.23 software.

Reference	Issue	Workaround
ASDHAL-10	Date format other than MM/DD/YYYY can cause synch to fail.	
ASDHAL-11	NAD83 GPS format does not display correctly in Halo Manager	
ASDHAL-17	It is not possible to upload new sample libraries to the Halo Instrument.	

MALVERN PANALYTICAL

Malvern Panalytical Ltd.
Groewood Road, Malvern,
Worcestershire, WR14 1XZ,
United Kingdom

Malvern Panalytical B.V.
Lelyweg 1, 7602 EA 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: © 2018 Malvern Panalytical. This publication or any portion thereof may not be copied or transmitted without our express written permission.

