

MSD[®] DISCOVERY WORKBENCH 4.0 RELEASE NOTES

Software Overview

DISCOVERY WORKBENCH[®] software is designed for acquiring and analyzing assay data using Meso Scale Discovery[®] instruments. Version 4.0 includes new features to increase efficiency and improve the quality of your data. The release notes below cover all updates since software launch, beginning with the most recent update.

DISCOVERY WORKBENCH 4.0 is also available in a desktop version for office PCs. The desktop edition includes all of the analysis, database management, and data integrity tools that are included in DISCOVERY WORKBENCH 4.0 for MSD instruments. The only difference is that the desktop version cannot be used to control an MSD instrument.

Released Software Versions

- DISCOVERY WORKBENCH 4.0.12.1
- DISCOVERY WORKBENCH 4.0.12
- DISCOVERY WORKBENCH 4.0.11
- DISCOVERY WORKBENCH 4.0.9 (original update from version 3.0)

DISCOVERY WORKBENCH 4.0.12.1

New Features

- Added support for the MESO[®] SECTOR S 600.

DISCOVERY WORKBENCH 4.0.12

Enhancements

- Prevents accidental installation of the desktop version of DISCOVERY WORKBENCH 4.0 onto MSD instrument computers. Installing the desktop version on the instrument computer can cause loss of data.
- Improved communications, particularly with respect to the stability of communications between the MSD instrument computer and the instrument.
- Updated contact information for customer support.

DISCOVERY WORKBENCH 4.0.11

New Features

- Added support for data acquisition using the new MESO QuickPlex[®] SQ 120 instrument.
- Enabled analysis of QuickPlex plate formats.

Enhancements

- Updated splash screen that appears on software initiation.
- Updated electronic version of the instrument manual and software user guide (see help menu).

Fixes

- Fixed miscellaneous bugs.



MSD[®] DISCOVERY WORKBENCH 4.0 RELEASE NOTES

DISCOVERY WORKBENCH 4.0.9

New Features

- A Configuration Manager profile was added with administrator rights. The Configuration Manager can control the access and permissions of other users and configure certain features that will customize DISCOVERY WORKBENCH for the needs of individual laboratories. Some of the features below are active only if enabled by a Configuration Manager.
- Users can be required to log in to the software using their Windows credentials before accessing the program. User access can be limited to a subset of valid Windows users.
- An audit log window tracks events, changes, and file imports.
- Data integrity tools add digital signatures to all exported text data files and quarantined data files from previously used plates and plates with unreadable or expired barcodes.
- A hide feature can be used to reduce the number of visible data sets, experiments, templates, and plate layouts.
- Kit layouts can assign assay locations automatically for multiplex plates and allow users to disable data acquisition from specified spots.
- A plate summary section and status window appears during plate runs.
- Easy sample import allows sample names to be added to a plate layout by copying/pasting or importing from a spreadsheet or text file.
- An experiment data table contains data for all plates, assays, and samples in the experiment.
- New experiment inter-plate statistics automatically calculate statistics for replicates across plates.
- QC acceptance criteria can automatically identify and flag samples for follow-up based on user-defined parameters for an experiment, assay, sample type, or individual sample.
- New curve fitting profiles include five parameter, linear, log-log, and exponential curve fitting.

Enhancements

- Data analysis software is compatible with multiple Microsoft operating systems: Windows 7, Vista, and XP Professional Edition.
- Plate data from different plate types (e.g., 96-well, single-spot and 96-well, 4-spot) can be analyzed in the same experiment.
- Users can create plate layouts on non-instrument computers without specifying an instrument type.
- Users can control the number of significant figures displayed in data tables.
- Data table customization improvements allow users to save default table profiles.
- Data tables can include additional information such as adjusted signal, plate name, sample description, and algorithm parameters.
- An excluded samples table displays samples that have been manually excluded from analysis.
- Plate data history can be saved to a text file or PDF for off-line viewing.
- Plate data history was expanded to include tabs for templates and plate layouts.
- A creation time column was added to the experiment tab in the plate data history.
- Cutting and pasting tables with merged cells into other programs has been improved.
- Experiment files can be exported from DISCOVERY WORKBENCH 4.0 into earlier software versions. Export experiments as Discovery Workbench 3.0 (*.zip) files to allow for import into DISCOVERY WORKBENCH 3.0.

MSD[®] DISCOVERY WORKBENCH 4.0 RELEASE NOTES

Fixes

- Corrected issues with using sample and experiment names with >10 numeric characters.
- Corrected issues with creating large PDF reports.
- Corrected other miscellaneous bugs.

System Requirements

Component	Minimum	Recommended
Operating System	Windows XP, Professional Edition	Windows 7, 32 bit, SECTOR [®] Imagers
		Windows 7, 64 bit, MESO QuickPlex SQ 120
		Windows 7, 32 or 64 bit, desktop version
Processor	INTEL PENTIUM 4, 1 GHz	INTEL Core 2 Duo, 2.1 GHz or higher
RAM	1 GB	2 GB or greater
Hard Disk Space	40 GB	160 GB or greater
CD-ROM/DVD-ROM Drive	DVD-ROM	16X CD/DVD+/-RW
Screen Resolution	1024 x 768	1280 x 1024
Screen Colors	256 colors	24-bit True Color
Other Software	Adobe Reader or PDF Viewing software	Adobe Reader or PDF Viewing software

Limitations

Plate layouts created in the desktop version of DISCOVERY WORKBENCH 4.0 cannot be imported into DISCOVERY WORKBENCH 4.0 for instrument computers. Layouts created on the instrument computer can be imported into DISCOVERY WORKBENCH 4.0 used on a desktop.

The desktop version of DISCOVERY WORKBENCH 4.0 is compatible with Microsoft operating systems in English, French, German, and Japanese. Operating systems in other languages are not supported.

Technical Support

Contact a technical representative at ScientificSupport@mesoscale.com for more information.

MESO SCALE DISCOVERY, MESO SCALE DIAGNOSTICS, MSD, DISCOVERY WORKBENCH, MULTI-ARRAY, MULTI-SPOT, QUICKPLEX, SECTOR, SECTOR PR, SECTOR HTS, SULFO-TAG, S-PLEX, V-PLEX, STREPTAVIDIN GOLD, MESO, www.mesoscale.com, SMALL SPOT (design), 96 WELL 1, 4, 7, & 10-SPOT (designs), 384 WELL 1 & 4-SPOT (designs), MSD (design), S-PLEX (design), V-PLEX (design), and SPOT THE DIFFERENCE are trademarks and/or service marks of Meso Scale Diagnostics, LLC.

©2013, 2014 Meso Scale Diagnostics, LLC. All rights reserved.