

Discovery Workbench[®] Data Analysis Tools (v3.0) Software CD

Overview: The MSD Discovery Workbench Data Analysis Tools (v3.0) software CD contains the full program for analysis of data collected by MSD SECTOR™ instruments. Upon installation, this software will function for 30 days without a permanent software seat license. After this demonstration period, the activation of a permanent software seat license is required. Software licenses are available and can be purchased from MSD before or after the end of the demonstration period. The process of activating a software license is described below.

Appropriate Use: *The MSD Discovery Workbench Data Analysis Tools (v3.0) software may only be used on a computer that is not attached to a MSD SECTOR instrument. Do not insert the MSD Discovery Workbench Data Analysis Tools (v3.0) disk into a computer that is connected to a MSD SECTOR instrument.*

Operating System Requirements: The software is compatible with the English language version of Windows 2000 or Windows XP operating systems.

Installation Instructions: Software installation occurs through a standard installer. Administrator level privileges are needed to install the software.

1. Insert the software disk into the CD drive.
2. The “auto run” feature will begin installing the software. (If the installer does not start, use the explorer to navigate to the CD and double click on the setup file.)
3. Enter the user information in the dialog box when prompted.
4. Select the default option “Install this application for anyone”.
5. Click “next” to proceed with installation.
6. When the installation wizard reports “installation complete”, the user has two options.
 - a. Reboot the machine now.
 - b. Reboot the machine later.The default option is “yes, I want to restart my computer now”. The computer must be restarted before using the software.

Installation note: One disk can be used to install software on multiple computers. However, each computer will require a computer specific license code to continue use of the software after the 30 day demo period.

License Activation Instructions: Upon each use of the software in demonstration mode, users will have an option to enter to enter a permanent software license code. To permanently activate a seat license, follow the instructions below.

1. Click on the “activate” icon on the Discovery Workbench start up licensing screen.
2. The system will generate “User Code 1” and “User Code 2”.
3. Report these user codes to MSD Customer Service by phone to +1 (301) 947-2025 (Maryland, USA) or email to CustomerSupport@Mesoscale.com. Also, provide your seat license purchase order number or MSD Sales Order confirmation number.
4. MSD will respond by providing your permanent seat license code. Enter this code into the field for “Activation Code 1”. Leave the field for “Activation Code 2” empty.
5. Click “continue” and the software will confirm successful activation.

Usage Note: The software CD contains the software user’s manual and training exercises.

