

Using Database Scripts to Install or Update the Microsoft® MSDE or Microsoft SQL Server™ Screening Database SR2 in Discovery-1 v6.1

Document ID
D50016

Product
Discovery-1, version 6.1

Created
30-May-2003

Last Reviewed
3-Mar-2004

Abstract

This document describes how to use the provided database installation scripts to install or upgrade your Microsoft SQL Server Desktop Engine (MSDE) or Microsoft SQL Discovery-1 screening database to Service Release 2 (SR2). The included scripts enable you to install the database for the first time, upgrade your existing database to SR2, or install additional databases to your system.

Notes

- ◆ This document assumes that Discovery-1 is installed with the included version of MSDE or that you already have a full version Microsoft SQL Server database server installed and running.
- ◆ The version requirement for MSDE is SQL Server 7.0 Service Pack 3 or above. The version requirement for Microsoft SQL is SQL Server 2000 Service Pack 1 or above.
- ◆ The installer must have Administrative privileges on the computer running the database server to use the scripts.
- ◆ Database names cannot contain spaces or special characters. The names screening 1 and screening 2 are not valid database names. The names screening_1 and screening_2 are valid database names. Limit the characters used for the database names to alphanumeric (0-9, a-z) and underscores (_).
- ◆ The scripts must be run on the computer that is running the database server. You cannot do a network installation of the scripts.
- ◆ The computer that is running the database server does not need to have Discovery-1 or MetaMorph software installed.
- ◆ The Scripts needed to perform the update are located on the Universal Imaging Corporation FTP site at: <ftp://www.universal-imaging.com/pub/uic/screening/scripts/microsoft/>

Instructions

Installing the Screening Database SR2

There are two stages in installing the database. In the first stage you run a script and answer questions about the database. The script generates a command that installs the database using the information you provided. Use the following procedure to generate the command and install the database:

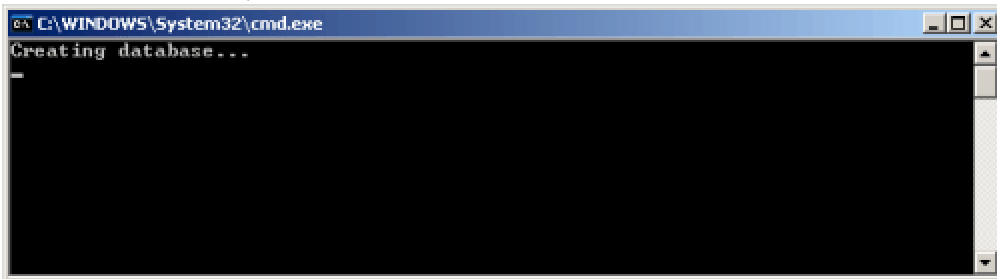
1. Open a web browser or ftp program and go to: <ftp://www.universal-imaging.com/pub/uic/screening/scripts/microsoft/>.
2. Copy the sr2 folder to your local computer running your database.
3. Using Windows Explorer, open the folder sr2 on your local computer.
4. Double-click generate-scripts.wsf. The Path dialog box opens.

Keywords: database msde microsoft v6.1 v6

Issue Type: dbs database admn

Discovery-1 Online Support

5. Select the default path in the text box and replace it with path of the folder that will contain the database.
6. Click *OK*. The database Name dialog box opens.
7. Type the name of the database in the text box.
8. Click *OK*. The username dialog box opens.
9. Type a user name in the text box.
10. Click *OK*. The Password dialog box opens.
11. Type a password for the user you created in Step 9.
Note: The username and password entered here have no connection with user names or passwords set up using the Meta Imaging Series Administrator.
12. Click *OK*. The password dialog box closes and the following three files are created in the sr2 folder:
 - o install-sr2.cmd
 - o create-sr2-job.sql
 - o create-sr2-schema.sql
13. Double-click `install-sr2.cmd`. An MS-DOS command prompt will open, run the command, and install the database, as shown below:



The command prompt will close when the installation is complete. No user interaction is required. If the install was not successful, read the `create-sr2-schema.out` file located in the sr2 directory. This is a text file containing pertinent information about the installation.

Note: After installation, you can safely delete the sr2 folder from your hard drive, or save it for future database installations.

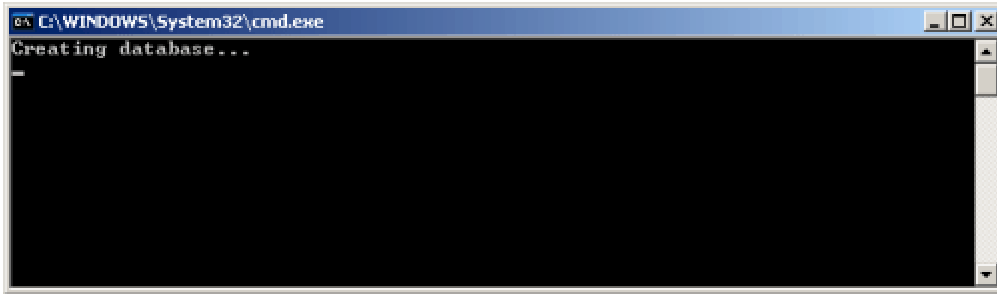
Updating the Screening Database to SR2

Use the following procedure to update the database to SR2:

1. Open a web browser or ftp program and go to: <ftp://www.universal-imaging.com/pub/uic/screening/scripts/microsoft/>.
2. Copy the sr1 to sr2 upgrade folder to the computer running your database.
3. Using Windows Explorer, open the folder sr1 to sr2 upgrade on your local computer.
4. Double-click `generate-scripts.wsf`. The oldDatabaseName dialog box opens.
5. Enter the current name of the database in the text box.
6. Click *OK*. The newDatabaseName dialog box opens.
7. Type either a new name or use the current name in the text box.
8. Click *OK*. The password dialog box closes and the following files are created in the sr2 folder:
 - o upgrade-sr1-to-sr2.cmd
 - o upgrade-sr1-to-sr2.sql

Discovery-1 Online Support

9. Double-click `upgrade-sr1-to-sr2.cmd`. An MS-DOS command prompt will open, run the command, and update the database, as shown below:



The command prompt will close when the update is complete. No user interaction is required. If the install was not successful, read the `upgrade-sr1-to-sr2.out` file located in the `sr1 to sr2` upgrade directory. This is a text file containing pertinent information about the upgrade.