

Using Database Scripts to Install or Update the Microsoft® MSDE or Microsoft SQL Server™ Screening Database SR3 for Discovery-1, Version 6.2

Document ID
D50017

Product
Discovery-1, version 6.2

Created
7-May-2004

Last Reviewed
15-Jun-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 3 (SR3). The included scripts enable you to install the database for the first time, upgrade your existing database to SR3, or install additional databases on your system.

Notes

- ♦ Version 6.2 of Discovery-1 requires the use of Discovery-1 Screening Database Service Release 3 (SR3). You cannot use SR2 with version 6.2 of Discovery-1.
- ♦ The Discovery-1 SR3 database can be run on the following three platforms:
 - ♦ MSDE SQL Server 7.0 Service Pack 3 or above
 - ♦ Microsoft SQL Server 2000 Service Pack 1 or above
 - ♦ Oracle 8i/9i Database Server
- ♦ This document does not cover the Oracle installation or upgrade procedure option. For information on installing the Oracle server, contact Universal Imaging Corporation Technical Support for assistance.
- ♦ 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 SR3

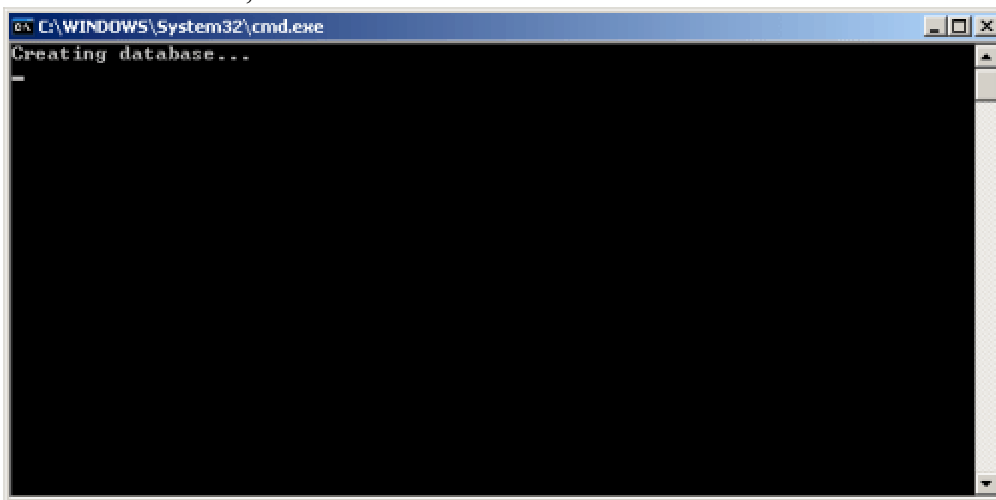
There are two stages in installing the database. In the first stage you run a script and answer questions about the database. The script generates a file that installs the database using the information you provided. Use the following procedure to generate the file and install the database:

1. Open a web browser or ftp program and go to: <ftp://www.universal-imaging.com/pub/uic/screening/scripts/microsoft/> .

Keywords: database msde microsoft v6.2 v6
Issue Type: dbs database admn

Discovery-1 Online Support

2. Copy the `sr3` folder to the local computer running your database.
3. Using Windows Explorer, open the folder `sr3` on your local computer.
4. Double-click `generate-scripts.wsf`. The Path dialog box opens.
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 `sr3` folder:
 - o `install-sr3.cmd`
 - o `create-sr3-job.sql`
 - o `create-sr3-schema.sql`
13. Double-click `install-sr3.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-sr3-schema.out` file located in the `sr3` directory. This is a text file containing pertinent information about the installation.

Note: After installation, you must configure the location of screening images and set the auto run timeout using the Database Utilities command in the Meta Imaging Series Administrator. You will not be able to save images to your database until this step is complete. For more information about the Database Utilities command, refer to the online help in the Meta Imaging Series Administrator.

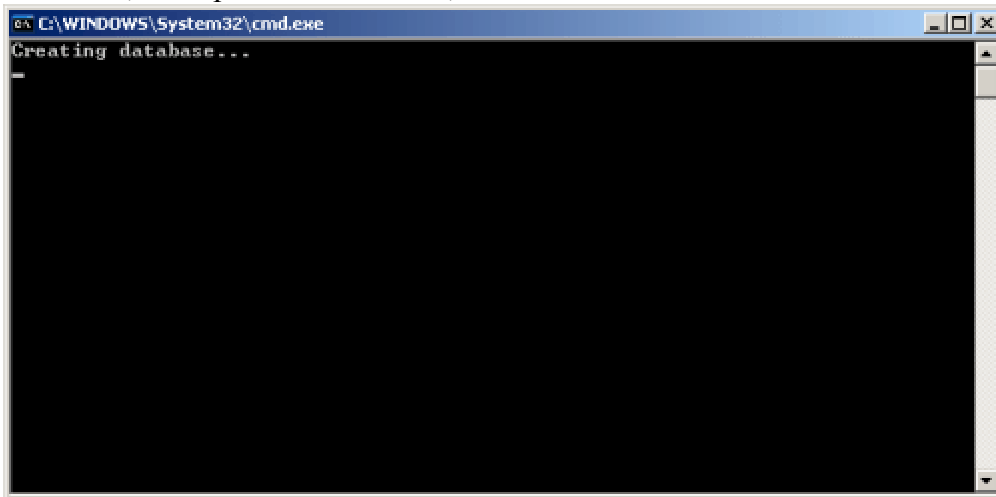
Updating a Screening Database to SR3

Use the following procedure to update a SR2 database to SR3:

1. Open a web browser or ftp program and go to: <ftp://www.universal-imaging.com/pub/uic/screening/scripts/microsoft/>.

Discovery-1 Online Support

2. Copy the `sr2 to sr3 upgrade` folder to the computer running your database.
3. Using Windows Explorer, open the folder `sr2 to sr3 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 password dialog box closes and the following files are created in the `sr2` folder:
 - o `upgrade-sr2-to-sr3.cmd`
 - o `upgrade-sr2-to-sr3.sql`
7. Double-click `upgrade-sr2-to-sr3.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-sr2-to-sr3.out` file located in the `sr2 to sr3 upgrade` directory. This is a text file containing pertinent information about the upgrade.

Note: After installation, you must configure the location of screening images and set the auto run timeout using the Database Utilities command in the Meta Imaging Series Administrator. You will not be able to save images to your database until this step is complete. For more information about the Database Utilities command, refer to the online help in the Meta Imaging Series Administrator.