

Installing Microsoft SQL Server Desktop Engine Service Pack 3

Abstract

This tech note describes the recommended way to install the Microsoft SQL Server Desktop Engine Service Pack 3. Installing this version of MSDE will protect against the SQL "slammer" worm.

Instructions

Recommended Procedure:

MetaMorph and MetaVue, versions 6.0 and later, use the Microsoft SQL Server Desktop Engine (MSDE) to manage the Image Browser thumbnail database. Microsoft released MSDE Service Pack 3 to solve the security hole exposed by the "Slammer Worm". Universal Imaging recommends the use of MSDE with Service Pack 3.

The Meta Imaging Series 6.0 CD installs the MSDE with Service Pack 1. For users with this version, we recommend you uninstall MSDE and then re-install using Service Pack 3. The following are the directions on how to accomplish this.

Please note that if you are running Meta Imaging Series applications earlier than version 6.0, you did not have MSDE installed on your system by the Meta Imaging Series installation CD.

Step 1: Determine the Version of SQL Server Installed on your Computer.

This can be done by following these instructions:

In the "Add or Remove Programs" Control Panel, look for "Microsoft SQL Server Desktop Engine."

If it is not present, MSDE is not in use -- consult with a person who is responsible for any SQL Server that is in use.

If "Microsoft SQL Server Desktop Engine" is present, highlight it. The words "Click here for support information" appear. Click on those words to display the "Support Info" dialog.

In that dialog, the Version information should be one of the following:

8.00.194 -- MSDE as Released To Manufacturing, also known as the RTM version. No Service Pack has been applied;

8.00.384 -- MSDE with Service Pack 1 (SP1);

8.00.534 -- MSDE with Service Pack 2 (SP2);

8.00.760 -- MSDE with Service Pack 3 (SP3).

Document ID
T20018

Product
MetaMorph
MetaVue
Discovery-1

Platform
Windows 2000

Created
28-Jan-2003

Last Reviewed
31-Jan-2003

Keywords: msde, sql,slammer, worm, virus, security, server

Issue Type: Application

If you have determined that your system is not already running SQL MSDE SP3, then you can use the following instructions to get MSDE SP3 up and running correctly on your system.

Uninstall the Microsoft SQL Server Desktop Engine by pressing the Remove button in Add or Remove Programs dialog if one of the previous versions exists on your computer.

Reboot the machine, only necessary after the previous uninstall.

Go to the link below and download the latest Service Pack (3) for SQL Server Desktop Engine. Specifically, the file you need from this site is SQL2KDeskSP3.exe

SQL Server 2000 SP3 Link:

<http://www.microsoft.com/downloads/details.aspx?FamilyId=9032F608-160A-4537-A2B6-4CB265B80766&displaylang=en#filelist>

Unpack the SQL2KDeskSP3.exe file. Choose the default directory, c:\sql2ksp3.

Using NotePad, add the following lines to the c:\sql2ksp3\MSDE\Setup.ini file:

```
SECURITYMODE=SQL  
INSTANCENAME=MSSQLSERVER  
BLANKSAPWD=1
```

Run the Setup.exe found in the c:\sql2ksp3 directory. This will install the Microsoft SQL Server Desktop Engine SP3.

In the Control Panel, go to Administrative Tools, Services and start the MSSQLSERVER and make sure the Startup Type is Automatic.

You can delete the files that were unpacked to c:\sql2ksp3

You can verify that you have the correct version of the Microsoft SQL Server Desktop Engine by viewing its Support Information in the Add or Remove Programs. Click on Microsoft SQL Server Desktop Engine -- the words "Click here for support information" appear. Click on those words to display the "Support Info" dialog.

In that dialog, the Version information should be MSDE SP3.