

# VWorks Automation Control 13.1.7

## Setup Guide Addendum

*For MySQL 8.0.18*

This guide contains the following topics:

- [“About this addendum” on page 2](#)
- [“Installing MySQL” on page 2](#)
- [“Installing and setting up a MySQL ODBC” on page 14](#)
- [“Connecting to the Inventory database” on page 18](#)

## About this addendum

This is an addendum to the VWorks 13 edition of the [VWorks Automation Control Setup Guide](#) (part number G5415-90067). This addendum is applicable if your system meets both of the following conditions:

- You are running VWorks 13.1.7 on Microsoft Windows 10 64-bit.
- Your system contains a device, such as a Labware MiniHub, that requires a database for the VWorks labware inventory management feature.

*Note:* The database may be configured already if the computer was provided by Agilent.

This addendum describes how to do the following:

- Install MySQL 8.0.18.
- Install and set up a MySQL 8.0.18 ODBC.
- Connect to the Inventory database for either of the following:
  - MySQL 5.7
  - MySQL 8.0.18

## Installing MySQL

### Before you start

#### Required installation files

Ensure you have the following installation files:

- VC\_redist.x64.exe
- VC\_redist.x86.exe
- mysql-installer-community-8.0.18.0.msi
- vworks\_mysql\_8.0.sql

This file is installed with the VWorks software in the following location:

C:\Program Files (x86)\Agilent  
Technologies\VWorks\Schema\vworks\_mysql\_8.0.sql

For assistance, contact Agilent Technical Support.

#### Required computer settings

Before you install the software:

- Log on to the computer as an administrator with full rights.
- In Microsoft Windows Explorer, select the following **View > Show** settings:
  - Select **Hidden items**.
  - Select **File name extensions**.

For detailed instructions, see the Microsoft Windows user documentation.

## Installation and setup workflow

Step	For this task...	See...
1	Uninstall any existing MySQL versions, and then restart the computer.	"Uninstalling previous versions of MySQL" on page 3
2	Install MySQL Server 8.0.18	"Installing MySQL 8.0.18" on page 6
3	Install MySQL ODBC and set up the data sources.	"Installing and setting up a MySQL ODBC" on page 14
4	Connect to the VWorks database.	<i>VWorks Automation Control Setup Guide</i> for VWorks 13.

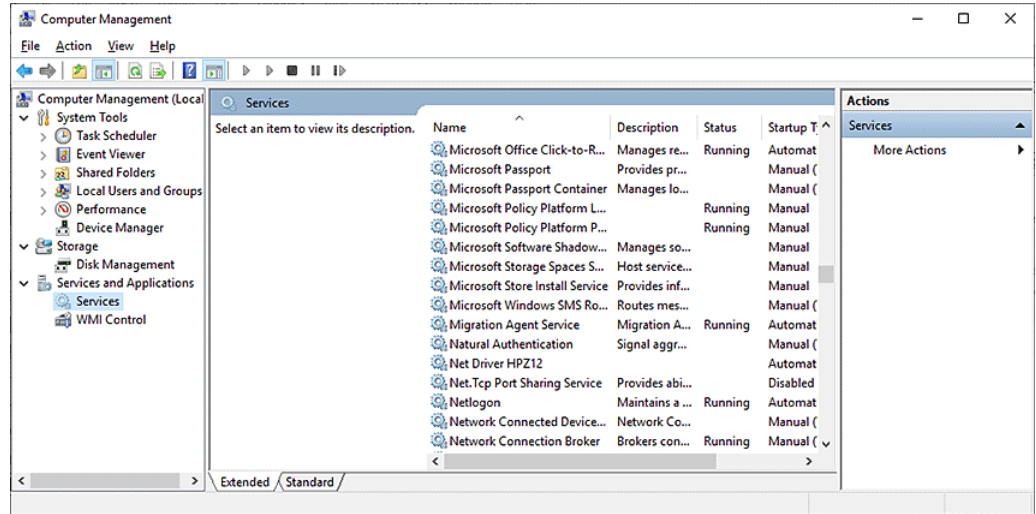
## Uninstalling previous versions of MySQL

### **Before you uninstall:**

Make a backup copy of your existing database before uninstalling it. To back up the Inventory database (velocity11), labware definitions, liquid classes, and pipetting techniques, see the [VWorks Automation Control User Guide](#).

### **To uninstall MySQL:**

- 1 Log on to the computer as an administrator.
- 2 In the Microsoft Windows desktop, go to **Settings > Apps & features**, and uninstall the following:
  - a MySQL Server
  - b MySQL Connector/ODBC
- 3 Restart the computer.
- 4 Verify that the MySQL service is completely removed:
  - a In Microsoft Windows Explorer, right-click the **This PC** icon. In the shortcut menu, click **Manage**. The Computer Management window opens.
  - b In the **Computer Management** window, click **Services and Applications > Services**.
  - c In the list of **Services**, ensure MySQL does not appear:
    - If MySQL does not appear, go to [step 8](#).
    - If MySQL appears in the list, continue to step 5.



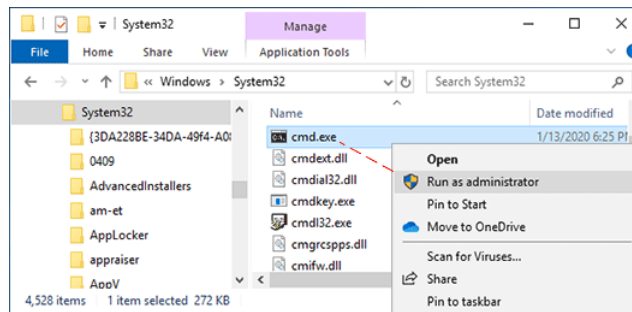
5 In the list of **Services**, right-click **MySQL** and click **Stop**.

*Note:* If an error message appears and states that it is unable to stop, ignore the message.

6 Open the Command Prompt window, as follows:

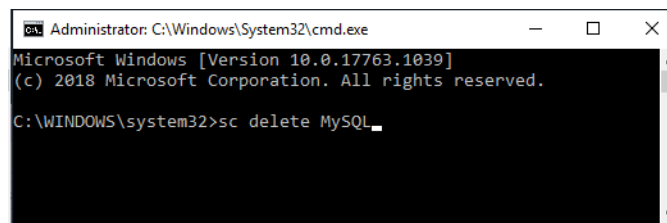
a In the **C:\Windows\System32** folder, right-click **cmd.exe**, and then select **Run as administrator**.

b Click **Yes** in the **User Account Control** dialog box that opens.



c Execute the following line in the Command Prompt window:

```
sc delete MySQL <enter>
```

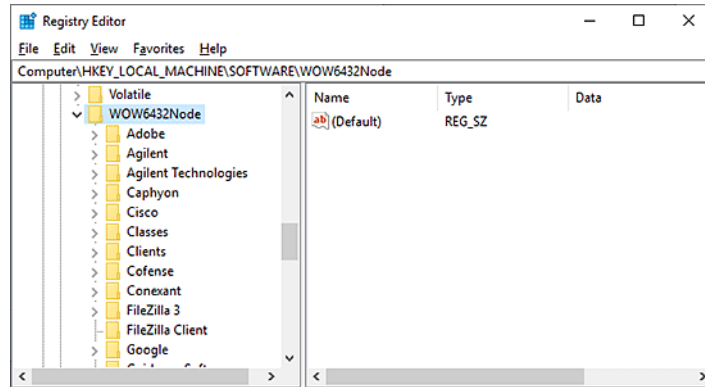


7 Check the registry keys for any references to old versions of MySQL as follows:

a Open the Registry Editor.

b Navigate to HKEY\_LOCAL\_MACHINE\SOFTWARE\WOW6432Node\MySQL AB\

c If present, delete MySQL Server 5.1 or MySQL Server 5.0.



8 Restart the computer.

### IMPORTANT

Ensure that you restart the computer after uninstalling the MySQL Server and MySQL Connector/ODBC.

## Installing MySQL 8.0.18

Before installing MySQL 8.0.18, install the Microsoft Visual C++ redistributable package.

### To install the Microsoft Visual C++ redistributable package:

- 1 Run the vc\_redist.x64.exe file.
- 2 Follow the Setup Wizard instructions.
- 3 Run the vc\_redist.x86.exe file.
- 4 Follow the Setup Wizard instructions.

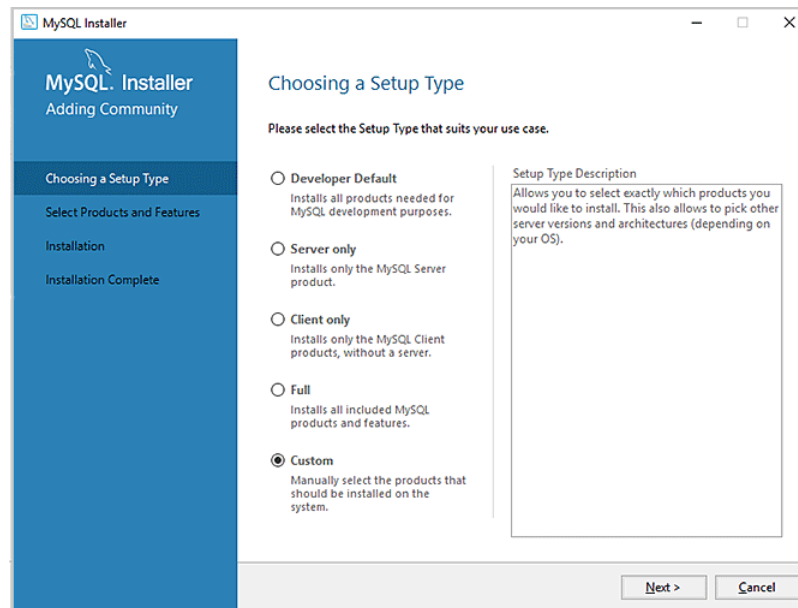
Note: If the computer already has Microsoft Visual Studio 2017 or Microsoft Visual C++ 2017 Redistributable installed, the MySQL installer might not recognize that the VC 2015 Redistributable is installed even if it is. To fix this issue, go to Settings > Apps. In the Apps & features window, locate Microsoft Visual C++ 2015, right-click, and then choose Modify. Click Repair when the installer window opens.

### To install MySQL 8.0.18:

- 1 Run the mysql-installer-community-8.0.18.0.msi. The MySQL Installer setup wizard opens.

Note: If Visual Studio 2017 or the Visual C++ 2017 redistributable package is installed on the computer, the MySQL installer might display an error message stating that Visual C++ redistributable is not installed even if it is. To fix this, go to Apps and Features, select Microsoft C++ Redistributable, right-click and then select Modify. In the Microsoft C++ Redistributable Setup screen, click Repair.

- 2 In the **Choosing a Setup Type** page, select **Custom**, and then click **Next**.

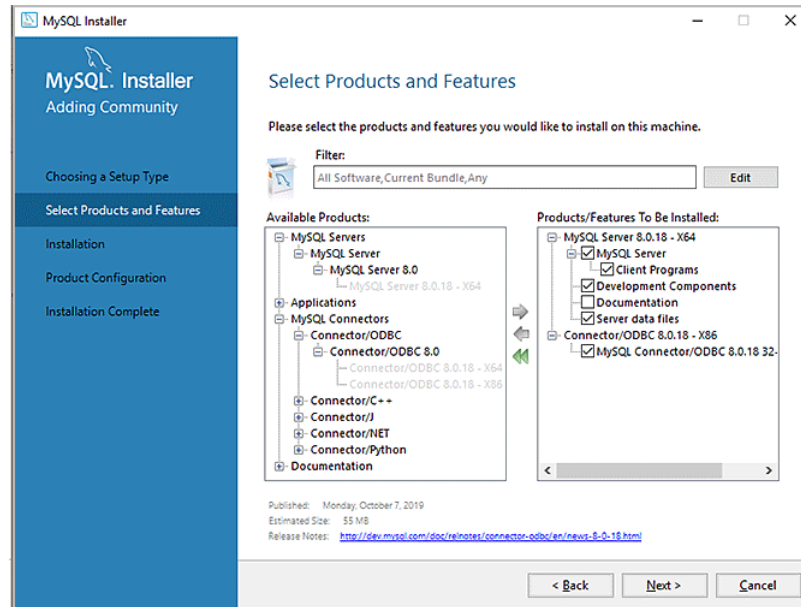


- 3 In the **Select Products and Features** page, select the following **Available Products** and then click the right arrow to move each selection to **Products/Features To Be Installed**:

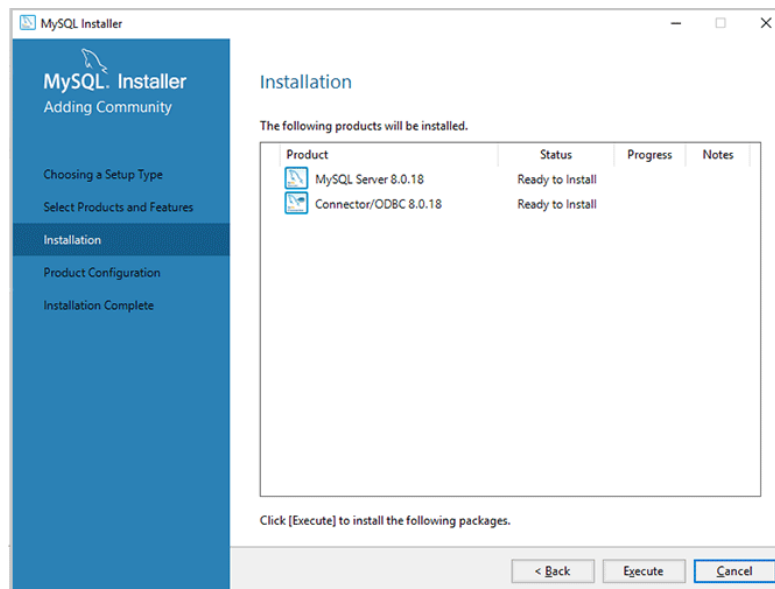
**MySQL Servers > MySQL Server > MySQL Server 8.0 > MySQL Server 8.0.0.1.8- X64**

**MySQL Connectors > Connector/ODBC > Connector ODBC 8.0.18 -X86**

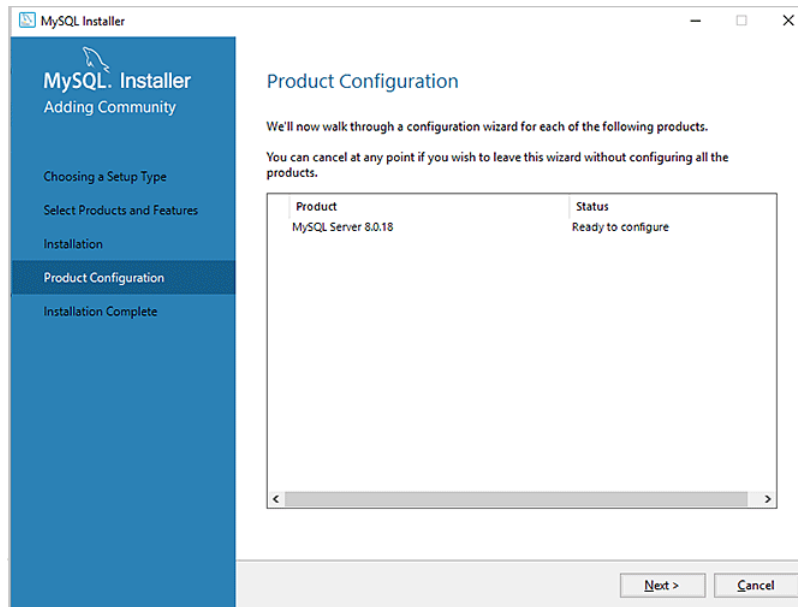
Click **Next**.



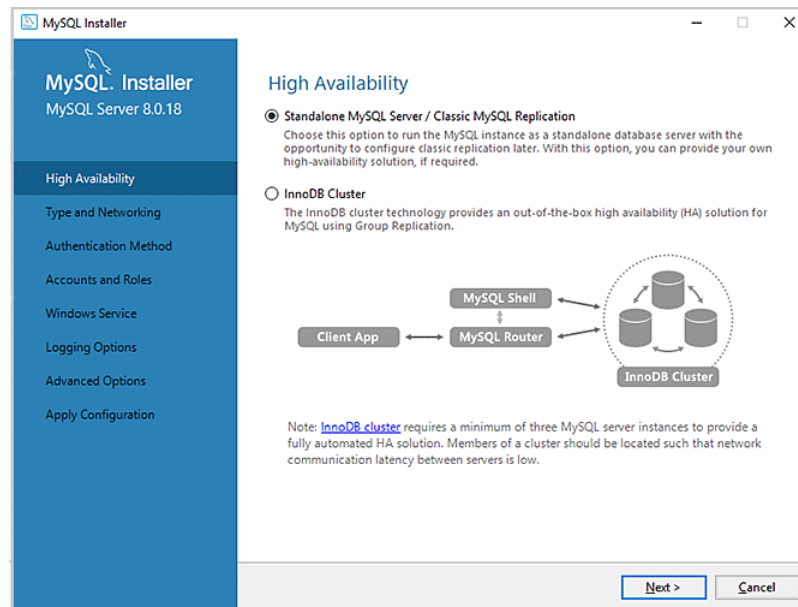
- 4 In the **Installation** page:
- Make sure that MySQL Server 8.0.18 and Connector/ODBC 8.0.18 are **Ready to Install**, and then click **Execute**.
  - When the installation is finished, click **Next**.



5 In the **Product Configuration** page, click **Next**.



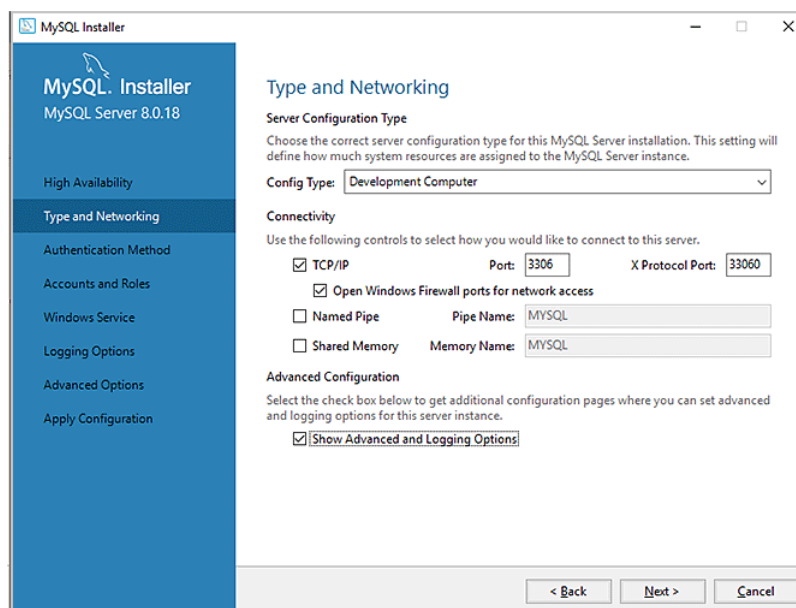
6 In the **High Availability** page, select **Standalone MySQL Server / Classic MySQL Replication**, and then click **Next**.





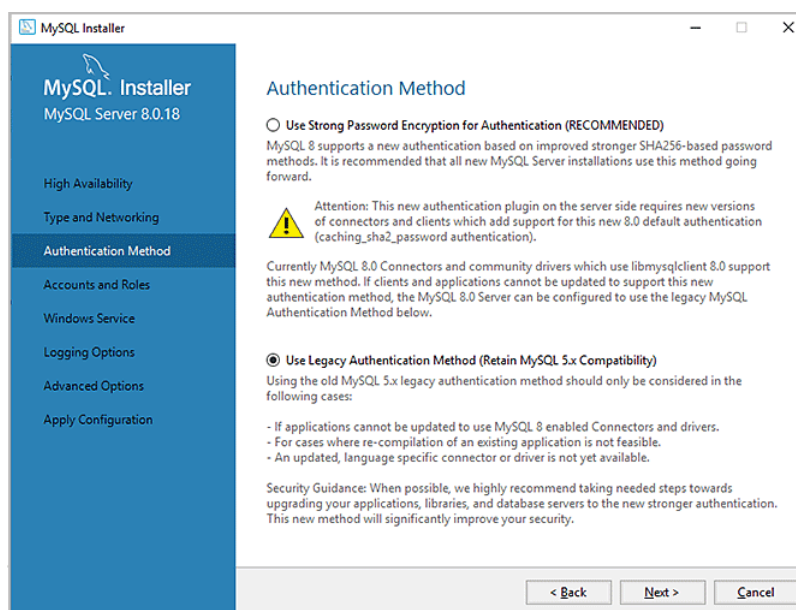
- 7 In the **Type and Networking** page, specify the following:
- **Config Type > Development Computer**
  - **TCIP/IP > Open Windows Firewall ports for network access**
  - **Show Advanced and Logging Options**

Click **Next**.

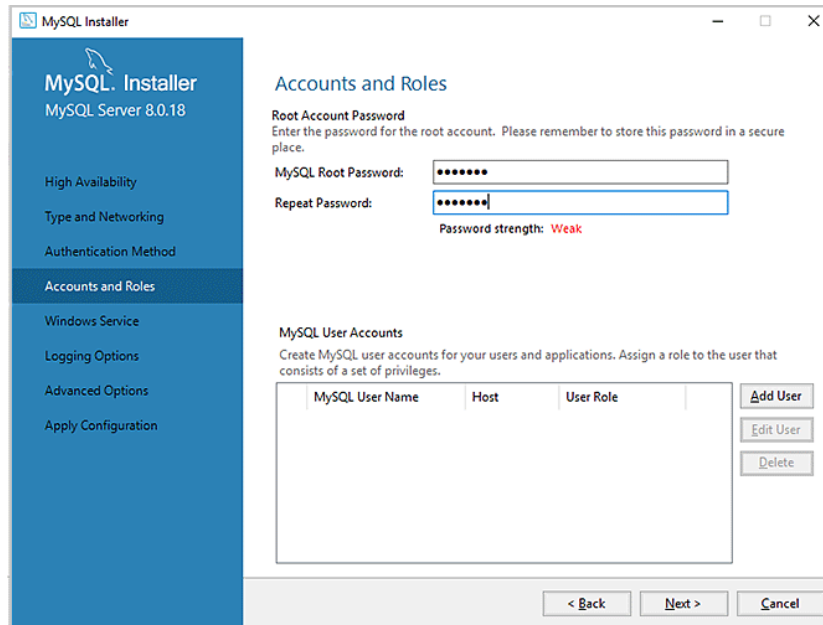


- 8 In the **Authentication Method** page, select **Use Legacy Authentication Method (Retain MySQL 5.x Compatibility)**.

Click **Next**.

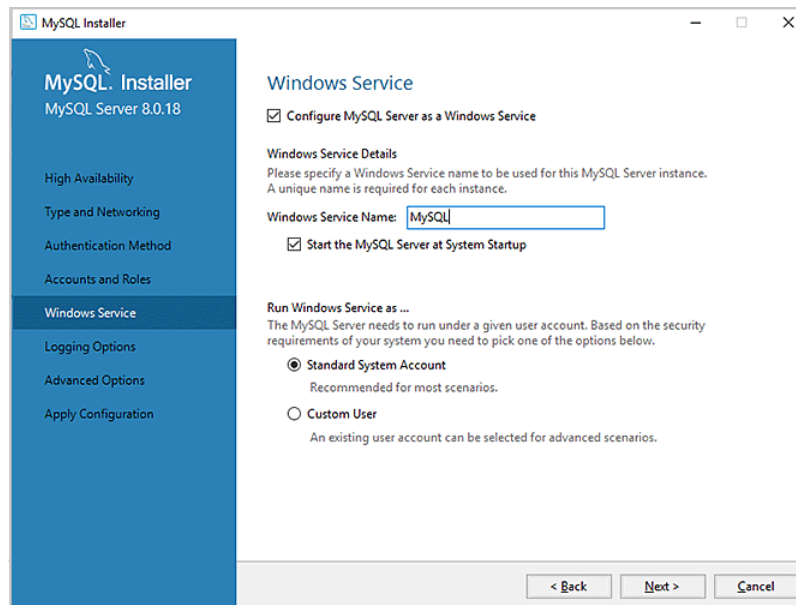


- 9 In the **Accounts and Roles** page, type `agilent` for **MySQL Root Password** and **Repeat Password** fields.  
Click **Next**.

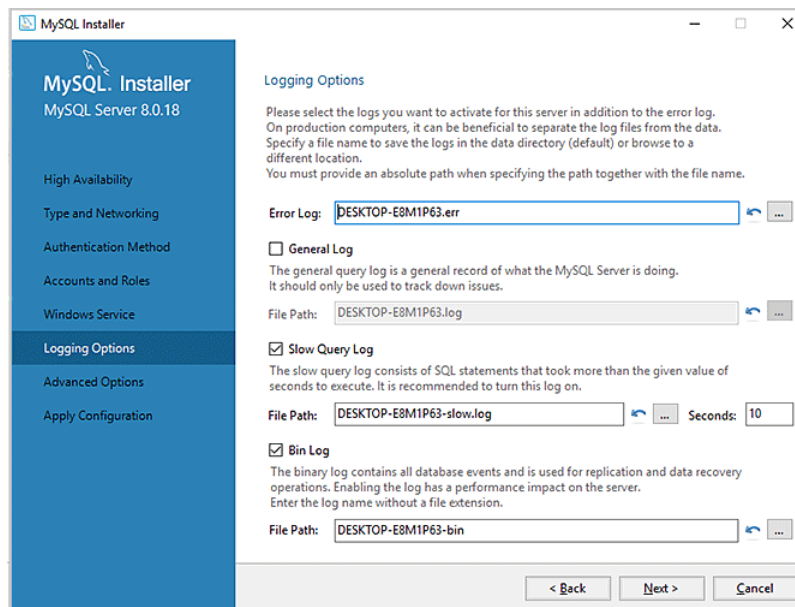


- 10 In the **Windows Service** page, specify the following:
- **Configure MySQL Server as a Windows Service**
  - Make sure you specify the following:
    - **Windows Service Name** > `MySQL`
    - **Start the MySQL server at System Startup**
    - **Run Windows Service as** > **Standard System Account**

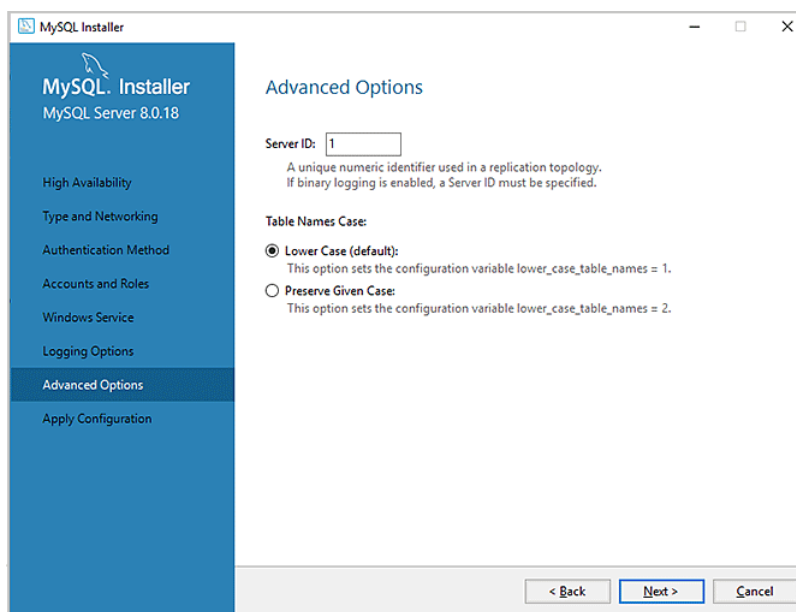
Click **Next**.



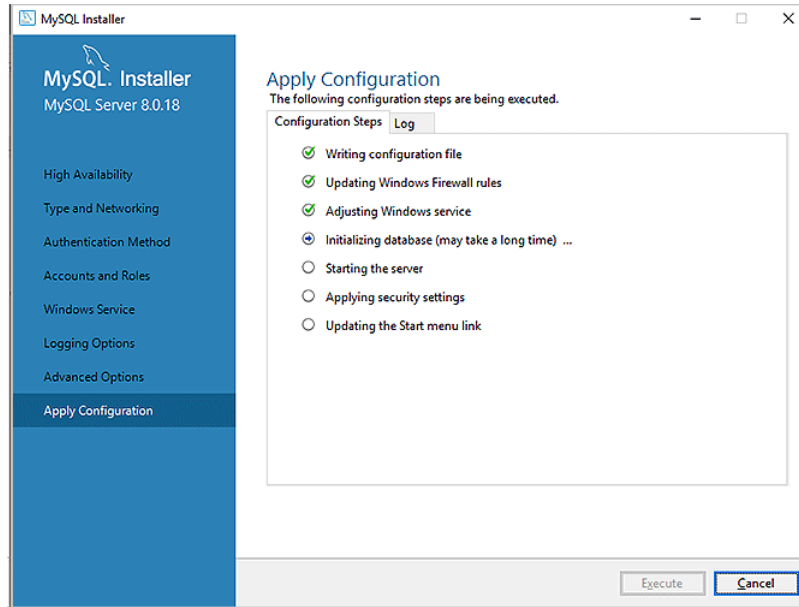
- 11 In the **Logging Options** page, accept the default settings. Click **Next**.  
Note: The file paths are based on the computer name.



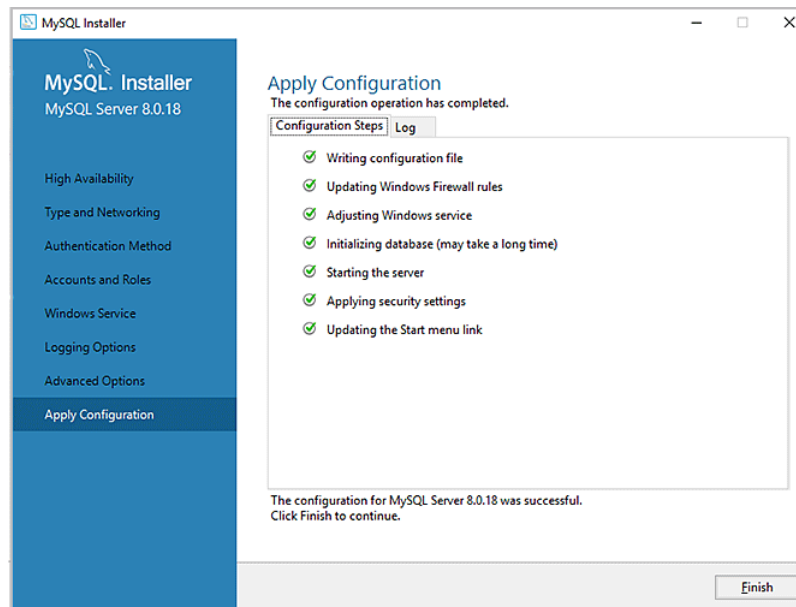
- 12 In the **Advanced Options** page, accept the default settings. Click **Next**.



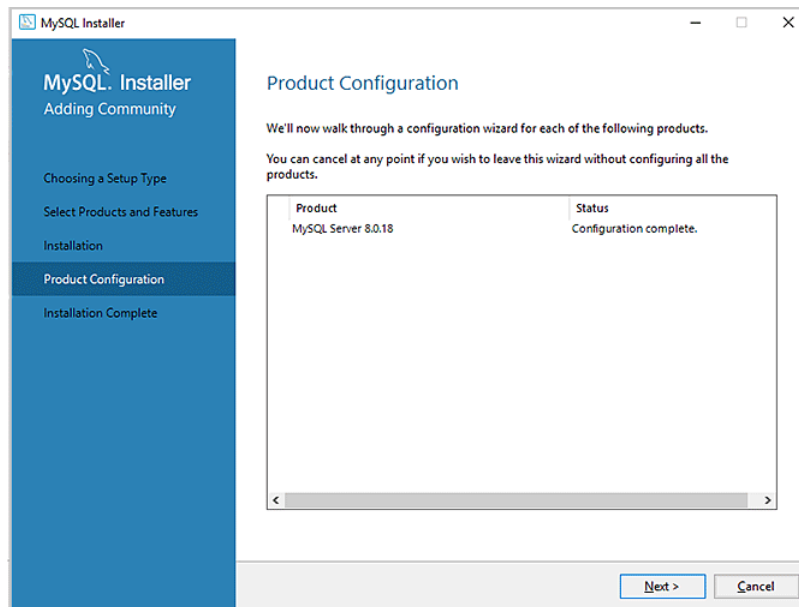
- 13 In the **Apply Configuration** page, click **Execute**.



14 When the screen displays the configuration was successful text, click **Finish**.



15 In the **Product Configuration** page, click **Next**.



16 Click **Finish** in the **Installation Complete** page.

# Installing and setting up a MySQL ODBC

## About this topic

This topic describes how to install and set up a data source in MySQL ODBC (open database connectivity) for the VWorks Inventory Editor.

During the MySQL ODBC connection setup you will:

- Import the VWorks database schema into MySQL. This sets up the database table structure for use by the VWorks Inventory Editor.
- Create, define, and configure new ODBC data source so that MySQL can access data in the database.

## Importing the database schema

### To import the database schema:

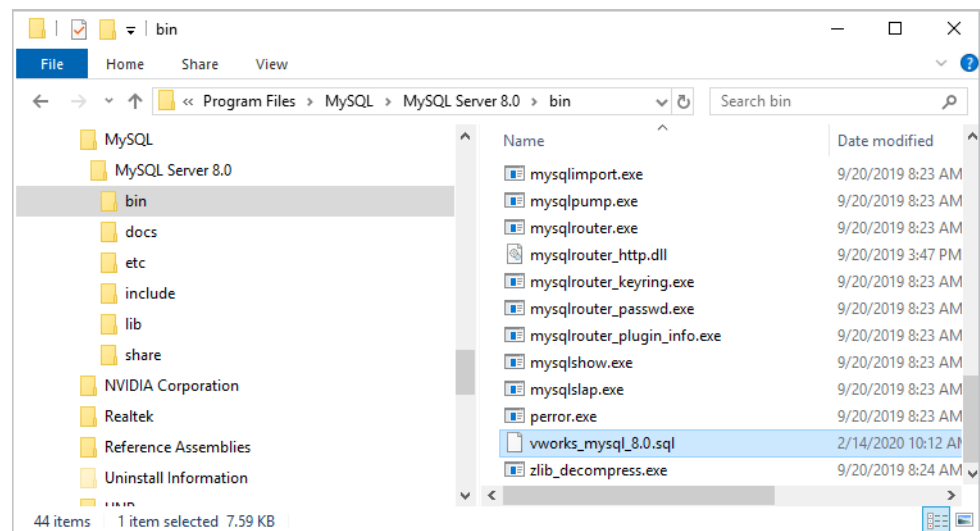
- 1 Copy the following database schema file:

vworks\_mysql\_8.0.sql

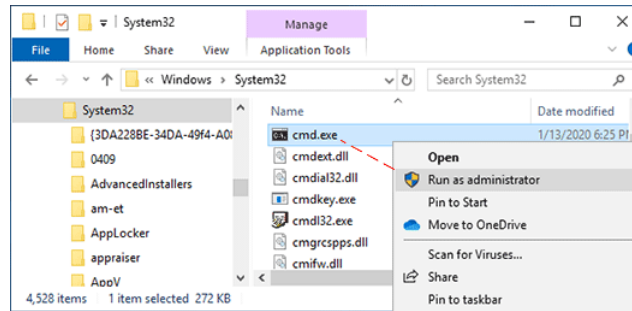
Note: This file is installed with the VWorks software in the folder:  
C:\Program Files (x86)\Agilent Technologies\VWorks\Schema\

Paste the copy into the following folder:

C:\Program Files\MySQL\MySQL Server 8.0\bin\



- 2 Open a Command Prompt window, as follows:
  - a In the C:\Windows\System32 folder, right-click **cmd.exe**, and then select **Run as administrator**.
  - b Click **Yes** in the **User Account Control** dialog box that opens.



- Execute the following lines in the Command Prompt window:
 

```
cd c:\program files\mysql\mysql server 8.0\bin <enter>
mysql -u root -p < vworks_mysql_8.0.sql <enter>
```

```
Administrator: C:\Windows\System32\cmd.exe - mysql -u root -p
Microsoft Windows [Version 10.0.17763.1039]
(c) 2018 Microsoft Corporation. All rights reserved.

C:\WINDOWS\system32>cd c:\program files\mysql\mysql server 8.0\bin

c:\Program Files\MySQL\MySQL Server 8.0\bin>mysql -u root -p < vworks_mysql_8.0.sql
Enter password: *****

c:\Program Files\MySQL\MySQL Server 8.0\bin>
```

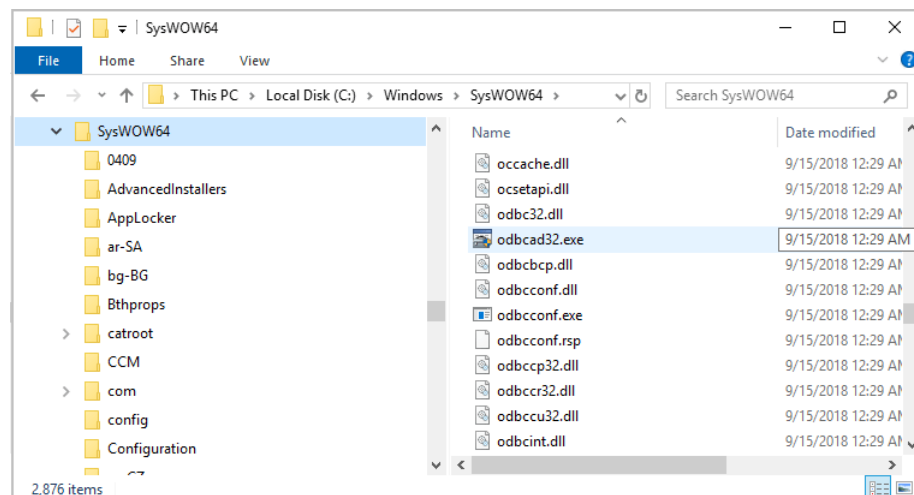
When prompted for password, enter `agilent`.

Make sure no errors occur while executing this command.

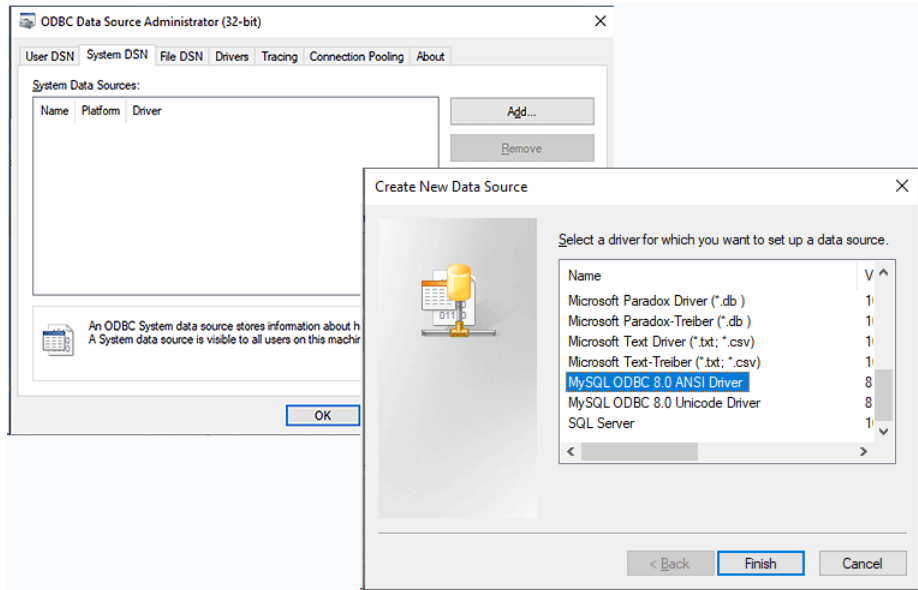
## Creating ODBC connections

### To create ODBC connections:

- Open the ODBC Data Source Administrator to create ODBC connections, as follows:  
In the `C:\Windows\SysWOW64` folder, right-click **odbcad32.exe**, and then select **Run as administrator**. The ODBC Data Source Administrator (64-bit) dialog box opens.



- 2 In the **ODBC Data Source Administrator** dialog box:
  - a Remove any previously defined MySQL ODBC Driver Sources as follows:
    - Click the **User DSN** tab, select any predefined MySQL ODBC Driver Sources named velocity11, and click **Remove**.
    - Click the **System DSN** tab, select any predefined MySQL ODBC Driver Sources named velocity11, and click **Remove**.
  - b In the **System DSN** tab, click **Add** to create a new data source.
  - c In the **Create New Data Source** dialog box, select **MySQL ODBC 8.0 ANSI Driver**, and then click **Finish**.



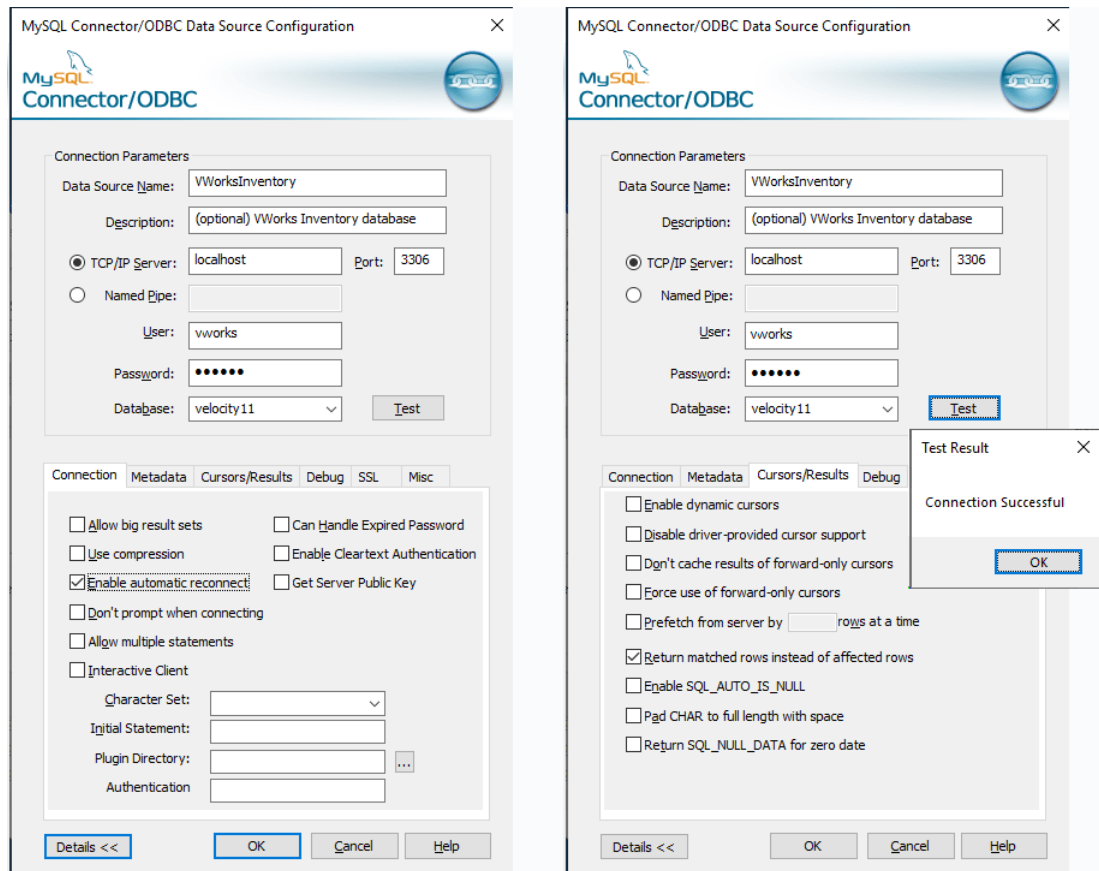
- 3 In the **MySQL Connector/ODBC Data Source Configuration** dialog box that opens, define your data source as follows:  
 Note: You do not have to provide a port number. Use the default number displayed.

Settings for VWorks Inventory database	
Connection Parameters	Settings
<b>Data Source Name</b>	VWorksInventory
<b>Description</b>	(optional) VWorksInventory database
<b>TCP/IP Server</b>	localhost
<b>User</b>	vworks
<b>Password</b>	vworks
<b>Database</b>	velocity11

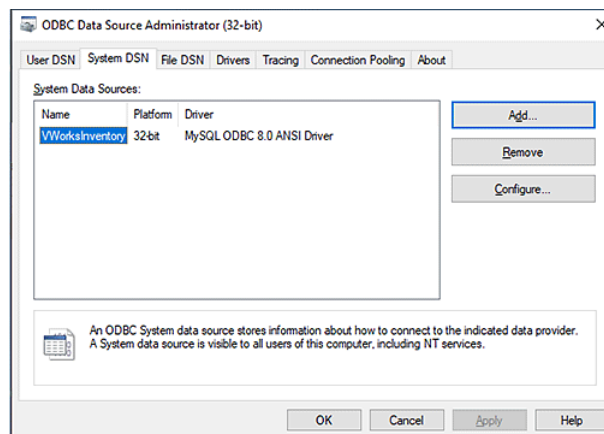
- a Click **Details** to expand the dialog box.
- b In the **Connection** tab, select the **Enable automatic reconnect** check box.
- c In the **Cursors/Results** tab, select the **Return matched rows instead of affected rows** check box.



4 Click **Test**. When the **Connection Successful** message appears, click **OK**.



5 In the **System DSN** tab, verify that the data source appears.



Click **OK** to close the dialog box.

For instructions on how to connect to the database in the VWorks software, see the [VWorks Automation Control Setup Guide](#) for VWorks 13.

# Connecting to the Inventory database

## Before you start

Before you can connect to the database, you must install and configure the inventory database. To install and configure the inventory database, see the appropriate instructions:

- For MySQL 8.0.18, see “Installing MySQL” on page 2 and “Installing and setting up a MySQL ODBC” on page 14.
- For MySQL 5.7, see the VWorks 13 edition of the *VWorks Automation Control Setup Guide* (part number G5415-90067).

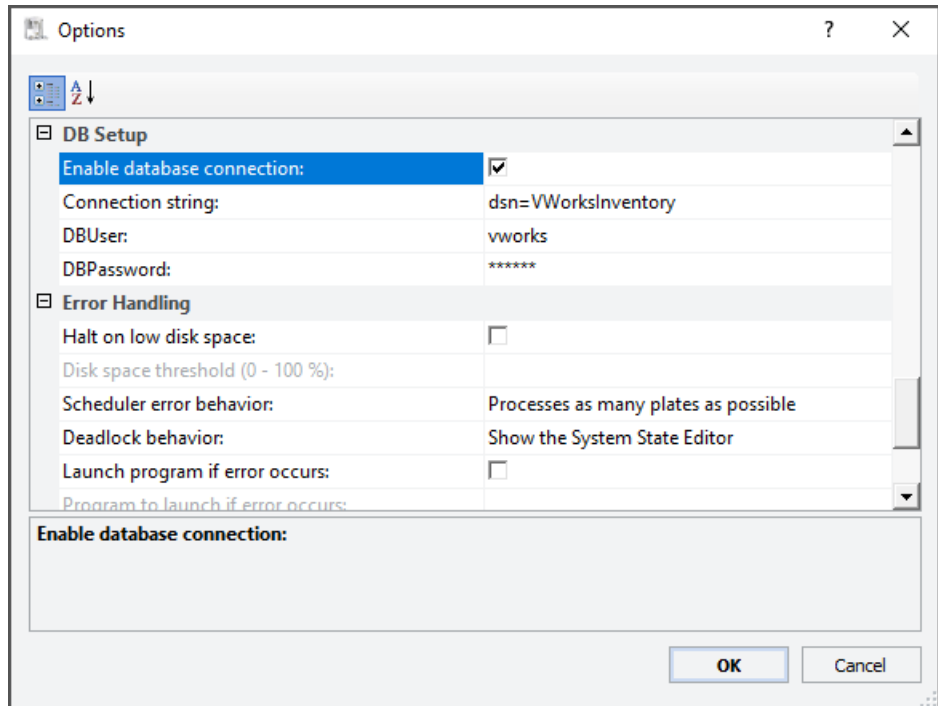
## Procedure

### To connect to the Inventory database:

- 1 In the VWorks window, click **Tools > Options**.
- 2 In the **Options** dialog box under **DB Setup**, verify the following, and then click **OK**.

Note: For MySQL 5.7, no entry is required for DBUser or DBPassword.

DB Setup parameters	MySQL 8.0.18	MySQL 5.7
Enable database connection	Selected	Selected
Connection string	dsn=VWorksInventory;	dsn=velocity11;
DBUser	vworks	
DBPassword	vworks	



## In This Book

This addendum to the VWorks Setup Guide describes how to install and configure MySQL 8.0.18 for VWorks 13.1.7.

[www.agilent.com](http://www.agilent.com)

© Agilent Technologies, Inc. 2021

February 2021



D0007558 Revision A

