

Berthold Detection Systems GmbH
Bleichstrasse 56–68
D-75173 Pforzheim/Germany



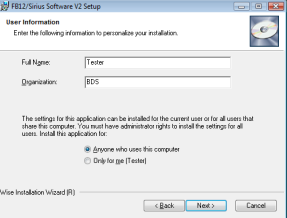

Phone: +49(0)7231/9206-0
Fax: +49(0)7231/9206-50
E-Mail: contact@berthold-ds.com
Internet: www.berthold-ds.com

Installing FB12/ Sirius Software V2.0 For USB Port of Sirius L

Product Information

Our tube luminometer Sirius L is equipped with a serial port as well as a USB port. For data connection via USB port, the installation of a USB driver is necessary. This driver software is part of the FB12/Sirius Software V2.0 package. It is therefore necessary to install the PC software prior to installation of the USB driver software.

Step 1: Installation of the FB12/ Sirius PC Software V2.0


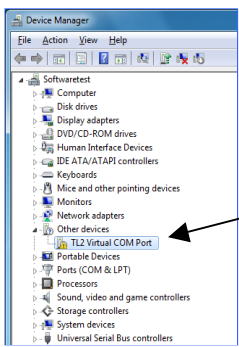
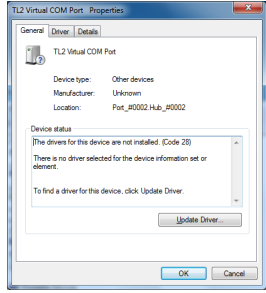
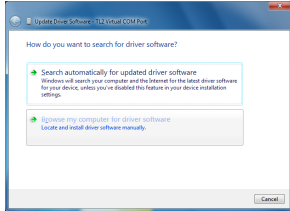
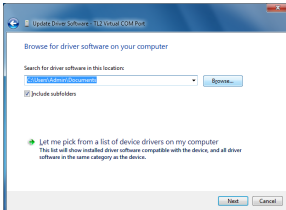
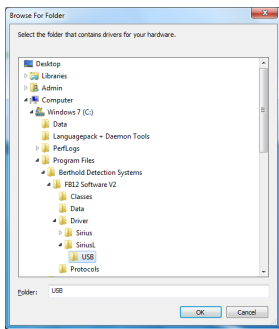
| | | | |
|--|---|--|---|
| <p>1. Insert your Product CD-ROM into your computer drive. The CD will automatically start.</p> <p>An installation wizard guides you through the software installation.</p> |  <p>Select Software. Click the FB12/ Sirius software icon. Double-click the Setup.exe file.</p> | <p>2. Follow the installation wizard instructions.</p> |  <p>Click <Next> to continue.</p> |
| <p>3. Enter user information and complete installation.</p> |  <p>Click <Next>, Follow remaining instructions</p> | <p>4. Finish Installation of Software.</p> |  <p>Click <Finish> The software is now ready installed. The FB12/Sirius Software icon is visible on your desktop.</p> |

Application Note 2010/02

Installing FB12/ Sirius Software V2.0 For USB Port of Sirius L

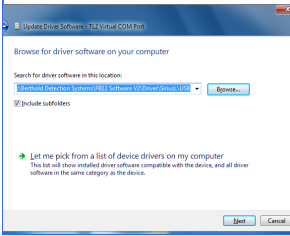
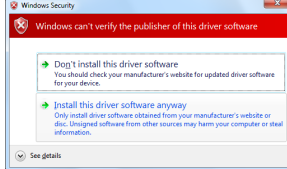
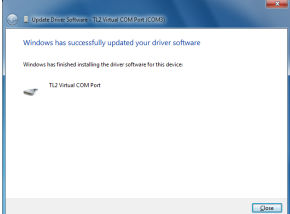
Step 2: Installation of USB driver

This description is valid for OS Win 7.

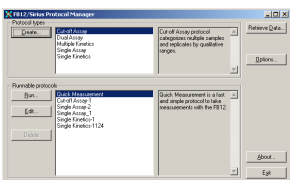
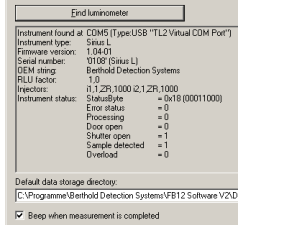

| | | | |
|--|--|---|---|
| <p>5. Switch on your luminometer and connect it to PC with the USB cable.</p> <p>The PC registers that there is a new device and starts to search for driver software, but does not find it automatically.</p> |  <p>Search for driver software in the following way:</p> | <p>6. Search for the driver software</p> <p>Click Start/Control Panel/System and Security/ System/Device Manager/Other Devices</p> <p>The TL2 Virtual COM Port Is present in the device manager.</p> |  <p>Click „TL Virtual COM Port“ with the right mouse-button</p> |
| <p>7. Update Driver installation</p> <p>A screen or a setting dialog appears.</p> |  <p>Click <Update Driver></p> | <p>8. Search driver software.</p> |  <p>Select „Browse my computer for driver software“</p> |
| <p>9. Browse for driver software on your computer</p> |  <p>Click <Browse></p> | <p>10. Browse for</p> <p>Computer /(C:) /Program Files /Berthold Detection Systems /FB12 Software V2 /Driver / Sirius L /USB</p> |  <p>Confirm your selection with <OK>.</p> |

Application Note 2010/02

Installing FB12/ Sirius Software V2.0 For USB Port of Sirius L

| | | | |
|--|---|---|---|
| <p>11. The selected location is assumed.</p> |  <p>Click <Next></p> | <p>12. A Windows Security Information appears</p> <p>The driver software is not licensed by Microsoft, but can be installed promptly.</p> |  <p>Select „Install this driver software anyway“</p> |
| <p>13. Complete the installation.</p> |  <p>Click <Close></p> | | |

Step 3: Verification

| | | | |
|---|--|--|---|
| <p>14. Double-click the FB12 Sirius Software V2.0 Icon on your desktop. The protocol manager will open.</p> |  <p>Click <Options></p> | <p>15. The options dialog apperas. Press the „Find luminometer“ button.</p> <p>Information about the instrument and the COM port is shown.</p> |  <p>Click <OK> to close the window. The protocol manager re-appears.</p> |
| <p>16. Start the Quick measurement protocol with <RUN>.</p> <p>The measurement window is opened.</p> |  <p>„RX“ appears on the right side of the status line. The system is now ready for use.</p> | | |

Best regards,
Berthold Detection Systems

March, 16th, 2010