

Sartorius **Wedge**

Getting Started

Demonstrates how to quickly capture measurements from a Sartorius device into Microsoft Excel.

Please visit taltech.com/sartoriuswedge for the latest information and product updates, including the most recent quick start guide, Excel files, WinWedge files, and troubleshooting tips.

Connect the Balance

First, let's connect your device to the PC. Most Sartorius balances use either a USB or serial cable, with USB being the preferred option. Some devices also have ethernet connectivity for accessing them over a network.

Additional steps may be required to properly connect your device to a PC. Please see your device's manual for complete steps.

USB & RS232 Serial

using WinWedge

USB: Plug the square end of the cable into the device and the rectangular end into a spare USB slot on your PC. Your PC should automatically install the necessary drivers.*

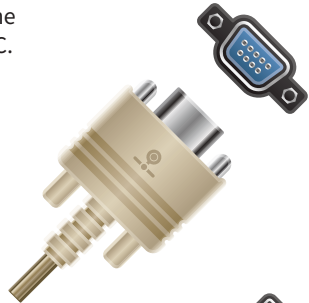
Serial: Plug the 25 pin female connector into the device and the 9 pin male connector into the PC.

For both Serial and USB devices, open the Device Manager on the PC and expand the "Ports" category. The available COM ports should be listed here. Keep note of the COM port number of your device.

TCP/IP

using TCPWedge

To set up the device for use over a network, connect an ethernet cable from the device to a switch or router on your network. The device may need to be configured to use the ethernet connection. See your device's manual or the complete quick start guide.



* If the PC does not find the drivers automatically, you may need to install them from your device's installation media. If you cannot locate the drivers, visit the Sartorius support web page for your device.

Open the Excel Workbook

There are several Excel workbooks included with the Sartorius Wedge package. Navigate to the WinWedge folder in the Start menu and select the Excel file for your device. The latest Excel files can be found online.

Initial Setup

The first time the Excel file opens, you'll need to configure a few basic settings. If the setup form did not automatically appear, press the "Setup button" on the first sheet. In this form, select the COM port number that the device is on. Also, type the timer interval (sec.) if you plan on using this feature. Click OK and WinWedge should automatically launch.

Capturing a Single Reading

To capture a single reading into Excel, press the "Print" button on your device or select the "Request Weight Reading" button on the Excel spreadsheet. This button sends a command to the device requesting the reading.

Capturing Multiple Readings on a Timer

If you'd like to collect multiple readings on a timed interval, click the Setup button to configure the interval and then click the "Enable Timer" button in Excel. This will enable the timer in WinWedge to send out a command to the device requesting data at the specified interval.

Troubleshooting Tips

1. Verify the Sartorius device is connected and powered on.
2. USB & Serial devices: verify the COM port is correct using Device Manager.
3. TCP/IP devices: disable the firewall or modify settings to allow local port.
4. Enable all macros in Excel macro security settings.

Read the PDF Quick Start Guide and the WinWedge Manual included on the CD and on taltech.com/sartoriuswedge for complete setup instructions.

Frequently Asked Questions

What is Sartorius Wedge?

Sartorius Wedge is a software package designed to capture data from Sartorius devices into Windows applications, specifically Microsoft Excel. Sartorius Wedge includes two of TALtech's award-winning data capture programs: WinWedge and TCPWedge. WinWedge connects to most serial or USB devices and TCPWedge connects to most networked devices.

Can WinWedge collect device data into any application?

This booklet shows how to capture data into Excel. However, you can also use WinWedge and TCPWedge to emulate the keyboard and send data to any Windows application. To do this, WinWedge must be configured in "Keystrokes Mode". More details can be found in the WinWedge manual PDF on the CD.

Why isn't any data coming into WinWedge?

Verify that all cables are securely plugged in and that the device is turned on. Then, open up WinWedge or TCPWedge, go to Port > Settings, and configure the settings EXACTLY as they appear on your device. To test the connection, go to Port > Analyze, press the "Print" button on your device. Data should populate in the Input Data field. If not, you either have the wrong settings for the device, the device isn't sending data, or there is a cable issue.

Why is data coming into WinWedge, but not Excel?

After WinWedge is activated, incoming data should appear in the WinWedge window and automatically transfer into Excel. If the data doesn't appear in Excel, verify that the Excel macro security setting is set to enable all macros. Also, make sure you're using the right spreadsheet for your device and that it is configured properly. If you get an error saying that the "COM port is not available", open up the Windows Device Manager and verify the COM port number. Also make sure there aren't any other instances of WinWedge running on that port.

Additional Troubleshooting

For more troubleshooting tips, visit taltech.com/sartoriuswedge