

Sartorius Wedge

Data Capture Software for Sartorius Balances

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.

Benefits of Sartorius Wedge:

- Eliminate manual data entry errors
- Log data from multiple devices simultaneously
- Capture real-time data to any Windows application
- Parse and filter the data to match your requirements
- 100% accurate, reliable, and compliant

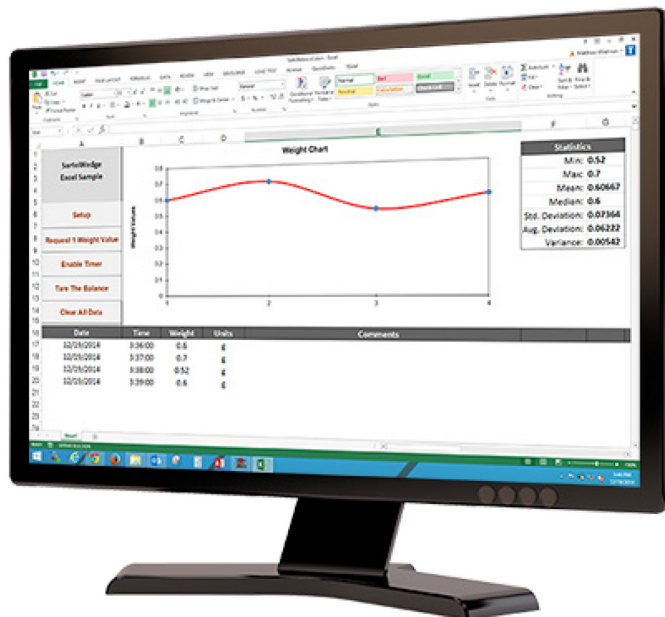


Excel Data Capture Spreadsheet Included

Everything you need to start capturing data into Microsoft Excel is included in the Sartorius Wedge package. The Excel files enable you to capture individual readings on demand or automatically capture readings on a timed interval. The readings are then charted in real-time.

System Requirements

Minimum: Windows 7, Excel 2016, Sartorius device and cables, USB port, serial port, or network card.



The Sartorius Wedge package includes software and files developed by TAL Technologies, Inc., specifically the WinWedge and TCPWedge software applications. © 2015 TAL Technologies, Inc. Sartorius is the distributor of Sartorius Wedge.

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

For further contacts, visit
www.sartorius.com