



SartoConnect.
The weigher-to-PC
data bridge.

sartorius

Why not transfer your data directly into your application program?

It's such a simple, professional way to bridge the data gap between your Sartorius weigher and your PC. You decide "what, when and where" – and let SartoConnect do the rest.

The SartoConnect Program

SartoConnect is a Windows 95/98/NT program for transmission of measured data. Data coming in through the serial interface can be inserted – together with additional information – directly into the desired PC application (e.g.: MS Word, MS Excel, Lotus 1-2-3, ..).

Alternatively, you can save the data in an ASCII file. Data can be transmitted either by hot-key functions or at time-based intervals. The data being transmitted can be freely formatted, filtered, and supplemented with additional information.

Above and beyond the weigher's own control system, SartoConnect lets you define your own data requests, and the weigher's replies can be sent to your receiving application. Once you have set your own specific parameters, SartoConnect saves them all in a configuration file.

Prerequisites

SartoConnect is suitable for all Sartorius laboratory balances and industrial scales, and IBM-compatible computers equipped with a serial interface.

Equipment Supplied

- 3.5" program diskettes in English/German
- RS-232 cable, optionally with D-sub 25 connector (YSC01L) or 12-pin round connector (YSC01I)



Example: MS Excel 5.0 diagram



Highlights

- Ease of operation
- User-definable formatting of measured data
- Remote control of the weigher
- Plausibility-limits check
- Time-based data acquisition
- Transmission of freely selectable text
- Supplementary identification, e.g., product ID
- 8 adjustable control commands
- Data acceptance via hot key
- ASCII-format storage

sartorius