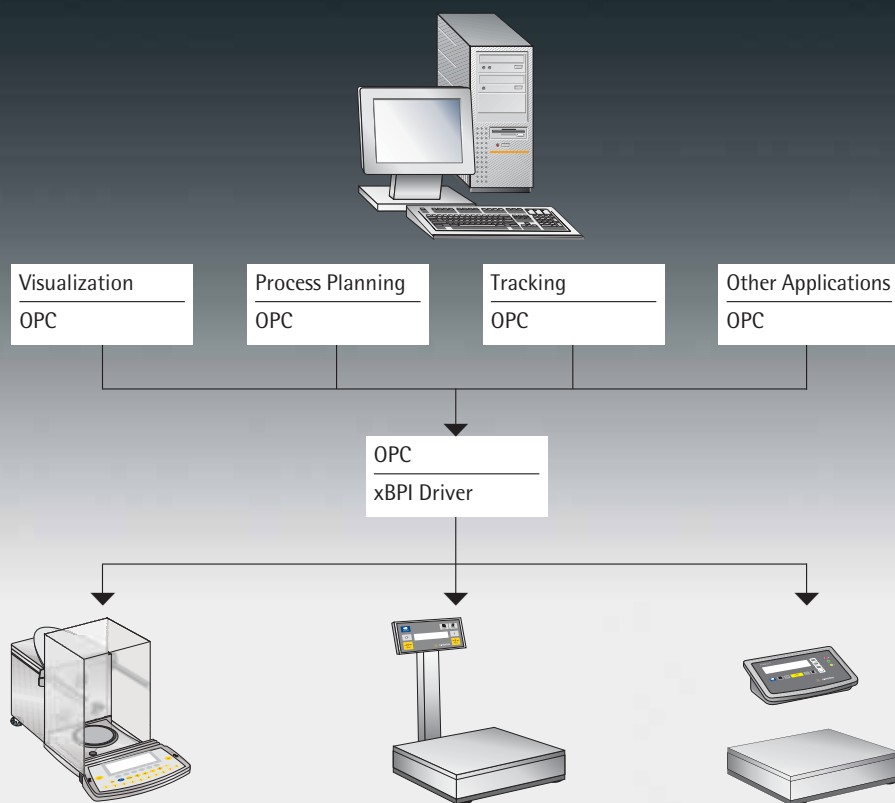




**sartorius**  
mechatronics



## Sartorius OPC Server

**A generous invitation\* to use  
Sartorius weighing equipment  
in higher-level electronic data  
processing systems.**

\* Visit our website at [www.sartorius.com/opc-server](http://www.sartorius.com/opc-server).  
There you will find your personal download version  
for free testing at no obligation.

Globalization, legal regulations (prepackage regulations and those on the net content of packages, FDA...), requirements (food safety laws, GMP, GLP) and standards (IFS, BRC, HACCP) entail comprehensive quality control and regular, 100% traceable data recording and logging. Without extensive networking of production machinery and software, it is hardly possible to integrate all information into a company's own data processing environment and to process it so that it is available to the particular programs used by this company. Sartorius now offers an OPC server for its xBPI weighing equipment – a standardized interface for simple connection of Sartorius balances and scales. The advantage is that it enables seamless communication of Sartorius equipment, which includes interface modules, components and transmission protocols.

### **Rely on standards**

The OPC server (OLE for Process Control) constitutes the bridge between hardware (Sartorius weighing equipment) and software. This server enables process and laboratory equipment to be easily integrated into users' particular applications. This means that users do not have to deal with a specific transmission protocol when they need to implement a Sartorius balance, scale, load cell or other equipment into their software. The OPC interface is subordinate to the application program and thus completely implemented by software.

Users of Sartorius process and laboratory instruments can now quickly and easily integrate them into and operate them in a network, without having to know much about programming. A further advantage: several client devices can simultaneously access process data.

# Sartorius OPC Server

62890PC | 62890PC-L –

**Our weighing equipment integrated into your app?  
No problem!**

- OPC server for all Sartorius laboratory balances and industrial scales that support the xBPI set of commands
- Easy use in nearly all of today's commonly used application development packages; OPC (OLE for Process Control) is the standard laboratory and process equipment interface
- Connection through RS-232 C | RS-485 P2P or through a serial | USB or serial | Ethernet adapter
- Suitable for Windows® 2000 | XP
- Is supplied with sample applications for simple conformance testing prior to use in your own application

## Equipment supplied

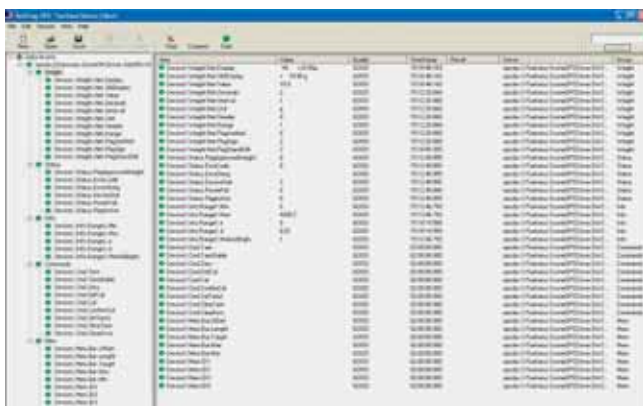
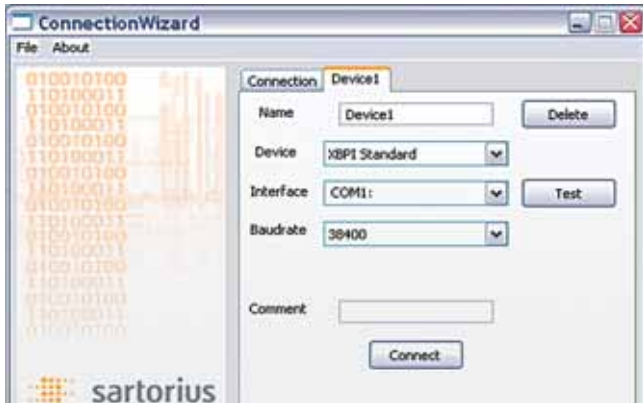
- Supplied by software download (here you will receive your personal 30-day free test version)...
- Sartorius OPC server for xBPI
- Installation | operating manual
- Website with access to weighing data as a sample application written in JavaScript
- Option: OPCScale sample application written in LabView 7.1 as a runtime version – requires LabView 7.1 runtime or a comprehensive full version with runtime and sources for use in your application (We will provide you with this module free of charge, without technical support.)

## Requirements

- IBM-compatible PC
- CPU at least 400 MHz
- RAM at least 256 MB
- Screen resolution at least 800 x 600 pixels
- Microsoft® Windows 2000 or XP operating system
- Sartorius balance or scale of the ME, MasterPro (LA), Combiics or Factory (FB, FC, FD) series, isi terminals, FCT indicators, IS digital platforms

## Advantages of the Sartorius OPC server

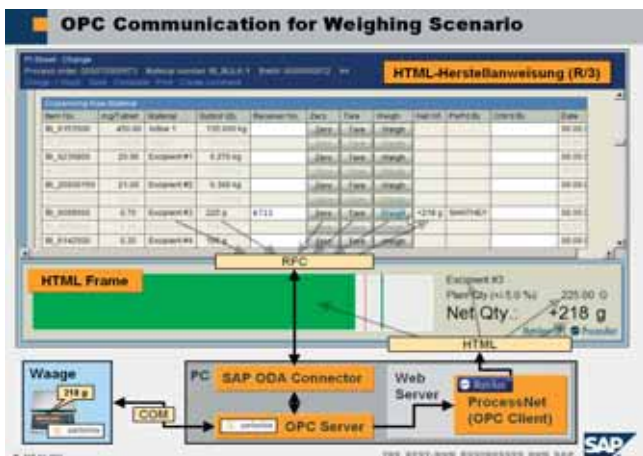
- Fast availability by download
- Fast installation as an application or Windows® service
- Easy use in application development software
- Supports OPC-compliant interface according to OPCDA and OPCXMLDA standards
- Up to ten balances or scales can be configured and up to three balances, scales or related weighing instruments operated simultaneously



The comprehensive set of commands for the OPC server is visible in the "softing" toolbox



Sartorius OPCScale sample application



Sartorius OPC server in the SAP R/3 weighing module

Sartorius AG  
Weender Landstrasse 94–108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

www.sartorius.com