



*Bluetooth*<sup>®</sup>  
Wireless Technology  
Putting the Bite on  
"Cable Spaghetti"

## Bluetooth® Wireless Technology\*. Wireless Weighing and Communication.

Both computer and measuring-instrument technologies are growing faster and more reliable all the time. The tools available for collection and evaluation of weighing data are becoming more powerful and, at the same time, easier to use. Only the data transmission path – the cable between weighing instrument and computer – has not kept pace with these rapid developments. This seems strange at a time when wireless communication has become so commonplace.

Bluetooth® wireless technology, widely used for laptops and mobile telephones, offers real advantages for both measurement and data storage processes. With a range of up to 100 meters, wireless connection of measuring stations, PCs and peripheral devices is now completely feasible for laboratory use.

No more cables to trip over, no more cable ducts collecting dust, no more inconvenient restrictions when positioning devices that have to be connected to one another.

Even the "line-of-sight" requirement imposed by infrared data interfaces is a thing of the past. Not only for mobile weighing, but also for clean room conditions or contaminated environments, Bluetooth® wireless technology presents a real alternative that eliminates connection problems before they occur.

Another major advantage of Bluetooth® wireless technology is the ability to connect multiple weighing stations in individual networks.

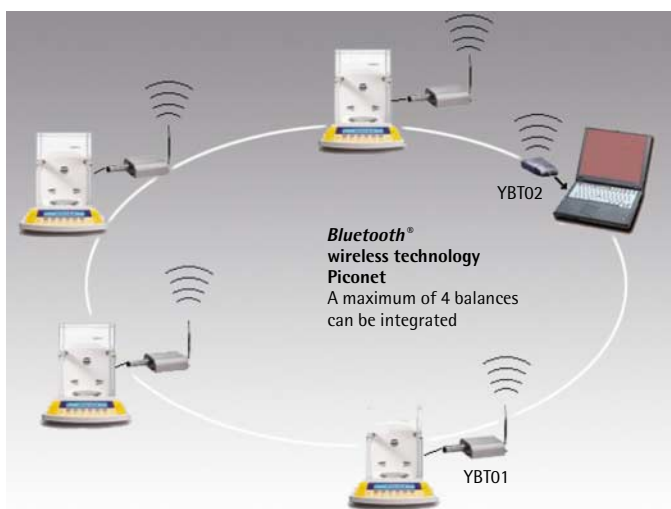
Installation of Sartorius communication modules featuring Bluetooth® wireless technology is as easy as can be. This technology uses the 2.45 GHz ISM band (for industrial, scientific, and medical usage). No fees are charged for this frequency, which means no added recurring costs for the user.

Data security is a high priority in Bluetooth® wireless technology. Data communication in both directions is protected by the use of frequency hopping, and other encryption techniques are also available. Thus, even sensitive areas are reliably secured.

With the YBT01 module for connection to the RS-232C port on the weighing instrument, and the YBT02 module for connection to the computer's USB port, Sartorius presents a solution that meets the most sophisticated requirements, with the same high quality as our Premium balances designed for use in the chemical and pharmaceutical industries.

The communication module has a stainless steel housing for optimal observance of the strictest standards of cleanliness. The data and data record transfer procedures will be familiar to anyone who has used RS-232C data interfaces. The new Sartorius network-capable SPEXXIS data acquisition program, is completely compatible with Bluetooth® wireless technology.

So put the bite on cable spaghetti with Bluetooth® wireless technology. The YBT01 and YBT02 modules are perfect for use with any of our Premium balances series, GENIUS ME, Master<sup>PRO</sup> LA or Expert LE.



### YBT01: RS-232C adapter with external antenna; for point-to-point connection only

Transmission power	In accordance with Class 1
Profiles supported	Serial port
Transmission speeds	1200 to 115,000 bit/s (configured by Sartorius Service)
Temperature range	0 to +40°C (+32 to +104°F)
IP rating	IP65
Dimensions (L × W × H)	121 mm × 84 mm × 32 mm (without antenna, cable, wall bracket)

### YBT02: USB adapter; for point-to-multipoint connections

Transmission power	In accordance with Class 1
Specification	Bluetooth® wireless technology V.1.1
Software	Bluetooth® device driver
Operating system	Windows® 98/2000, XP

Sartorius AG  
Weender Landstrasse 94-108  
37075 Goettingen, Germany

Phone +49.551.308.0  
Fax +49.551.308.3289

[www.sartorius.com](http://www.sartorius.com)

\* The brand name and logo for Bluetooth® wireless technology are owned by Bluetooth SIG Inc., USA. The use of this trademark by Sartorius is under license.

Specifications subject to change without notice.  
Printed in Germany on paper that has been bleached without any use of chlorine.  
W/sart-168 – W  
Publication No.: W-1096-e04032  
Order No.: 98649-002-21