



Sartorius YC013



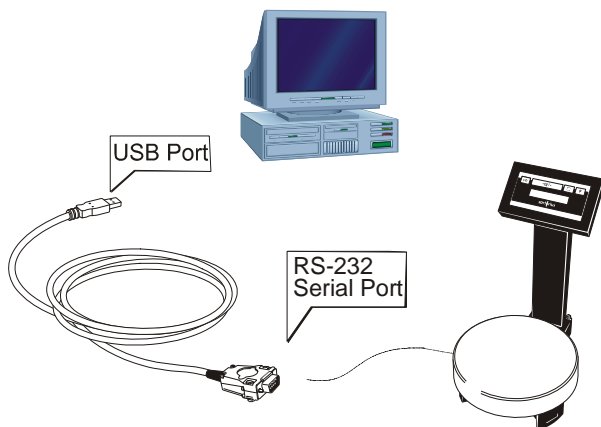
Product features:

- USB-connection cable to connect a PMA scale to a PC with USB-port
- For data communication using extended mode (BPI)
- USB 2.0 full speed device
- Can be used under Windows 98SE® / Windows ME® / Windows 2000® / Windows XP®

Standard accessories:

- RS232/USB connection cable
- Installation-CD with software drivers

Connection of a PMA to a PC



USB connection:	USB A-plug
RS232 connection:	9-pin D-SUB plug
Baud rate	300 Baud – 923 KBAud
Signal lines	RxD, TxD, CTS, DTR
Operating temperature:	0 – 40 °C
Cable length:	approx. 1,8 m
Net weight:	approx. 130 g
