



Signum® Complete Scale

Option A1 - UniCOM: RS-232 / YD001SW-232

Option A2 - UniCOM: RS-422 / YD001SW-485422

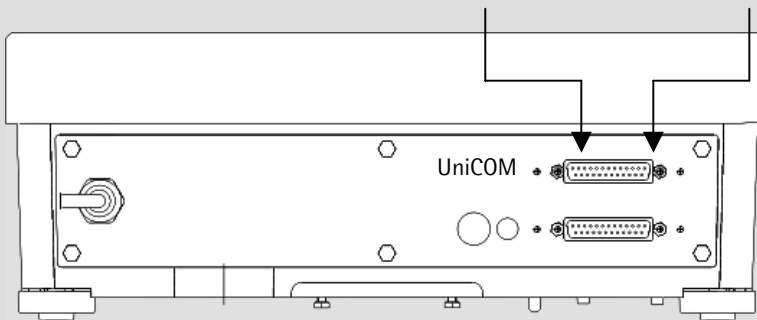
Option A3 - UniCOM: RS-485 / YD001SW-485422



RS-485/422



RS-232



- Options for a second interface on the scale (optional for all application levels) for data transfer to printers or external displays, or for data exchange with computers or controllers:
- The bidirectional RS-485 interface can be used to network up to 32 scales over an XBPI bus, or data exchange with a computer/controller or additional weighing platform with an RS-485 data output.
- The RS-422 and RS-485 interfaces are electrically isolated.
- The bidirectional RS-232 interface enables data transfer with different printers, data exchange with a computer/controller (SBI, XBPI or SMA operation) and connection of other accessories (see overview).
- The bidirectional RS-422 interface enables data exchange over distances of up to 1000 m.

Distances and protocols for the interfaces

Interface	Protocol	Maximum distance
RS-232	SBI/XBPI	20 m at 57 kbit/s (shorter distances required for faster speeds)
RS-422	SBI/XBPI	1000 m in accordance with the RS-422/485 specification
RS-485	XBPI	1000 m in accordance with the RS-422/485 specification

Connectable devices

RS-232 (Option A1)	Computer with serial RS-232 input port SBI/XBPI protocol YAM01IS external Alibi memory YBT01 external Bluetooth adapter YRD02Z second display USB adapter cable for connecting a computer over USB YCC01-USBM2 Second weighing point: scale with RS-232 data output Second weighing point: IS platform with optional RS-232 data output
RS-422 (Option A2)	Point-to-point connection over SBI/XBPI protocol. Important: interface converter for RS-422 to RS-232 required
RS-485 (Option A3)	Network up to 32 scales over XBPI bus Additional IS platform with standard RS-485 data output