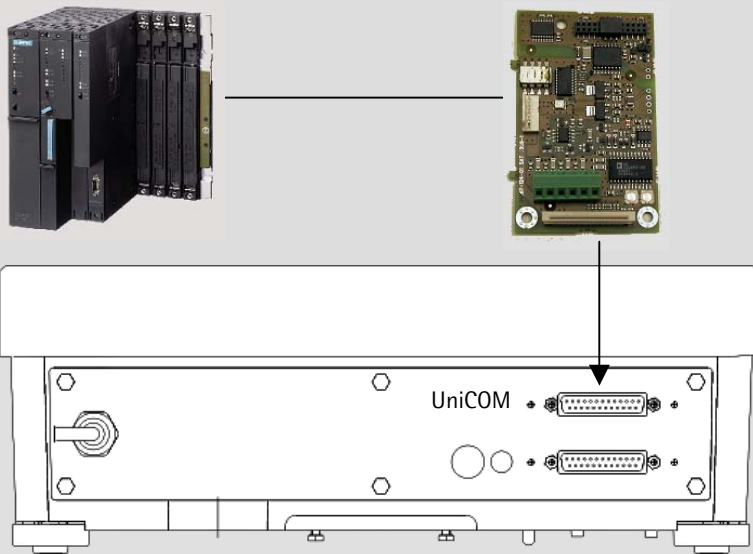




Signum® Complete Scale

Option A9 - UniCOM: Analog output port / YD001SW-A0



- The analog output port enables connection of a programmable logic controller or an external analog display unit

Data on Analog Interface

Electrically isolated data output

Can be used as voltage interface: standard: 0 to 10 V, or as current interface: 0 to 20 mA / 0 to 24 mA / 4 to 20mA (set by a switch)

- : 16 bit resolution
- : Current output:
 - Output error (at 25°C): 0.15% (max)
 - Drift of the output error: 25 ppm/°C (typical)
- : Voltage output:
 - Output error (at 25°C): 1.00% (max)
 - Drift of the output error: 100 ppm/°C (typical)

Connectable devices

Controllers with analog input