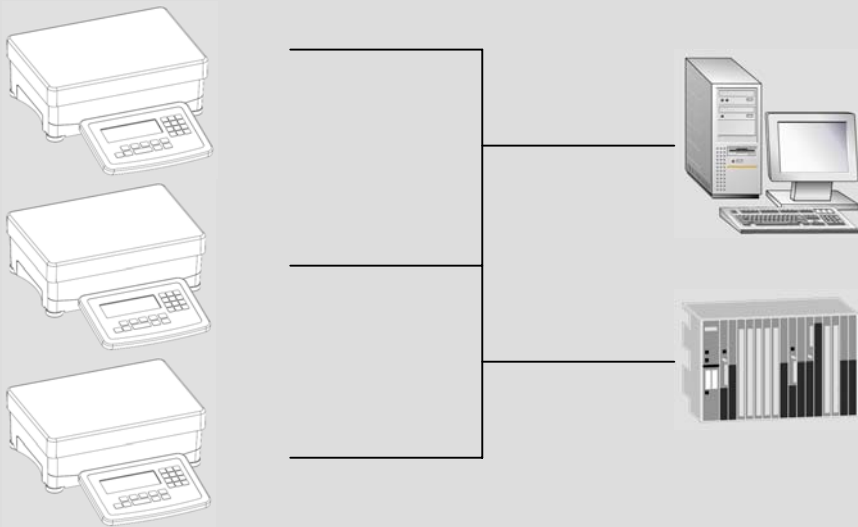




Signum® Complete Scale

Option B9: Ethernet Interface Port



- The interface supports Internet protocols:
TCP (data exchange with acknowledgement)
UDP (data exchange without acknowledgement of receipt)
DHCP (to use host names rather than IP addresses)
- Data speeds are automatically detected and selected:
10 Base T (10 Mbit/s) standard for embedded devices, such as scales
100 Base T (100 Mbit/s)

Connection

- : RJ45
- : CAT5 twisted pair cable

User data

The following protocols can be tunneled:

- : SBI (as client (master) or server (slave))
- : XBPI (as server (slave))
- : SMA (as server (slave): US standard scale protocol)

The Modbus/TCP protocol over Ethernet (as slave (server)) can also be used

The Unicon software tool can be used for data acquisition with a computer (order number: 6289-UniCon).

Furthermore, a "virtual COM port" (Com Port Redirector) can be set up to make the customer's existing serial controller software Ethernet-compatible for use with the SBI and SMA protocols.

Download:

<http://www.lantronix.com/device-networking/utilities-tools>