



## Signum® Complete Scale

### Option A5 - UniCOM: Digital 5 x IN/5 x OUT / YD001SW-DIO



- This optional data output enables direct control of signal transmitters and controllers.
- Electrically decoupled switch outputs can activate signals according to the specifications listed below (5 inputs and 5 outputs)

Digital Input Function	Level 1	Level 2	Level 3	Digital Output Function	Level 1	Level 2	Level 3
Zero/Stop calibration/Calibrate	•	•	•	Stable weight value (scale stability)	•	•	•
Tare	•	•	•	Large minimum sample quantity for application		•	•
Print	•	•	•	Small minimum sample quantity for application		•	•
Quit application/Delete one character (CF key)		•	•	Underflow L (underload)	•	•	•
"FN" key function (configurable in the the operating menu: toggle between weight units, gross and net values, or normal and 10-fold higher resolution)	•	•	•	Overflow H (overload)	•	•	•
Start function/application ("OK" key)		•	•	Control output "<"		•	•
"Info" key function (view application info or manual tare values)		•	•	Control output "="		•	•
Scale toggling		•	•	Control output ">"		•	•
Combined zero/tare function	•	•	•	Control output "Set"		•	•
Toggle between applications	•	•	•	Scale ready to operate	•	•	•
"isoTEST" key function	•	•	•	Data in tare memory			
Print GMP footer ("Print" key held > 2 sec)	•	•	•				
Switch scale on/off	•	•	•				

#### Digital inputs (electrically isolated by opto-electronic coupler)

Input voltage range -3 to -30VDC

Input voltage: -3V to 5V (logical 0)  $I_{(5V)}=100\mu A$

Input voltage: 11-30V (logical 1)  $I_{(15V)}=2.6mA$

#### Digital outputs (electrically isolated by opto-electronic coupler)

Output voltage range 0 to 30 VCD  $I_{max}=100mA$