









Quick Start - Used for Initial Start Up Only









Perform steps 9 to 20 for all units.

<p>NOTE: Once any Setup step is completed, meaning you pressed the  key a second time, you cannot repeat the step to make corrections. You can make corrections after the unit is started, refer to the manual Section IV.</p> <p> Press  to continue the setup procedure.</p>	<p>9  UnitS UnitS is the temperature display. Range: °C or °F Factory Default: °C</p> <ul style="list-style-type: none"> Press  The display will flash between UnitS and °C If desired, use  to change the scale to °F Press  to sequence to the next display 	<p>10  Hi t Hi t sets the fluid's High Temperature Alarm Limit. Range: +4°C to +42°C Factory Default: +42°C</p> <ul style="list-style-type: none"> Press  The display will flash between Hi t and 42 If desired, use  to adjust the value Press  to sequence to the next display
<p>11  Lo t Lo t sets the fluid's Low Temperature Alarm Limit. Range: +4°C to +42°C Factory Default: +4°C</p> <ul style="list-style-type: none"> Press  The display will flash between Lo t and 4 If desired, use  to adjust the value Press  	<p>12  Hi P1 Hi P1 sets the Pump's High Pressure Discharge Alarm Limit. Range: 4 to 100 psi Factory Default: 100 psi</p> <ul style="list-style-type: none"> Press  The display will flash between Hi P1 and 100 If desired, use  to adjust the value Press  	<p>13  dELAY dELAY is the length of time the pump can exceed the Hi P1 Alarm Limit before shutting down. Range: 0 to 30 seconds Factory Default: 0 seconds</p> <ul style="list-style-type: none"> The display will flash between dELAY and 0 If desired, use  to adjust the value Press 
<p>14  Lo P1 Lo P1 sets the Pump's Low Pressure Discharge Alarm Limit. Range: 4 to 100 psi Factory Default: 4 psi</p> <ul style="list-style-type: none"> Press  The display will flash between Lo P1 and 4 If desired, use  to adjust the value Press  	<p>15  dELAY dELAY is the length of time the pump can exceed the Lo P1 Alarm Limit before shutting down. Range: 0 to 30 seconds Factory Default: 10 seconds</p> <ul style="list-style-type: none"> The display will flash between dELAY and 10 If desired, use  to adjust the value Press  	<p>16  ALr ALr configures the unit's reaction to temperature, pressure, and flow (optional) alarm limits - either shut down (fLt) or continue to run (indC). See manual for more information. Range: fLt* or indC** Factory Default: fLt</p> <ul style="list-style-type: none"> The display will flash between ALr and fLt If desired, press  to display indC Press  *fLt = fault (shut down) **indC = indicate (continue to run)
<p>17  Sound Turns the unit's audible alarm on or off. Range: on or OFF Factory Default: on</p> <ul style="list-style-type: none"> Press  The display will flash between Sound and on If desired, press  to display OFF Press  	<p>18  StArT StArT enables/disables auto restart. Range: on or OFF Factory Default: on</p> <ul style="list-style-type: none"> Press  The display will flash between StArT and on If desired, press  to display OFF Press  	<p>19  CArE CArE is used to set the preventative care cleaning frequency reminder for the unit's air and fluid filters. Range: L1 - 1000 hours, L2 - 2000 hours, L3 -3000 hours Factory Default: L1</p> <ul style="list-style-type: none"> Press  The display will flash between CArE and L1 If desired, use  to change display to L2 or L3 Press 




















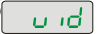



(If applicable, see boxes on right to set up options)

<p>20  StorE</p> <ul style="list-style-type: none"> Press  to save all settings The unit will automatically start. Press  to abort all settings The display will go blank Press  to restart the procedure 	<p>The Setup procedure is now complete.</p> <p>When the unit starts the controller will display the process fluid temperature.</p> <p>If desired, you can change/verify the unit's setpoint by pressing </p>	<p> SP SP is used to adjust the setpoint. Range: +5°C to +40°C Factory Default: +20°C</p> <ul style="list-style-type: none"> The display will flash between SP and 20 If desired, use  to change the setting Press  to return to the temperature display
--	---	--

Option - Flow Transducer (Steps A and B)

<p>A  Hi FLo HiFLO sets the high flow alarm limit. Range: 0.5 to 5.0 GPM Factory Default: 5.0 GPM</p> <ul style="list-style-type: none"> Press  The display will flash between HiFLO and 5.0 If desired, use  to adjust the value Press  	<p>B  LoFlo LoFLO sets the low flow alarm limit. Range: 0.5 to 5.0 GPM Factory Default: 0.5 GPM</p> <ul style="list-style-type: none"> Press  The display will flash between LoFLO and 0.5 If desired, use  to adjust the value Press  <p>If your unit does not have serial communications see Step 20.</p>
---	--

Option - Serial Communications (Steps C to H)

<p>C  SEr SEr is used to enable/disable and to configure serial communications mode. Range: off, rS232, rS485 Factory Default: off</p> <ul style="list-style-type: none"> Press  The display will flash between SEr and OFF If desired, use  to change the mode Press  	<p>D  BAud BAud is used to select the speed rate for serial communication. Range: 9600, 4800, 2400, 1200, 600, or 300 bits per second. Factory Default: 9600</p> <ul style="list-style-type: none"> Press  The display will flash between BAud and 9600 If desired, use  to change the rate Press 
<p>E  dAtA dAtA is used to display the number of bits per communication. Display: 8</p> <ul style="list-style-type: none"> Press  The display will flash between dAtA and 8 Press  	<p>F  StoP StoP is used to indicate the end of a communication's packet. Range: 2 or 1 Factory Default: 2</p> <ul style="list-style-type: none"> Press  The display will flash between StoP and 2 If desired, use  to change the setting Press 
<p>G  PAr PAr is used as a means to check for communication errors. Range: even, odd, or none Factory Default: none</p> <ul style="list-style-type: none"> Press  The display will flash between PAr and none If desired, use  to change the setting Press  	<p>H  u id u id (user id) is used in RS485 only. It sets the number of devices connected to the RS485 port. Range: 1 to 99 Factory Default: 1</p> <ul style="list-style-type: none"> Press  The display will flash between u id and 1 If desired, use  to change the setting Press  <p>See Step 20.</p>