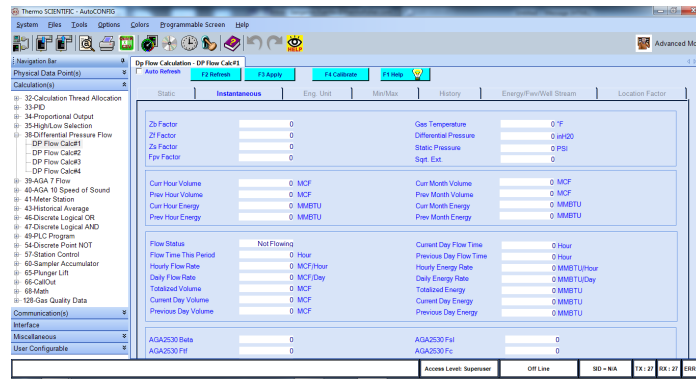


Frequently Asked Questions



Common questions for AutoCONFIG software for use with our Flow Computers

Which Thermo Scientific instruments have AutoCONFIG software integrated?

- AutoXP Gas and Liquid Flow Computer
- AutoEXEC 32-Run Expandable Gas and Liquid Flow Computer
- AutoPILOT PRO Multi-Run Hydrocarbon Flow Computer
- AutoXP-SMV Smart Multivariable Transmitter
- SOLA iQ On-line Sulfur Analyzer

Are you able to monitor multiple wells simultaneously?

Thermo Scientific AutoCONFIG™ software is capable of recording multiple wells including tubing and casing from each well, along with other critical process variables to create trend reports allowing the user to take corrective action where needed throughout the process.

Can you consolidate measurements?

Configured for multiple primary devices, the software can consolidate oil, gas and water measurements for maximum control and efficiency.

Can you calculate net tank production?

By integrating with a variety of tank gauges AutoCONFIG will record and archive net tank production and withdrawals.

Are you able create user-configurable screens?

The AutoCONFIG software is designed to empower the user through the capability to create custom user-configurable screens.

What type of communication ports connect to the AutoCONFIG software?

Connect to AutoCONFIG software via RS232, RS485 or Ethernet ports.

Is it possible to save a connection list?

For simplicity, a connection list can be created, saved, and shared amongst technicians.

Can you open multiple sessions?

Yes, you can open multiple sessions of the software at the same time.

To simplify the view, can features/functions not active be hidden?

There are two modes, advance and standard. Standard mode will hide all non active/disabled functions. A user can switch between these modes, but the standard mode hides all functions that are not enabled.

Are separate tools or software required for custom views?

No, the configurable screen development tool is accessible when you download the AutoCONFIG software.

Can custom displays and configurations be shared with other users?

Easily email or store files on a corporate sharepoint for access by other users.

Do PIDs require programming?

No, firmware based PIDs provide all control and switchover logic that is needed. In addition, AND/OR logic and math tables provide increased configurable control.

What resources are available for training?

Request a live demo or view the user guide at www.thermofisher.com/autoconfig

Find out more at thermofisher.com/autoconfig