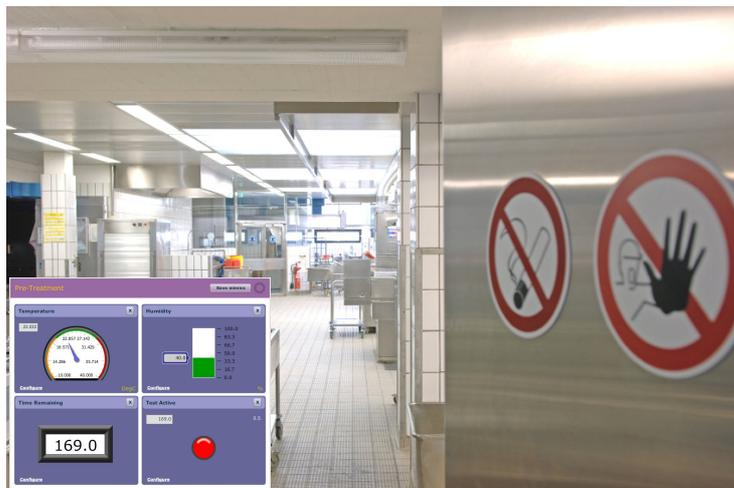




Customer Requirements

A decontamination and sterilisation facility uses Ethylene Oxide (EtO) preconditioning rooms as part of the sterilisation process. The environment within these rooms needs to be monitored for both safety and quality assurance purposes. The huge range of products sterilised require a minimum temperature of 22 degrees Celsius and 40% humidity level. The two stage preconditioning process takes approximately 12 hours.

The logging solution used must be easy to use, provide alarms for unsuitable conditions and simple to retrieve data from.



Decontaminate it!: A food sterilisation facility uses *dataTaker* and *dEX* to monitor the EtO preconditioning of food products.

dataTaker DT80

- 1 A cost effective data logger expandable to 100 channels, 200 isolated or 300 single-ended analog inputs
- 2 Built-in web and FTP server allows for remote access to logged data, configuration and diagnostics
- 3 Modbus slave and master functionality allows connection to Modbus sensors and devices and to SCADA systems
- 4 Smart serial sensor channels capable of interfacing to RS232, RS485, RS422 and SDI-12 sensors
- 5 Rugged design and construction provides reliable operation under extreme conditions
- 6 Includes USB memory stick support for easy data and program transfer



dataTaker Solution

Equipment

dataTaker DT80 data logger

Sensors

Humidity
Temperature

Implementation Notes

The data logging system comprises a *dataTaker* DT80 data logger and a workstation running a web browser. Through the web browser, the *dataTaker* web interface (*dEX*) can be accessed. *dEX* allows users to configure the data logger, monitor the connected sensors, collect data and check the logging and memory status, all from within a web browser. The logger can be scheduled to sample and record the temperature and relative humidity at predetermined intervals such as every second or minute from up to 15 single-ended sensors placed in the EtO preconditioning room(s).

Mimics can be used to visually monitor the progress of tests and to clearly identify to the facility staff when a test is in progress for a specific chamber.

Within the logger, alarms can be configured to automatically alert staff when the preconditioning environmental conditions are outside the acceptable levels.

The data can be later downloaded from the logger and saved for quality assurance purposes in either CSV or DBD formats either via *dEX* or using a USB stick.