Process NIR Instrument













Reliable, Accurate, Real-time





Modern on-line process sensor.

The DA 7440 measures moisture, fat, sugar, flavorings and more in real-time. It's widely used in various food and agricultural processing applications and is used to monitor incoming ingredients and raw materials, optimize production processes and verify finished product quality. It's typically installed over conveyor belts but its flexibility allows for other options as well.

Results are easily integrated into an existing automation system. Thanks to its use of modern diode array technology the DA 7440 gives accurate and stable readings even when your products change.



Optimize your production

The DA 7440 is successfully used to measure and control moisture and more in various processes. By accurate, real time measurements, the user can automatically or manually control drying, cooling, ingredient additions, blending or segregations, depending on current needs.

The modern DA technology automatically detects if there is product on the belt or not and handles process changes such as varying product height, colour and temperature without any undue effects on the measurement results.

Information integration

By integrating the instrument into the SCADA system, the operator can see the results in a familiar operating environment. Further, the results are available for automated controls and reporting through a central reports system. The DA 7440 offers a range of communication options including.

- OPC DA
- Fieldbus including Profibus DP, Modbus, etc.
- Ethernet TCP/IP based Modbus/TCP & PROFINET
- Analogue outputs

Further, data can easily be downloaded from the web-based user interface for evaluation.



Industrial Design

The DA 7440 is available in different models for different processes. The DA 7440 General Purpose is typically used in dry processes, whilst the DA 7440 Sanitary Design version is used in food production areas or other environments that require higher ingress protection. A cooled version of the instrument is also available for installation in areas with elevated temperatures. The Smart Sensor is equipped with an Air-purge window that ensures the sensor window is kept clean and free from contamination.

The sensor communicates over Ethernet and can be plugged straight into an existing plant network or connected to a wireless network.

This connectivity reduces the need for additional cabling to be run separately.

The light indicator on the sensor indicates operating status and alarm conditions with different colors, making it easy for operators to see that everything is working as expected.





Process Plus

Process Plus is the dedicated process instrumentation software from Perten Instruments. Process Plus is a modern and easy to use software that is built with the specific demands of the process industry in mind. The software includes a web-based user interface with both numerical and graphical displays as well as functions for setting up instrument and product specific features. Process Plus can be installed either on a dedicated PC in the junction box or on a server with automated back-up and redundancy. Process Plus is part of the Perten Plus software suite.

Operator At-line display

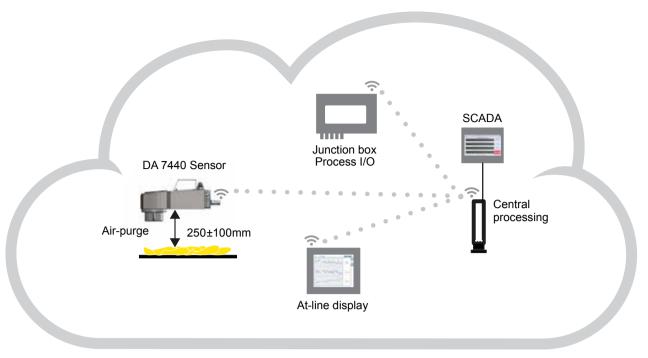
Further enhancing the functionality of the system, we offer an at-line display. The large and clear screen provides at-a-glance information of actual measurement results in graphical or numerical form. The instrument status is shown with signaling and status messages, clearly showing the current status. The touch-sensitive display is also used when taking grab samples. When the sample button is pushed, the instrument tags the corresponding measurements in the instrument database for use at later validation or calibration enhancements.

Junction box To ease the installation, we offer a junction box that includes:

- 24V DC power supply & UPS
- 4 port switch
- Wireless access point
- Industrial computer for Process Plus
- Gateway for Analogue output and Fieldbus interfaces
- External sample button input

The junction box is typically mounted in the vicinity of the sensor and provides a convenient way to quickly install the required supplies to the sensor.





Multiple sensors with one processing unit is possible.

Modern communication infrastructure

Communication between the various system components is over Ethernet. The system is available with wired or wireless interface between sensor, central processor and peripherals. The user interface is web-based, meaning that any user can access real-time measurement information from a web-browser anywhere on the network. For example, whilst the operator is using the At-line display, the production manager can view measurements in his/her office in real time. The laboratory can, simultaneously and independently, add reference analysis results to the manual sampling report, allowing a quick and continues verification of sensor performance. This multi-user capability requires NO additional software to be installed on any computer.

Advantages

- Accurate and precise results for all products
- Web-based user Interface
- Easily integrated into SCADA



Diode Array 7440
Process NIR Instrument

Manufactured by Pertentiation of the makes

Calibrations In difference to instruments that use only a few NIR wavelengths to make the measurement, the Diode Array technology captures a complete part of the NIR spectra. This facilitates the use of modern calibration techniques such as PLS, the proprietary HR or ANN. Perten offers ready-to-use calibrations for all important parameters and continued support for the calibrations. Alternatively, the customer can chose to develop and maintain PLS calibrations, using third party software, the Unscrambler® from Camo.

Support Managing ones assets in order to ensure optimum performance becomes more and more profitable. Perten offers a range of affordable support packages to ensure your instrument keeps performing as new, year after year. The support includes calibration maintenance and up-dates, replacement of certain consumables as well as software updates.

Specifications

Temperature range: 0°C - 40°C, 0-70° with optional cooler

Ingress protection: IP65 or IP69k

Humidity: 50% RH ±30% Mounting: 38 mm diameter

Power: 230 V AC 50-60 Hz to the junction box.

Network: Ethernet/TCP-IP 100/100 Mbps network to sensor, cabinet and at-line display

Compressed air for air purge: 3 bar, 6 mm diameter.

Flow rate: 30-60 l/min.

