

Diode Array 7250

At-line & Lab NIR Analysis System



Fast



Easy-to-Use



At-line Ruggedness



Web Interface Software

Accurate Analysis – Anything, Anywhere, Anytime, Anybody.



Industries: Grain and Oilseed Processing, Dairy, Meat Processing, Animal Feed, Flour Milling, Petfood, Snackfoods, Starch Production, Sugar Production, Bakeries, Ethanol Production, Wood-based Products.

Parameters: Moisture/Dry Matter, Solids, Protein, Fat/Oil, Ash, Fiber, Starch, Amino Acids, Fatty Acids, Sugars, Brix, Lactose, and more.

Fast, Accurate, Easy-to-use, Versatile, Robust.

Key features

- 6-second analysis of all types of samples
- Modern technology and a wide range of available calibrations
- Easy sample handling and touch-screen operation
- IP65 certified for dust- and water resistance
- Anytime access to results via web browser



Diode Array 7250

At-line & Lab NIR Analysis System

Accurate Analysis -

The DA 7250 simply makes sense. It's fast, accurate, easy-to-use, versatile and robust. Our years of experience and expertise have guided its design – optics, housing, sample presentation, workflow and software simplification. It comes in two versions – one for general use and a stainless steel, sanitary design version, for use in the food industry.



Robust and reproducible Diode array technology allowed us to create an instrument with no moving optical components. This improves accuracy and stability, and means that hardware can be standardized, minimizing instrument differences. The DA 7250 sets new standards in calibration transferability. It also removes sources of potential failure as there are no expensive monochromators to wear out and no wiggling mirrors that can be subject to vibration.



Fast and accurate spectra The DA7250 uses a sensor with excellent signal/noise ratio and sensitivity. In combination with the outstanding energy throughput, the DA7250 collects more than 20 spectra per second, as the sample moves. The instrument is equipped with advanced automatic standardizations of the wavelength and absorbance scales. This improves accuracy, corrects for changes in ambient conditions, and alerts users to potential problems.



Connectivity The DA 7250 instrument and software platform is designed for optimal connectivity and data is readily exchanged with other programs. Ethernet and USB ports are provided and Windows peripherals such as a Wi-Fi dongles, cellular/mobile modem, bar code readers and printers are supported. The innovative cloud-based NetPlus software provides measurement result viewing from anywhere using web browsers. The instrument can be remotely administered through NetPlus or Team Viewer.



Anything, Anywhere,

Grains, powders, pellets, pastes, slurries and liquids

Anything

Grains, powders, pellets, pastes, slurries and liquids – nearly any type of sample – can be analyzed on the DA 7250, with no requirement for expensive, cumbersome modules. A large, representative sample area is analyzed meaning little or no sample preparation is required. The use of open-faced dishes removes tedious and time consuming cleaning between analyses. Available factory calibrations cover a wide variety of products and parameters and are built from our global database which encompasses hundreds of thousands of samples. The DA 7250 is ISO 12099 compliant, ensuring analysis that follows international standards.

Anywhere

The DA 7250 is as at home in the lab as it is at-line in processing environments. It uses no moving optical components and employs a solid state drive (SSD) making it vibration resistant. It is IP65 certified (dust proof/water resistant) and temperature stabilized for consistent results. The DA 7250 excels in dusty, dirty and otherwise challenging environments. In food production areas the stainless steel Sanitary Design model meets hygienic requirements. It is designed with an open analysis area and without sharp corners, making it easily cleaned.

In the Lab & At-line



Anytime, Anybody.



Anybody

Plant personnel to experienced chemists find the DA 7250 easy-to-use yet powerful. Its large touch screen with intuitive user interface and straightforward sample handling ensures rapid operator confidence. Operating procedures are designed to eliminate sources of operator error producing accurate results regardless as to who performs the analyses. Configuring and validating instrument measurements is easily accomplished using Netplus Remote web based maintenance software. Accurate analysis is available at all times, enabling you to optimize and verify product quality 24/7, improving efficiency and reducing scrap.

Anytime

The DA 7250 is fast; test samples anytime, with optimized workflow, little or no sample preparation, in just 6 seconds. When analysis results are needed, the DA 7250 provides them immediately. It automatically performs background standardization tests ensuring accuracy. Not only can samples be analyzed in essentially real-time, but analysis results are available anytime from any location around the world using NetPlus Reports Web based software.





Lab, At-line and Process

The DA 7250 can be used in the lab and at-line. It is available in two models; General Purpose (GP) and Sanitary Design (SD). For continuous process analysis we offer two sensor models. The DA 7300 series is used for in-line analysis and the DA 7440 series for over-the-belt on-line analysis. As they are based on the same NIR technology, many calibrations can be transferred between the models.



Web interface reporting and instrument management

The DA 7250 is supported by the NetPlus web based cloud service. Instruments are connected through a local network or internet and are accessed from anywhere using web browsers.

NetPlus Reports provides access to analysis results. Monitor production, verify quality of ingredient shipments or get an update on latest analyses. See the results in tables and charts on your laptop or tablet.

NetPlus Remote lets you configure instruments, monitor performance and update calibrations. Whether you manage one instrument or an NIR network of one hundred instruments, NetPlus Remote streamlines your tasks.

Accessories

A wide range of sample presentation accessories are available for the DA7250. Nearly every type of sample, solids to liquids, can be analyzed with little to no sample preparation.



Specifications

Analysis time: 6 seconds

Sample types: Grains, Pellets, Powders, Pastes, Slurries, Liquids and more

Sample area analyzed: Up to 108 cm²

Sample volume: Flexible 1-400 ml, depending on sample type

Display: 12" color touch screen

Language: Multiple languages included. English, Chinese, Thai, Russian and many more

Interfaces: USB and Ethernet

Operating temperature range: 5 to 40°C

Ingress Protection per IEC60529: IP65 certified

Wavelength Range: 950-1650 nm

Detector: Indium Gallium Arsenide (InGaAs)

Wavelength Accuracy: < +/-0.05 nm

Power requirements: 115 or 230V, 50 or 60 Hz (specify by order)

Dimensions (HxWxD): 517 x 370 x 390

Net Weight: 13 kg