

# Inframatic 8800

NIR Grain Analyzer



On-farm



Take Anywhere



GPS



Protein Maps

Accurate Portable Grain Analyzer

**Perten**  
INSTRUMENTS  
a PerkinElmer company

# The take-anywhere answer to your grain quality questions.



*Near-infrared (NIR) grain testing has been a fixture at large elevators and grain processors for years. With the use of newly available technology, Perten Instruments is excited to introduce a tester designed specifically for on-farm use.*



**Designed** for farm conditions, this rugged meter is ideal for testing grain for moisture, protein and oil. An automated shutter protects the tester from light, insects and vermin to ensure reliable, low maintenance operation for years to come.

**Portable** – Compact, light enough to hand carry and powered by 12V, or battery (up to 2 hours), the Inframatic 8800 goes with you where you need it most: in the pickup truck, in the cab of the header and to the silo. The optional carrying bag protects the IM 8800 and is useful for transporting the instrument to inspection sites, fields etc. It's also useful for long-term, protective storage of the instrument after harvest.

**GPS** – The Inframatic 8800 is equipped with a GPS so you can create protein maps of your fields. Make fast harvesting and binning decisions to manage your grain quality.

**Protein maps** allow farmers to extract additional profit by identifying pockets of premium grain. Topography, fertilization, and run-off are known to cause variation in protein content within a field. The Inframatic 8800 helps you get the most value from the grain you are already producing.



Measure in the Field



## Why the Inframatic 8800 is different

An on farm grain analyzer has been desirable for some time. The demands on such an instrument are high, however, and must balance many requirements. It must be rugged and robust. It must be portable to tote around the field. It must be simple to operate, but sophisticated enough to provide accuracy similar to elevator and lab instruments. And it must do it all in a cost-effective manner.

Recent technological developments have allowed us to develop just such an instrument. The Inframatic 8800 uses solid state components and diode array technology meaning no moving optical components. The lack of moving parts allows us to align and match instruments at the factory – one instrument is exactly like the next. This means the instruments are accurate, repeatable, and reproducible.

## Measure in the field

It's all about portability and fit-for-purpose with the IM 8800. Place it in the cab of your pickup or harvester. It's small enough to carry around the field for spot testing allowing you to determine optimal harvest time. The 2 hour battery life and GPS allow you to create protein maps of fields.

Variability of protein in wheat within a single row has been shown to be up to 6%. The maps allow you to capture that value. Maximize profitability by better controlling drying, identifying your highest quality grain, and selling it for a premium.

## Quality Control before delivery

Deliver the right grain to the right buyer. Test your grain as it's going into storage and as it's going into the truck. Send the highest quality grain to those willing to pay premiums. Make sure you get the most for your grain.

*"Get the most from your grain with fast, accurate analysis."*

Before delivery





## Standardized and Matched

The Inframatic 8800 is standardized to our elevator/lab instrument – the Inframatic 9500. The Inframatic 9500 is well established in the market and is an approved for payment instrument in many countries. It has approvals in the USA, Germany, France and Australia to name a few.

## Perten Instruments

With 50+ years in agri-food business, we have the experience and expertise to make your use of the IM 8800 successful.

## Perten Support

Through our extensive group of trained distributors and direct offices, Perten provides unparalleled support of our instruments. It starts with instruments engineered to meet the needs of our customers and the demands of the environments in which they are used. It extends to the software provided for routine operation, networking and remote administration, and – ultimately – installation and training. Your purchase of an instrument is the beginning of our relationship – not the end.

## Specifications

**Products:** Grains and Oilseeds

**Parameters:** Moisture, Protein, Oil, Wet Gluten and more

**Available Calibrations:** Wheat, Durum, Barley, Canola/Rape seed, Sorghum, Corn and Soybean

**Analysis Time:** ~90 s

**Sample Size:** ~400 ml

**Subsamples:** Up to 10 per sample

**Analysis Principle:** Diode-array detector, Transmittance

**Wavelength Range:** 850-1050 nm

**Size (W x D x H):** 349 x 265 x 274 mm

**Weight:** 7 kg

**Interfaces:** 4 x USB-A ports, 1 x Ethernet port (RJ45)

**Display:** 5.7" color touch screen

**Protection:** Dust and humidity protected

**Battery Operation:** ~2 hrs.

**Positioning:** GPS module, connected through USB port

**Ambient Temperature:** 5-45 °C

