## Grain Moisture Meter









Oilseeds



Grain Intake

Revolutionary Grain Moisture Accuracy



Design a Quality of Perfect Particular of the AM experience in every a moisture in The higher Unif

Aquamatic 5200 Grain Moisture Meter

The AM 5200 is Perten's second-generation high frequency moisture meter. Drawing on experience from more than 5,000 installed units, we have designed a moisture meter improved in every aspect – accuracy, repeatability, reliability, and user interface. It can be a stand-alone moisture meter for use at terminals or can be integrated into automated testing systems.

The higher frequency used penetrates deep into samples for improved results. The updated Unified Grain Moisture Algorithm (UGMA) provides a more accurate analysis regardless of temperature or crop type. The improved mechanisms provide better repeatability. Together these make the AM 5200 the most accurate and repeatable moisture meter available. It analyzes grains, oilseeds, pulses, beans, lentils, seeds and more for moisture, Test Weight/Hectoliter Weight, and temperature.

## Features and Benefits

**Superior Accuracy:** Superior accuracy and repeatability are obtained using a 150 MHz measurement cell made of robust, high precision die-cast aluminum. The new cell is slightly larger providing a more representative sample. The large contact area of the temparature sensor with the grain provides a superior temperature accuracy from larger representation of grain in comparison to other moisture meters.

**Accurate Results on Fresh Field Samples:** Older moisture meters read up to 1% low in moisture when analyzing grain straight from the field (known as the rebound effect). The higher frequency measurement employed by the AM 5200 penetrates deeper into the grain providing more accurate results and shorter return on investment.

**Easy To Use:** The large color touch screen with intuitive user interface makes the AM 5200 simple to use. Input sample IDs, view results on a remote screen, and update through USB memory stick all add to its flexibility and ease of use. The touchscreen uses resistance technology and even responds to gloved hands.

**Rapid Analysis:** Just pour the sample and the AM 5200 begins analysis automatically. Results are displayed in less than 10 seconds. An optional flow-through provides for automated testing when the sample can be discarded.

**Modern Technology:** Using higher frequency, advanced computational power, USB data transfer, and the updated UGMA, the AM 5200 is designed for today's grain trade.

**USDA Certified:** The USDA certified the AM 5200 as UGMA compatible and for use for official moisture measurements.

**Officially Approved:** The AM 5200 is officially approved in a number of countries including the USA, Germany, France, Romania, Serbia, Croatia and Tunisia.

## Specifications

**Products:** Grains, oilseeds, pulses, lentils and more **Parameters:** Moisture, specific weight (hectoliter

weight/test weight) and temperature Analysis Time: ~ 10 seconds Display: 5.7" color touch screen Connectivity: Ethernet, USB

Measurement Technique: RF dielectric

constant at 150 Mhz

**Sample Temperature:** -20 to  $45^{\circ}$  C / -4 to 113 F (moisture <18%); 0 to  $45^{\circ}$  C / 32 to 113 F (moisture >18%)

Sample Size: 700 ml

Power Requirement: 115/230 V, 50/60 Hz

Dimensions HxWxD (mm/inches):

415/16.3 x 424/16.7 x 353/13.9 **Net Weight (kg/lb):** 18.3/40.3



