PERTEN FALLING NUMBER WORLD STANDARD METHOD



Perten Falling Number (FN 1000)

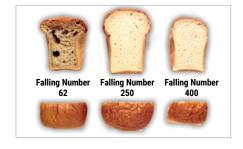
Approved Methods

- AACC No. 56-81.04
- ICC No. 107/1
- ISO No. 3093

The Only Validated Instruments for the Approved Methods

Falling Number Method

Alpha-amylase activity has great influence upon the quality of baked goods, pasta and noodles. Sprout damage is caused by alpha-amylase – a naturally occurring enzyme in grain that increases in concentration during wet harvests. Perten Falling Number instruments by PerkinElmer are the only validated instruments capable of running the World Wide Standard method for measuring alpha-amylase activity in flour and meals of wheat, durum, rye, barley, other grains and malted cereals.



The Falling Number method is a fast and easy test to help protect your organization from the effects of sprout damage. The FN method provides a common international quality-language along the whole production and process chain - from wheat growers to bakers - to ensure quality, efficiency and fair trade.

Perten Falling Number® 1000 instrument

The Perten Falling Number 1000 is an automatic dual sample analysis system designed for simple operation. The system includes functions for automatic water level control, automatic start and automatic stop at operator set FN target. It also includes functions for registration of sample ID, calculation of moisture corrected sample weight, mean value calculation, moisture corrected results, altitude correction, and calculation of blends and malt addition. The user can also select the optional Fungal Falling Number method. With its 5.7" touch screen, operation is simple and intuitive – including local language options. The FN 1000 has USB and Ethernet ports for printer and bar code reader and results can be readily copied to a memory stick or sent for external collection e.g. in a central LIMS system.



Features and Benefits		
Faster	User set min FN result auto-stop option for time saving during high sample load periods.	
Safer	Isolated water bath and stirring mechanism, and reduced steam.	
Easy to Use	Auto-start, automated water level control, temperature and atmospheric pressure sensing. Confidently used by non-technical operators.	
Reliable	Simple, robust design provides exceptional instrument life and low cost of ownership.	
Calibration-Free	Saving users time and ensuring correct, reliable measurements.	
Altitude Correction	Automatic re-calculation of FN results. Built-in atmospheric pressure sensor alerts operator if altitude correction is required.	
World Standard	Uniform reporting for grain growers, traders and processors.	
The Only Approved Instruments	Use for trade and export purposes. The Perten Falling Number models are the only validated instruments covered by the approved International standards: AACC/No. 56-81.04, ICC/No. 107/1, ISO/DIS 3093.	



Uses		
Segregation	Save money by avoiding mixing sound and sprouted grain.	
Blend Optimization	Blend grains or flours to create a product with specific characteristics.	
Quality Assurance	Ensure deliveries meet end-user specifications and purchase agreements.	
Fungal FN	Verify total enzymatic activity in flours supplemented with fungal enzymes.	

Recommended Accessories		
Easily and accurately dispenses 25 ml of water.		
Saves water and environment by re-circulating cooling water		
Automatic shaker for fast and uniform sample mixing.		
Rapid Falling Number tube cleaner.		
Approved hammer mills for preparation of grain.		
Compact USB printer for hard copy results.		













Laboratory Mill 3100

Printer

PerkinElmer, Inc. 940 Winter Street Waltham, MA 02451 USA P: (800) 762-4000 or

(+1) 203-925-4602 www.perkinelmer.com

