

Mark 3 LTE Moisture Analyzer

Simple. Dependable. Accurate. Fast.



Out of the box solution at your fingertips!

Fast and accurate moisture or solids measurement with the Mark 3 LTE allows adjustments to be made in the processing cycle in minutes rather than hours. One button operation to perform a test. No pre-programming required on the part of the operator.

Thousands of material program settings already established for the following markets:

- | | |
|------------------|---------------------|
| - Pharmaceutical | - Food |
| - Chemical | - Wastewater |
| - Cosmetics | - Ethanol |
| - Pulp and paper | - Paint and coating |

In-House Testing:

Sartorius Mechatronics will provide program development, free of charge.

Free Trial:

Call for a free two week trial of a Sartorius Mechatronics product at your location.

info@sartorius-omnimark.com
www.sartorius-omnimark.com

Tel (800) 835-3211



Specifications

Method of Measurement	Loss on Drying
Components Analyzer	Modular: Control and Heater modules
Optional Configuration	Up to 2 Heater modules per Control module
Heat Source	Four parallel infrared quartz cylinders
Temperature Steps	Programmable one or two
Temperature Ramp Rates	Two: Normal and Rapid
Temperature Adjustment	Optional fixture for internal electronic two point with programmable prompt
Standby Temperature	30 – 160° C in one degree increments
Equilibration Time	Programmable setting of time at Standby Temperature
Temperature Display	Display of both target and current temp at each stage of drying and standby
Conserve Mode Clock	Programmable time by day of week to shut down and wake up analyzer
Balance Capacity	40 Grams
Balance Resolution	0.001 grams
Balance Repeatability	+/- 1 mg
Balance Calibration	Internal calibration weight of external manual with prompt
Balance Module Design	Thermally isolated with inner and outer stainless steel containers
Balance Spill Protection	Reservoir under draft shield to collect liquids, powders or pellets
Ideal Weight	Audio and visual display for programmable target sample weight addition
Display of Results	0.01%
Recommended Working Range	0.1 to 99.9%
Target Limit Alert	Minimum and maximum programmable limits and screen alert when test done
Program Storage	60 programs with unique settings, alphanumeric naming
Result Storage	999 individual results
Statistical Evaluation	Average, SD, RSD, high and low
Units of Measure of Results	Weight, % Moisture, % Solids, % Volatiles, Fuel Moisture, PPM
Endpoint Determination	Actual slope, Calculation and Time modes
Sample ID	Numeric, Numeric incrementing, alphanumeric or turned off
Modes of Operation	Standard or Concentration modes
Security Mode	Programmable passcode to lock Setup menu
Graph Mode	Graph of Temperature/Time and Percent/Time on display or optional printer
Display	Quarter VGA with contrast adjustment (black on white or white on black)
Keypad	Sealed membrane with audible tactile feedback
Keys	Start/Stop, Paper Feed, numeric, decimal, Enter, 4 navigation and 4 soft keys
Optional Internal Printer	Internal thermal, 40 character graphic
Printout Format	Selectable header (3 lines), analyzer info, operator, Sample ID, Calibration info, program name, program parameters, interval print, weight info, result and signature line
External Input/Output	Serial, USB and Ethernet
Internal Clock	Date and Time: International, Military or AM/PM
Audible Tone	Key stroke, ideal weight met and end of test, selectable On or Off
System Status Indicator Lights	Two tri-color LEDs (balance & heater) on front of heater module
Power Requirements	90-250 VAC 50/60 Hz self adjusting power supply
Dimensions (inches)	16 1/4 x 19 1/2 x 9 1/4 (L x W x H) combined modules
Warranty Period	2 years

The Mark 3 LTE is backwards compatible to many of the standard Mark 2 drying programs



Sartorius Mechatronics Corporation

5 Orville Drive

Bohemia, NY 11716

Phone: 800.645.3108

Fax: 631.254.4252

www.sartorius-mechatronics.com