

Mark 3 LTE Moisture Analyzer

Modular Infrared Moisture Analyzer with Quartz Heater

Advantages

- Economical and reliable alternative to the drying oven
- Intuitive operation for fast learning
- Easy one-button measurement
- Robust and fast, optimal for routine operation
- Multi-module operation for parallel measurements



Product Description

Mark 3 is a flexible moisture analyzer system for labs and production facilities. With its resolution of 0.01%, the Mark 3 LTE has been optimized to achieve excellent reproducibility in routine measurements and meets even highest demands because of its built-in Sartorius weighing technology.

Fast

Quick measuring times give you greater flexibility and let you adapt your production processes fast.

Flexible

Whether your samples are solid, liquid or pasty, the Mark 3 LTE provides optimal flexibility in analyzing a variety of samples with moisture contents ranging from 1.00% to 99.90%. With its optional modular add-ons of up to two heating units, it provides the best solution for every measuring task – regardless of whether you need to perform three or 300 measurements a day.

Easy

Even inexperienced users can reliably operate the Mark 3 intuitively and easily. Thanks to effective password protection, access can be restricted to strictly measuring functions. One touch of a button is all it takes to start a measurement.

Areas of Application

- Foods
- Chemical products
- Biomass & water purification
- Building materials
- Paper, pulp and composites
- Cosmetics
- Paints and coatings
- Agricultural products

Technical Specifications

Display of results	1 mg; 0.01%
Weighing capacity	40 g
Typical repro- ducibility of the weighing system	Starting at an initial sample weight of approx. 1 g: \pm 0.2% Starting at an initial sample weight of approx. 5 g: \pm 0.05%
Method of measurment	Thermogravimetric; loss on drying
Recommended moisture range	1.00 – 99.90%
Heat source	Four parallel infrared quartz cylinders
Modular extension	Optional; up to 2 heater modules connectable
Data printer	Optional internal thermal printer; 40 characters
Temperature range and adjustment	30 – 210°C in one-degree increments
Temperature steps	Programmable; one or two
Standby temperature	Standby temperature selectable from 30 – 160°C
Temperature display	Display of target temperature and actual temperature during the measurement and in standby mode
Display mode	Weight loss in mg % moisture % dry weight % volatiles
Modes of operation	Standard or concentration mode
Endpoint determination	Timer mode; when value falls below a defined weight loss per time unit (%/min); trend shutoff
Program storage	60 programs; battery-backed memory
Data storage	Results with statistics of the last 999 measurements: average, SD, RSD, min and max
Temperature adjustment	Optional set for internal electronic two-point adjustment
Adjustment of the weighing system	Manual calibration and adjustment using an external weight
ldeal initial weight	Audio and visual display for programmable weighing of a sample to an initial target
Initial sample weight lock	Lockable initial sample weight for optimal repeatability
Pass-fail function	Programmable target value, including limits; pass-fail information at the end of a measurement

Sample ID	Numeric, consecutive numbering, alphanumeric or turned off
User management	20 users can be stored with password protection
Password protection	Password protection against undesired adjustment of parameters
Languages	English, French, German, Italian, Portuguese, Spanish
Date and time settings	dd mm yy, mm dd yy, 12 h- or 24 h-format
Saving mode	Integrated timer to switch the analyzer on and off Programmable sleep mode after a longer period of non-use
Display	Backlit VGA display
Parameter and data input	Sealed membrane foil overlay with audible tactile feedback
Real-time curves	Real-time curves of the loss in weight (% or grams) as well of the temperature curves during and after the measurement
Acoustic signals	Key stroke, ideal weight met and end of test; selectable (on or off)
System status indicator lights	Two tricolor LEDs (balance & heater) on front of heater module
Data input output	Serial and Ethernet; USB for service
Printout format (Printer optional)	Configurable printout: selectable header (3 lines), analyzer info, operator, sample ID, calibration info, program name, program parameters, interval print, weight info, result and signature line
Weighing system	Monolithic weighing system for high-precision results
Balance spill protection	Reservoir under draft shield to collect liquids, powders or granules
Operating range	15-30°C ($59-86$ °F) $10-85$ % rel. humidity; non-condensing
Power requirements	90-250 VAC 50 60 Hz, self-adjusting power supply
Housing dimensions $(W \times D \times H)$	$41.3 \times 49.5 \times 23.5$ cm (164" \times 191" \times 94"), modules combined
Weight	10 kg (22 lbs.), modules combined
Warranty	2 years

Order Numbers

LMA110-SET-P	 Mark 3 LTE moisture analyzer with printer, consisting of – LMA100PA control module, incl. printer – LMA110SQ heater module for measurement (1 mg, 40 g)
LMA110-SET	Mark 3 LTE moisture analyzer without printer, consisting of – LMA110SA control module – LMA110SQ heater module for measurement (1 mg, 40 g)
Add-ons: LMA110SQ-000U	Mark 3 LTE heater module (1 mg, 40 g) as a modular add-on to a Mark 3 LTE (max. 2 heater modules)

Equipment Supplied

- Inter-module cable
- Power module cable
- Power cord
- Tweezers
- Sample pans
- Operating instruction manual

Addition for Mark 3 LTE with printer:

- Roll of paper

Accessories

A variety of consumables, cables and inspection, measuring and test equipment are available for the Mark 3 moisture analyzer along with other accessories. We will be glad to advise you in the selection of appropriate accessories.



Mark 3 LTE in parallel operation with 2 measuring modules

Europe

Germany Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen

Phone +49.551.308.0

France & Suisse Romande

Sartorius France 2. rue Antoine Laurent de Lavoisier ZA de la Gaudrée 91410 Dourdan

Phone +33.1.70.62.50.00

Austria

Sartorius Austria GmbH Modecenterstrasse 22 1030 Vienna

Phone +43 1 7965760 0

Belgium

Sartorius Belgium N.V. Rue Colonel Bourg 105 1030 Bruxelles

Phone +32.2.756.06.90

Finland & Baltics

Sartorius Biohit Liquid Handling Oy Lainnatie 1 00880 Helsinki

Phone +358.9.755.951

Hungary

Sartorius Hungária Kft. Kagyló u. 5. 2092 Budakeszi

Phone +3623.457.227

Ireland

Sartorius Ireland Ltd. Unit 41, The Business Centre Stadium Business Park Ballycoolin Road Dublin 11

Phone +353.1.8089050

Italy

Sartorius Italy S.r.l. Via Torino 3/5 20814 Varedo (MB)

Phone +39.0362.5557.11

Netherlands

Sartorius Netherlands B.V.

Phone +31.30.60.53.001

info.netherlands@sartorius.com

Sartorius Poland sp.z o.o. ul. Wrzesinska 70 62-025 Kostrzyn

Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS" Vasilvevsky Island 5th line 70, Lit. A 199178 St. Petersburg

Phone +7.812.327.53.27

Spain & Portugal

Sartorius Spain, S.A. Avda, de la Industria, 32 Edificio PAYMA 28108 Alcobendas (Madrid)

Phone Spain +34.913.586.095 Phone Portugal +351.800.855.800

Switzerland

Sartorius Mechatronics Switzerland AG Ringstrasse 24a 8317 Tagelswangen (ZH)

Phone +41.44.746.50.00

U.K.

Sartorius UK Ltd. Longmead Business Centre Blenheim Road, Epsom Surrey KT19 9QQ

Phone +44.1372.737159

Ukraine

LLS "Sartorius RUS" Post Box 440 "B" 01001 Kiev, Ukraine

Phone +380.44.411.4918

Americas

IISA

Sartorius Corporation 5 Orville Drive, Suite 200 Bohemia, NY 11716

Phone +1.631.254.4249 Toll-free +1.800.635.2906

Argentina

Sartorius Argentina S.A. Int. A. Ávalos 4251 B1605ECS Munro **Buenos Aires**

Phone +54.11.4721.0505

Brazil

Sartorius do Brasil Ltda Avenida Senador Vergueiro 2962 São Bernardo do Campo CEP 09600-000 - SP- Brasil

Phone +55.11.4362.8900

Canada

Sartorius Canada Inc 1173 North Service Road West, D4 Oakville, ON L6M 2V9

Phone +1.905.569.7977 Toll-Free +1.800.668.4234

Mexico

Sartorius de México, S.A. de C.V. Libramiento Norte de Tepotzotlan s/n, Colonia Barrio Tlacateco, Municipio de Tepotzotlan, Estado de México, C.P. 54605

Phone +52.55.5562.1102 leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C. Avenue Alberto del Campo 411 Floor 12 - The Office 15076 - San Isidro, Lima

Phone +51.1.441 0158

Asia | Pacific

Australia

Sartorius Australia Ptv. Ltd. Unit 5, 7-11 Rodeo Drive Dandenong South Vic 3175

Phone +61.3.8762.1800

China

Sartorius (Shanghai) Trading Co., Ltd. 3rd Floor, North Wing, Tower 1 No. 4560 Jinke Road Zhangjiang Hi-Tech Park **Pudong District** Shanghai 201210, P.R. China

Phone +86.21.6878.2300

Hong Kong

Sartorius Hong Kong Ltd. Unit 1012, Lu Plaza 2 Wing Yip Street Kwun Tong Kowloon, Hong Kong

Phone +852.2774.2678

Sartorius Weighing India Pvt. Ltd. #69/2-69/3, NH 48, Jakkasandra, Nelamangala Tg 562 123 Bangalore, India

Phone +91.80.4350.5250

Japan

Sartorius Japan K.K. 4th Fl., Daiwa Shinagawa North Bldg. 8-11, Kita-Shinagawa 1-chome Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.3740.5408

Malaysia

Sartorius Malaysia Sdn. Bhd Lot L3-E-3B, Enterprise 4 Technology Park Malaysia Bukit Jalil 57000 Kuala Lumpur, Malaysia

Phone +60.3.8996.0622

Singapore

Sartorius Singapore Pte. Ltd 10 Science Park Rd The Alpha #02-13/14 Singapore Science Park II Singapore 117684

Phone +65.6872.3966

South Korea

Sartorius Korea Ltd. 8th Floor, Solid Space B/D, PanGyoYeok-Ro 220, BunDang-Gu SeongNam-Si, GyeongGi-Do, 463-400

Phone +82.31.622.5700

Thailand

Sartorius (Thailand) Co. Ltd. 129 Rama 9 Road, Huaykwang Bangkok 10310

Phone +66.2643.8361-6