



Mark 3 HP Moisture Analyzer

Modular Infrared Moisture Analyzer
with Quartz Heater

Advantages

- The alternative to the Karl-Fischer titration and oven drying methods
- Intuitive operation for fast learning
- Easy one-button measurement
- Robust and fast; optimal for routine operation
- Extended multi-module operation for parallel measurements



Product Description

Mark 3 is a flexible moisture analyzer system for labs and production facilities. Its high-resolution weighing system also makes the Mark 3 HP suitable for precisely analyzing the moisture and water content of very dry samples. As a result of this feature, the system is an economical and reliable alternative to Karl-Fisher titration.

Fast

Quick analysis times give you greater efficiency and let you adapt your production processes fast. On top of this, the Mark 3 will even meet the highest requirements of difficult materials. Its integrated two-phase drying allows the combination and separate evaluation of two different temperatures to determine, for example, the surface and total moisture content of a sample in just one measurement.

Reliable

The fully automatic calibration function ensures optimal measurement conditions – every time. In addition, an extensive range of security settings guarantees reliable results and 100% traceability.

Flexible

Whether your samples are solid, liquid or pasty, the Mark 3 HP provides optimal flexibility in analyzing a variety of samples with moisture contents ranging from 0.001% to 99.999%. With its optional modular add-ons of up to four heating units, it provides the best solution for every measuring task, as it enables even parallel

measurements of different samples – 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
- Pharmaceutical products
- Plastics
- Chemical products
- Sugar
- Biomass & water purification
- Building materials
- Paper, pulp and composites
- Cosmetics
- Paints and coatings

Technical Specifications

Display of results	0.1 mg; 0.001%	Sample ID	Numeric, consecutive numbering, alphanumeric or turned off
Weighing capacity	100 g	User management	20 users can be stored with password protection
Typical reproducibility of the weighing system	Starting at an initial sample weight of approx. 1 g: $\pm 0.1\%$ Starting at an initial sample weight of approx. 5 g: $\pm 0.02\%$	Password protection	Password protection against undesired adjustment of parameters
Method of measurement	Thermogravimetric; loss on drying	Languages	English, French, German, Italian, Portuguese, Spanish
Recommended moisture range	0.001 – 99.999%	Date and time settings	dd mm yy, mm dd yy, 12 h- or 24 h-format
Heat source	Four parallel infrared quartz cylinders	Energy-saving mode	Integrated timer to switch the analyzer on and off; programmable sleep mode after a longer period of non-use
Modular extension	Optional; up to 4 heater modules connectable	Display	Backlit VGA display
Data printer	Integrated thermal printer; 40 characters	Parameter and data input	Sealed membrane keypad overlay with audible tactile feedback
Temperature range and adjustment	30 – 210°C in one-degree increments	Real-time curves	Real-time curves of the loss in weight (% or grams) as well of the temperature curve during and after the measurement
Temperature steps	Programmable; one or two with selectable ramp rate	Acoustic signals	Key beeps; signal when ideal sample weight is reached; and end-of-measurement signal; selectable as On Off
Standby temperature	Standby temperature selectable from 30 – 160°C	System status indicator lights	Two tricolor LEDs (balance & heater) on front of heater module
Temperature display	Display of target temperature and actual temperature during the measurement and in standby mode	Data input output	Serial and Ethernet; USB for service
Display modes	Weight loss in mg % moisture % dry weight % volatiles ppm moisture ppm dry weight concentration mg/L	Printout format	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
Modes of operation	Standard, syringe, concentration or 2-phase drying mode	Weighing system	Monolithic weighing system for high-precision results
Endpoint determination	Timer mode; when value falls below a defined weight loss per time unit (%/min); trend shutoff	Balance spill protection	Reservoir under draft shield to collect liquids, powders or granules
Program	300 programs; battery-backed memory	Operating range	15 – 30°C (59 – 86°F) 10 – 85% rel. humidity; non-condensing
Data storage	Results with statistics of the last 999 measurements: average, SD, RSD, min and max	Power requirements	90 – 250 VAC 50 60 Hz, self-adjusting power supply
Temperature adjustment	Optional set for internal electronic two-point adjustment	Housing dimensions (W × D × H)	41.3 × 49.5 × 23.5 cm (164" × 191" × 94"), modules combined
Adjustment of the weighing system	Internal calibration weight or manual calibration and adjustment using an external weight	Weight	10 kg (22 lbs.), modules combined
Ideal initial weight	Audio and visual display for programmable weighing of a sample to an initial target	Warranty	2 years
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		

Order Numbers

LMA100-SET-P	Mark 3 HP moisture analyzer with printer, consisting of <ul style="list-style-type: none">- LMA100PA control module, incl. printer- LMA100PQ heater module for measurement (0.1 mg, 100 g)
Add-ons: LMA100PQ-000U	Mark 3 HP heater module for measurement (0.1 mg, 100 g) as a modular add-on to a Mark 3 HP (max. 4 heater modules)

Equipment Supplied

- Inter-module cable
- Power module cable
- Power cord
- Tweezers
- Sample pans
- Operating instruction manual
- 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 HP in parallel operation with 3 measuring units

Sales and Service Contacts

For further contacts, visit www.sartorius.com

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
Modectcenterstrasse 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
Laippatie 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

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830

Russian Federation

LLC "Sartorius RUS"
Vasilyevsky 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

USA

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 Pty. 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

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
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



◀ www.sartorius.com