

The New MA160 Moisture Analyzer Manages Your Sophisticated Tasks



Fast, Precise and Easy Moisture Analysis

To be able to perform fast and accurate analysis on the widest variety of product samples, you need a reliable instrument that gives you maximum flexibility.

The MA160 uses the thermogravimetric method to determine the moisture content of liquids, pastes and solids – conveniently, reliably and in minimum time. It delivers outstandingly fast, absolutely reproducible results and features an intuitive Method Development Assistant that enables you to develop new methods in just three easy steps. The MA160 gives you the functions you need for professional and secure management of the methods you have created.

High-speed measurements

Reliable Method Development Assistant function

High memory capacity for results and methods

User-friendly operation

Effortless cleaning

Display of the measurement status

Reliable performance testing

Compatibility with Sartorius MA150

MA35 mode

These methods can be stored via an SD card and transferred to other MA160 analyzers. During a measurement, a status light indicates the current progress of a measurement.

At the touch of a button, you can run the built-in performance test with a ReproEasy pad to check the functionality of the MA160 – ensuring flawless performance over the long term.



Applications

With its optimized AURI heating unit, the MA160 delivers high-speed measurements.

The MA160 is ideal for moisture analysis of the widest variety of samples under different conditions. Its Method Development Assistant function enables you to create your own methods for different samples and efficiently manage the methods in a library. This supports your work in the QC lab and in process monitoring.

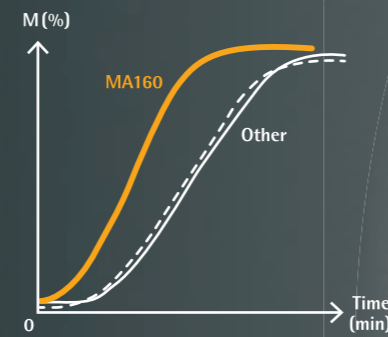
Typical areas of application for the MA160 include moisture analysis of foods, beverages, pharmaceuticals, chemicals, paper materials and products for environmental protection.



Features

High-speed measurements

Both the high-performance AURI heating unit and the sample chamber's geometry ensure rapid heating and fast uniform drying of your sample with up to 30% quicker analysis time.



User-friendly operation

The MA160's intuitive user interface, including a touch screen, and easy-to-understand menu prompts for guidance enable you to easily operate the analyzer – without having to study the manual first.

High memory capacity

MA160 features a high memory capacity, letting you store up to 100 different methods and efficiently manage them in a library. In addition, the results of the last 999 measurements are automatically saved in the data memory, ensuring traceable GMP-compliant documentation at any time.

Fast and precise development of methods

The MA160's Method Development Assistant function solves the time-consuming problem of having to determine and evaluate the right parameters. It enables you to develop new methods quickly in just three simple steps.



Display of the measurement status

The light indicates the current status of a measurement in progress: "running | START," "analysis completed | STOP," "analyzer OFF" or "error."

Reliable performance testing

A ReproEasy Pad in combination with the built-in performance test lets you check the functionality of the entire MA160 moisture analyzer within just 5 minutes. This test guarantees reliable results, saving you from going to the trouble of performing extensive temperature adjustment.

Compatibility with Sartorius MA150

The downward compatibility of the analyzer lets you readily use MA150 method parameters. The additional MA35 mode even enables you to run moisture analysis with parameters developed for the MA35.

Effortless cleaning

The BetterClean design lets you easily and thoroughly clean the instrument parts. Additionally, the heating unit and the sample chamber plate are removable for cleaning in a laboratory washer.



Technical Specifications

Max. weighing capacity	200 g
Reproducibility, typical	Starting at an initial sample weight of approx. 1 g: $\pm 0.2\%$, Starting at an initial sample weight of approx. 5 g: $\pm 0.05\%$
Readability	1 mg, 0.01%
Typical sample quantity	5 g – 15 g
Display modes for results	Moisture content in % M and % g dry weight in % S and g ATRO (ratio) in % M/S
Temperature range and settings	40°C – 200°C, in increments of 1°C Standby temperature selectable from 50°C – 120°C
Sample heating	Infrared heating using an AURI heating unit, 600 W
Heating programs	Standard drying, gentle drying, MA35 mode
Shutoff parameter	Fully automatic, semi-automatic, manual and with timer settings
Forceps for samples	For easy handling of sample pans
Data interface	Mini USB port; automatic printer detection, direct transfer to Microsoft® Office
Data transfer	SD card; function for importing and exporting methods and results
Housing dimensions (W x D x H)	215 x 400 x 210 mm (8.5" x 15.7" x 8.3")
Weight	Approx. 6.2 kg (approx. 13.6 lbs.)

Accessories

6965542	Disposable sample pans, 80 units, aluminum, 90 mm Ø
6906940	Glass fiber pads for analysis of pasty and fatty samples, hard quality; 80 units; 90 mm Ø
6906941	Glass fiber pads for analysis of liquid and fatty samples, soft quality; 200 units; 90 mm Ø
YHP01MA	ReproEasy pads, 10 units for performance testing to check the functionality of the analyzer's heating unit and weighing system
YCW512-AC-02	External calibration weight, 100 g (E2) with DAkkS calibration certificate*
YDP40	Standard laboratory printer
YDP30	Premium GLP laboratory printer
YCC03-D09	Adapter cable for connecting a YDP20-OCE printer
YDP20-OCE	Data printer
YST01MA	Forceps for samples
YMD02B (incl. short report)	Customer-specific service for individual method development at the Sartorius Application Laboratory in Goettingen, Germany, or in the USA

* DAkkS = German accreditation body recognized throughout Europe



Disposable sample pans



ReproEasy pad

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
Fax +49.551.308.3289

France & Suisse Romande

Sartorius France
2, rue Antoine Laurent de Lavoisier
ZA de la Gaudrée
91410 Dourdan
Phone +33.1.70.62.50.00
Fax +33.1.64.59.76.39

Austria

Sartorius Austria GmbH
Modecenterstrasse 22
1030 Vienna
Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Rue Colonel Bourg 105
1030 Bruxelles
Phone +32.2.756.06.90
Fax +32.2.481.84.11

Finland & Baltics

Sartorius Biohit Liquid Handling Oy
Laippatie 1
00880 Helsinki
Phone +358.9.755.951
Fax +358.9.755.95.200

Hungary

Sartorius Hungária Kft.
Kagyó u. 5.
2092 Budakeszi
Phone +3623.457.227
Fax +3623.457.147

Ireland

Sartorius Ireland Ltd.
Unit 41, The Business Centre
Stadium Business Park
Ballycoolin Road
Dublin 11
Phone +353.1.8089050
Fax +353.1.8089388

Italy

Sartorius Italy S.r.l.
Viale A. Casati, 4
20835 Muggiò (MB)
Phone +39.039.4659.1
Fax +39.039.4659.88

Netherlands

Sartorius Netherlands B.V.
Phone +31.30.60.53.001
Fax +31.30.60.52.917
info.netherlands@sartorius.com

Poland

Sartorius Poland sp.z o.o.
ul. Wrzesinska 70
62-025 Kostrzyn
Phone +48.61.6473830
Fax +48.61.6473839

Russian Federation

LLC "Sartorius RUS"
Uralskaya str. 4, Lit. B
199155 St. Petersburg
Phone +7.812.327.53.27
Fax +7.812.327.53.23

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
Fax Spain +34.913.589.623
Fax Portugal +351.800.855.799

Switzerland

Sartorius Mechatronics Switzerland AG
Ringstrasse 24a
8317 Tagelswangen (ZH)
Phone +41.44.746.50.00
Fax +41.44.746.50.50

U.K.

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road, Epsom
Surrey KT19 9QQ
Phone +44.1372.737159
Fax +44.1372.726171

Ukraine

LLS "Sartorius RUS"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918
Fax +380.50.623.3162

Americas

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1.631.254.4249
Toll-free +1.800.635.2906
Fax +1.631.254.4253

Argentina

Sartorius Argentina S.A.
Int. A. Avalos 4251
B1605ECS Munro
Buenos Aires
Phone +54.11.4721.0505
Fax +54.11.4762.2333

Brazil

Sartorius do Brasil Ltda
Avenida Senador Vergueiro 2962
São Bernardo do Campo
CEP 09600-000 - SP- Brasil
Phone +55.11.4362.8900
Fax +55.11.4362.8901

Canada

Sartorius Canada Inc.
2179 Dunwin Drive #4
Mississauga, ON L5L 1X2
Phone +1.905.569.7977
Toll-Free +1.800.668.4234
Fax +1.905.569.7021

Mexico

Sartorius de México, S.A. de C.V.
Libramiento Norte de Tepetzotlan s/n,
Colonia Barrio Tlacateco,
Municipio de Tepetzotlan,
Estado de México,
C.P. 54605
Phone +52.55.5562.1102
Fax +52.55.5562.2942
leadsmex@sartorius.com

Peru

Sartorius Peru S.A.C.
Av. Emilio Cavenecia 264 San Isidro
15073 Lima, Perú
Phone +51.1.441 0158
Fax +51.1.422 6100

Asia | Pacific

Australia

Sartorius Australia Pty. Ltd.
Unit 5, 7-11 Rodeo Drive
Dandenong South Vic 3175
Phone +61.3.8762.1800
Fax +61.3.8762.1828

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
Fax +86.21.6878.2882

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwun Tong
Kowloon, Hong Kong
Phone +852.2774.2678
Fax +852.2766.3526

India

Sartorius Weighing India Pvt. Ltd.
#69/2-69/3, NH 48, Jakkasandra,
Nelamangala Tq
562 123 Bangalore, India
Phone +91.80.4350.5250
Fax +91.80.4350.5253

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
Fax +81.3.3740.5406

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
Fax +60.3.8996.0755

Singapore

Sartorius Singapore Pte. Ltd
1 Science Park Road,
The Capricorn, #05-08A,
Singapore Science Park II
Singapore 117528
Phone +65.6872.3966
Fax +65.6778.2494

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
Fax +82.31.622.5799

Thailand

Sartorius (Thailand) Co. Ltd.
129 Rama 9 Road,
Huaykwang
Bangkok 10310
Phone +66.2643.8361-6
Fax +66.2643.8367



◀ www.sartorius.com

Microsoft® is a registered trademark of Microsoft Corporation, USA.