



MA37

Fully Automatic Infrared Moisture Analyzer
for Daily Routine Operation

Advantages

- High-speed measurements
- Effortless cleaning
- User-friendly operation
- Status light for display of measurement progress



Product Description

The MA37 is the new Sartorius basic analyzer for moisture determination. It operates even faster and is easier to use than its predecessor MA35. In addition, it features a compact footprint and delivers highly accurate results. Moreover, its BetterClean design also enables effortless cleaning, and the convenient status light indicates the current progress of a measurement. The MA37 is therefore ideal for operations involving identical routine measurement tasks, guiding you step by step throughout moisture analysis.

Applications

The MA37 moisture analyzer is designed for quick and reliable determination of the moisture content of liquids, pastes and solids using the thermogravimetric method. The MA37 is perfectly suited to daily recurring moisture analysis tasks, such as in production or incoming goods inspection.

Performance

Fully automatic endpoint determination eliminates the inconvenience of time-consuming screening of shutoff parameters. The moisture analyzer monitors the drying process and ends the measurement once it detects a constant sample weight. The built-in weighing system with 1 mg resolution has been specially optimized for use in higher temperature ranges, delivering the required measurement accuracy. A high-performance metal tube heater heats the samples in the MA37. This heater unit is fast, extremely rugged and durable. Unlike glass heating lamps, such as infrared lamps or halogen heaters, it is especially resistant to dirt and vibration. The functionality of the MA37 can be verified at any time by running the built-in performance test with a ReproEasy pad.

Technical Specifications

Max. weighing capacity	70 g
Reproducibility, typical	For initial sample weights of approx. ≥ 1 g: $\pm 0.2\%$ For initial sample weights 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 metal heating elements
Heating programs	Standard drying, gentle drying, MA35 mode
Shutoff parameter	Choice of modes: – Fully automatic – Semiautomatic mg (1 mg – 50 mg 5 sec. – 300 sec.) – Semiautomatic % (0.1% – 5.0% 5 sec. – 300 sec.) – Timer settings (02:00 – 99.59 min.) – Manual
Access to sample chamber	Removable hood with wide opening angle; SoftClose mechanism
Measuring program	1 program saved in the non-volatile memory (user-definable method parameters)
Memory for data storage	Results are saved until the start of the next measurement
Operator guidance features	Intuitive user interface, including touch screen and easy-to-understand menu prompts
Languages selectable in menu	English set as the default; can be changed in the menu to Chinese, French, Italian, Japanese, Polish, Portuguese, Russian, Spanish and Turkish
Status light	Displays the status "analysis running/START," "analysis completed/STOP" or "analysis error"
Sample inspection	LED-illuminated sample chamber; inspection window with grid above the hood
Draft shield	Integrated draft shield
Cleaning	Removable hood, grid and sample chamber plate for easy cleaning in a laboratory dishwasher
Printout	– Printout using the optional, internal printer YDP30; alternatively, the printer YDP20-OCE can be used with an adapter (YCC03-D09) – GLP-compliant, inalterable standard for printout of the results paper-saving mode for short record
Performance test	Menu function for testing the analyzer's performance using a ReproEasy pad
Monitoring and control of the accuracy of the analyzer as inspection, measuring and test equipment	External calibration using optional calibration weights; Temperature adjustment with optional adjustment set
Data interface	Mini USB port – Automatic detection of the Sartorius printers YDP30 and YDP40 – Direct data transfer to Microsoft® Office programs without any additional software – Programmable data output interval – SBI data transfer protocol
Frequency	50 60 Hz
Power consumption	Max. 640 VA
Temperature range	10°C – 30°C
Housing dimensions (W × D × H)	215 × 400 × 210 mm (8.5" × 15.7" × 8.3")
Weight	Approx. 6.2 kg (approx. 13.6 lbs.)

Available Models

MA37-1	Standard model without country-specific additions and with automatic voltage recognition of 115 V 230 V
MA37-1US	Standard model with country-specific additions for the USA

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
YCW452-AC-02	External calibration weight, 50 g (E2) with DAkkS* calibration certificate
YDP40	Standard printer
YDP30	Premium GLP laboratory printer
YCC03-D09	Adapter cable for connecting the printer YDP20-OCE
YDP20-OCE	Data printer
YTM15MA	Temperature adjustment set
YHP01MA	ReproEasy pads, 10 units for performance testing to check the functionality of the heating unit and weighing system of the analyzer
YST01MA	Forceps for samples
YMD01B (incl. short report)	Customer-specific service for individual method development at the Sartorius Application Laboratory in Goettingen, Germany, or in the USA

Equipment Supplied

Moisture analyzer, power cord, user manual, aluminum sample pans, glass fiber pads – hard quality

* DAkkS = German accreditation body recognized throughout Europe

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.
Kagyló 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