

MA37

Fully Automatic Infrared Moisture Analyzer
for Daily Routine Operation

Advantages

- High-speed measurements
- Effortless cleaning
- User-friendly operation
- Display of process status



Product Description

The MA37 is the new Sartorius basic model for moisture determination. It works 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 allows for effortless cleaning, and the status light indicates the current process status. The MA37 is ideal for operations characterized by consistently routine measurement tasks.

Applications

The MA37 moisture analyzer can be used for quick and reliable determination of the moisture content of liquid, pasty and solid substances using the thermogravimetric method. Typical applications of the MA37 include the analysis of foods, beverages, pharmaceuticals, chemicals, paper materials and environmental protection products. For daily recurring moisture analysis tasks, such as in production or incoming goods inspection, the MA37 is perfect!

Performance

Fully automatic endpoint determination eliminates the cumbersome capture of termination criteria. The moisture analyzer monitors the drying process and stops the measurement once the sample weight is constant. The built-in weighing system with 1 mg resolution is specifically optimized for use in higher temperature ranges and delivers the required measurement accuracy. Two powerful metal tube heating elements heat the samples in the MA37. These heating elements are fast, extremely rugged and durable. Compared to glass heating lamps, e.g. infrared lamps or halogen heaters, they are especially resistant to dirt and vibrations.

Technical Specifications

Max. weighing capacity	70 g
Repeatability, average	for initial sample weight approx. > 1 g: ± 0.2 % for initial sample weight approx. > 5 g: ± 0.05 %
Readability	1 mg, 0.01 %
Typical sample quantity	5 – 15 g
Result display	Moisture content in %M and g dry matter in %S and g ATRO in %M/S
Temperature range and settings	40°C–160°C, stand-by temperature selectable from 40–100°C in increments of 1°C
Sample heating	Infrared metal heating elements
Heating programs	Standard drying, gentle drying
Shutoff parameter	Optional: – Fully automatic – Semiautomatic mg (1–50 mg 5–300 sec.) – Semiautomatic % (0.1–5.0 % 5–300 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 a non-volatile memory (freely selectable 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 guidance
Language selection	English, German, French, Spanish, Italian, Russian, Chinese, Polish, Japanese, Portuguese, 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, inspection window with grid and sample chamber plate for easy cleaning in the dishwasher
Log 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 results printout paper-saving short record
Monitoring of inspection and testing equipment	External calibration using optional calibration weights
Data interface	Mini USB – Automatic detection of the Sartorius printers YDP30 and YDP40 – Direct data transfer to Microsoft® Windows programs without any additional software – Programmable data output interval – Data transfer protocols SBI, table format, text format
Frequency	50/60 Hz
Power consumption	Max. 640 VA
Temperature range	10°C–30°C
Housing dimensions (W × D × H)	215 × 400 × 210 mm
Weight	Approx. 6.2 kg

Available Models

MA37 115V/230V Automatic voltage detection 115V/230V

Accessories

6965542	Disposable sample pans, 80 pcs., aluminum, Ø 90 mm
6906940	Glass fiber pad for analysis of pasty and fatty samples, hard quality, 80 pcs., Ø 90 mm
6906941	Glass fiber pad for analysis of liquid and fatty samples, soft quality, 200 pcs., Ø 90 mm
YCW452-AC-02	External calibration weight, 50 g (E2) with DKD certificate
YDP40	Standard printer
YDP30	Premium GLP laboratory printer
YCC03-D09	Adapter cable for connecting the printer YDP20-OCE

Equipment Supplied

Moisture analyzer, power cord, user manual, aluminum sample pans, glass fiber filter

Sales and Service Contacts

For further contacts, visit www.sartorius.com

Europe

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Weender Landstrasse 94-108
37075 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
Franzosengraben 12
1030 Vienna

Phone +43.1.7965760.0
Fax +43.1.7965760.24

Belgium

Sartorius Belgium N.V.
Leuvensesteenweg, 248/B
1800 Vilvoorde

Phone +32.2.756.06.71
Fax +32.2.253.45.95

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 & Ukraine

LLC "Sartorius ICR"
and LLC "Biohit"
Uralskaya str. 4, Lit. B
199155, Saint-Petersburg

Phone +7.812.327.5.327
Fax +7.812.327.5.323

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine

Phone +380.44.411.4918
Fax +380.50.623.3162

Spain & Portugal

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

Phone Spain +34.902.123.367
Phone Portugal +351.800.855.800
Fax Spain +34.91.358.96.23
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

America

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. Ávalos 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.
Circuito Circunvalación Poniente
No. 149
Ciudad Satélite
53100, Estado de México
México

Phone +52.5555.62.1102
Fax +52.5555.62.2942

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 Scientific
Instruments (Beijing) Co., Ltd.
33 Yu An Road, Airport Industrial Park
Zone B,
Shunyi District, Beijing 101300,
P.R.China

Phone +86.10.8042.6300
Fax +86.10.8042.6486

Hong Kong

Sartorius Hong Kong Ltd.
Unit 1012, Lu Plaza
2 Wing Yip Street
Kwung 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 Stedim Japan K.K.
4th Fl., Daiwa Shinagawa North Bldg.
8-11, Kita-Shinagawa 1-chome
Shinagawa-ku, Tokyo, 140-0001 Japan

Phone +81.3.4331.4300
Fax +81.3.4331.4301

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