

The New Moisture Analyzer – MA37

Fast – Compact – Accurate



Your Reliable Partner for Routine Tasks

MA37 is a next-generation moisture analyzer and continues the success story of the MA35. The MA37 delivers fast, accurate results and is easy to operate. This compact device has a BetterClean design which offers effortless cleaning. The status light makes it easy to track the current status of the measurement.

Therefore, the MA37 is ideal for users who require a fast, reliable moisture analyzer which is easy-to-operate and to clean.

Fast measurements

Display of measurement status

User-friendly operation

Effortless cleaning





Your Reliable Partner for Routine Tasks

MA37 is a next-generation moisture analyzer and continues the success story of the MA35. The MA37 delivers fast, accurate results and is easy to operate. This compact device has a BetterClean design which offers effortless cleaning. The status light makes it easy to track the current status of the measurement.

Therefore, the MA37 is ideal for users who require a fast, reliable moisture analyzer which is easy-to-operate and to clean.

Fast measurements

Due to its high-performance heating elements, the MA37 heats quickly which results in reduced measuring time.

Display of measurement status

The status light indicates the current measurement status – "running | START", "process finished | STOP", "analyzer OFF" or "error". It is visible from a distance of up to 10 meters.

User-friendly operation

The intuitive user interface, including touch-screen and easy-to-understand menu, considerably simplifies operation of the MA37.

Effortless cleaning

BetterClean design enables the user to easily and thoroughly clean the instrument parts. Additionally, the heating module and the sample chamber plate are removable and dishwasher safe.

Fast measurements

Display of measurement status

User-friendly operation

Effortless cleaning



Technical Specifications

Max. weighing capacity	70 g
Repeatability, typical	Starting at an initial sample weight of approx. 1 g: $\pm 0.2\%$ for initial sample weight approx. > 5 g: $\pm 0.05\%$
Readability	1 mg, 0.01%
Typical sample quantity	5 – 15 g
Reading 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, in increments of 1 degree Standby temperature selectable from 40 to 100°C
Sample heating	Infrared heating using a metal tube heater
Heating programs	Standard drying, gentle drying
Shutoff parameter	Fully automatic, semi-automatic, manual and with timer settings
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
Sample inspection	LED-illuminated sample chamber, inspection window with grid above the hood
Interface	Mini USB, automatic printer detection, direct transfer to Microsoft® programs
Housing dimensions (W × D × H)	215 × 400 × 210 mm
Weight	Approx. 6.2 kg

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
YHP01MA	ReproEasy performance test pads, 10 pcs. for performance testing to verify the repeatability of the analyzer
YCW452-AC-02	External calibration weight, 50 g, E2 with DKD certificate
YDP40	Standard data printer
YDP30	Premium GLP laboratory printer
YCC03-D09	Adapter cable for connecting the YDP20-OCE printer

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
Modecenterstrasse 22
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

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

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

Ukraine

LLC "Biohit"
Post Box 440 "B"
01001 Kiev, Ukraine
Phone +380.44.411.4918
Fax +380.50.623.3162

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
Surrey KT19 9QQ
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
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.

Specifications subject to change without notice. Printed in the EU on paper bleached without chlorine. | W
Publication No.: W-1513-e150303 - Order No.: 98649-015-97 - Ver. 03 | 2015