

MA37 Moisture Analyzer

For Daily Routine Operation

Simplifying Progress

SARTURIUS

# Your Reliable Partner for Routine Tasks



#### Fast measurements

The new, high-performance heating unit integrated into the MA37 ensures fast and uniform drying of the sample.

#### Display of the measurement status

The light indicates the status of the analyzer. As a result, you will have everything under control – even from a distance.

#### User-friendly operation

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

#### Effortless cleaning

The BetterClean design makes it so easy to thoroughly clean the parts of the sample chamber. You can even remove the chamber hood and the sample chamber plate to clean them in a laboratory washer.

#### MA35 mode

This mode lets you readily use method parameters already developed for the MA35.

## Technical Specifications

Max. weighing capacity	70 g
Reproducibility, typical	Starting at an initial sample weight of approx. 1 g: ± 0.2% Starting at an initial sample weight of approx. 5 g: ± 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 a metal tube heater
Heating programs	Standard drying, gentle drying, MA35 mode
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 the non- volatile memory (user-definable 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 port; automatic printer detection; direct data transfer to Microsoft® Office programs
Housing dimensions (W × D × H)	215 × 400 × 210 mm (8.5" × 15.7" × 8.3")
Weight	Approx. 6.2 kg (approx. 13.6 lbs.)

#### Accessories

6965542	Disposable sample pans, 80 units, aluminum, Ø 90 mm
6906940	Glass fiber pads for analysis of pasty and fatty samples, hard quality; 80 units; Ø 90 mm
6906941	Glass fiber pads for analysis of liquid and fatty samples, soft quality; 200 units; Ø 90 mm
УНРО1МА	ReproEasy pads, 10 units for perfomance testing to check the functionality of the analyzer's heating unit and weighing system
YCW452-AC-02	External calibration weight, 50 g (E2) with DAkkS calibration certificate*
YDP40	Standard data printer
YDP30	Premium GLP laboratory printer
YCC03-D09	Adapter cable for connecting the YDP20-OCE printer
YDP20-0CE	Data printer
YST01MA	Forceps for samples
YMD01B (incl. short report)	Customer-specific service for individual method development for the MA37 at the Sartorius Application Laboratory in Goettingen, Germany, or in the USA

\*DAkkS = German accreditation body recognized throughout Europe



### Germany

Sartorius Lab Instruments GmbH & Co. KG Otto-Brenner-Strasse 20 37079 Goettingen Phone +49 551 308 0 Sartorius Corporation 565 Johnson Avenue Bohemia, NY 11716

**USA** 

Phone +1 631 254 4249 Toll-free +1 800 635 2906

For further contacts, visit www.sartorius.com