

MA150Q Moisture Analyzer

Fully Automatic Infrared Moisture Analyzer with Coiled Quartz Heating Element (CQR)

Moisture Analyzer



- Time-saving and convenient alternative to a drying oven
- Utmost measurement accuracy and durability "made in Germany"
- 20 program memory modules for quick method changing
- Quick heating and particularly even heat distribution

| | |
|--------------------------------|--|
| Max. weighing capacity | 150 g |
| Repeatability, 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 – 15 g |
| Reading display | % moisture content % dry weight RATIO value mg weight loss g residue g/kg residue g/l residue |
| Temperature range and settings | 40°C–220°C, stand-by temperature selectable from 40–100°C in increments of 1 degree |
| Sample heating | Infrared heating using a coiled quartz heating element (CQR) |
| Heating programs | Standard drying, gentle drying |
| Shutoff criteria | Optional: – Fully automatic – Semi-automatic (1–50 mg/5–300 sec.) – Semi-automatic (0.1–5.0%/5–300 sec.) – Timer setting (0.1–99.9 min.) – Manual |
| Access to sample chamber | Hinged cover |
| Measuring program | 20 programs saved in a non-volatile memory (freely selectable parameters) |
| Data storage | Results are saved until the start of the next measurement |
| Operator guidance features | Menu-driven, alphanumeric dialogue text (5 languages selectable) and symbols |
| Parameter data entry | Via softkeys, menu-specific key allocation |
| Report printout | – Printout using the optional, external printer YDP20-OCE – Configurable GLP-report for measuring results calibration adjustment (5 languages selectable) |



Open sample chamber



Version without glass

| | |
|--------------------------------|---|
| Test equipment monitoring | <ul style="list-style-type: none"> – External calibration using optional weight sets – Automatic temperature calibration with optional temperature adjustment set YTM03MA |
| Safety | Parameter settings password-protected against unauthorized access |
| Voltage frequency | 48-60 Hz |
| Power consumption | Max. 700 VA |
| Operating temperature range | 10°C...30°C |
| Housing dimensions (W × D × H) | 214 × 320 × 180.5 |
| Weight | Approx. 5.5 kg |

Available Models

| | Supply voltage | Design |
|-----------------|------------------|--|
| MA150C-000115V1 | 115 V +15%, –20% | Due to the design of the coiled quartz heating element, this model is not available as a version without glass |
| MA150C-000230V1 | 230 V +15%, –20% | Due to the design of the coiled quartz heating element, this model is not available as a version without glass |

Accessories

| | |
|------------|--|
| 6965542 | Disposable sample pans, 80 units, aluminum, round, Ø 90 mm |
| 6906940 | Fiberglass pad, Ø 90 mm for pasty and fatty samples, hard quality, 80 units |
| 6906941 | Fiberglass pad, Ø 90 mm for liquid and fatty samples, soft quality, with high suction force, 200 units |
| YAT01MA | Single-use pipettes for weighing liquid samples, 500 units |
| YDS04MA | Panel replacement set, aluminum panels for replacing glass panels, to meet FDA HACCP regulations (conversion kit) |
| YSC02 | SartoCollect, software for data communication between moisture analyzer and PC (connecting cable must be ordered separately) Prerequisite: MS Windows 2000 Windows XP Professional, Vista Ultimate, Windows 7 Ultimate. |
| YCW5128-02 | External calibration weight, 100 g (E2) with DKD certificate |
| YTM03MA | Temperature adjustment set with manufacturer certificate for adjustment of the heat source |
| YDP20-OCE | Data printer, suitable for external verification, with date, time, statistics, numerator including cable |
| 6906918 | Ink ribbon for data printer |
| 6906937 | Paper rolls for data printer, 5 rolls, 50 m each |

Equipment Supplied

- Moisture analyzer with data interface
- Power cord
- User manual
- 80 alu sample pans
- In-use dust cover for keypad
- 1 pair of forceps

Sartorius Weighing Technology GmbH
Weender Landstrasse 94–108
37075 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289

info.mechatronics@sartorius.com
www.sartorius-mechatronics.com

Subject to change without notice.
Printed in Germany on paper that has been bleached
without any use of chlorine. | W
Publication No.: W--2031-e11081
Order No.: 98649-012-26