

Cubis[®] MCM1004

Manual Mass Comparator

User Benefits

- Complete mass standard laboratory in a single unit
- Integrated climate sensors for recording all data relevant for determining measurement uncertainty
- Integrated workflow control for efficient and error-free mass comparison
- Fast measurement cycles according to the ABA, ABBA or AB_n...B_nA method

Highlighted Performance Features

- Cubis[®] MSA color touch screen for fast and simple configuration of parameters and workflows
- Sensor-equipped climate module integrated into the draft shield for recording the temperature, humidity and air pressure
- Integrated calibration workflows for ABA, ABBA, AB_n...B_nA cycles to ensure efficient, error-free mass comparison
- Fully integrated function for determining the measurement uncertainty in accordance with OIML and ASTM recommendations
- Filters for optimal adaptation of the mass comparator to ambient conditions
- Monolithic weighing technology
- Additional applications for density determination, statistics and individual identifiers are integrated as standard programs
- Built-in SD card slot for storage and transfer of all data and settings
- For display and evaluation, complete electronics and power supply separated from the weighing system to prevent heat from affecting the results
- Graphical level indicator for interactive user guidance during levelling
- Easy logging of reference weight data
- Continuous weighing range display: any weight between 0 g and the maximum capacity can be displayed
- USB, RS-232C and Ethernet interface ports to integrate the mass comparator into networks or to enable it to communicate with external software via third-party protocols, standardized communication protocols or web services



Technical Specifications

Metrological Specifications

Maximum capacity	1110 g
Application range	0 – 1110 g
Readability	0.1 mg
Repeatability, optimal ¹⁾	0.05 mg
Repeatability, standard E ²⁾	0.07 mg
Repeatability, E 1/10 load ²⁾	0.05 mg
Repeatability standard, F ³⁾	0.2 mg
Electronic weighing range and tare range	610 g
Substitution weights	500 g
Linearity	0.1 mg 600 g
Eccentric load deviation	30 µg mm
Stabilization time	3 s
Cycle time, ABBA in s	90 s

Basic Equipment

Interfaces	RS232C USB LAN
isoCAL	✓
Draft shield	✓
Application programs	Basic weighing, mass unit conversion, individual identifiers, density determination, statistics
Below-comparator weighing port	✓
Air temperature sensor	✓
Air humidity sensor	✓
Air pressure sensor	✓
PC connecting cable	USB

Ambient Conditions

Permissible operating temperature range	10 – 30 °C
Recommended operating temperature	22 °C
Temperature fluctuations	0.3°C/h 0.5°C/12h
Max. air current	< 0.2 m/s
Humidity range	40 – 70 %
Humidity fluctuations	5% 4 h
Power supply	100 – 240 V AC/50 – 60 Hz
Power consumption	< 35 VA

Dimensions

Weighing pan diameter	90 mm
Sample size (D × H)	135 × 140 mm
Weigh cell (W × D × H)	222 × 431 × 301 mm
Electronic unit (W × D × H)	239 × 320 × 56 mm
Net weight	16.3 kg
Gross weight	31.5 kg
Number of packages	1
Packaging data 1	87 × 60 × 96 cm
Pallet	84 × 60 × 95 cm
Optimal height for setup	800 mm

The standard deviation "s" is the repeatability calculated from 5 ABA cycles under the following conditions:

- 1) Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone no drafts from above
- 2) Standard conditions E: measured by hand in a laboratory under E1 conditions, on a decoupled weighing stone; no drafts from above
- 3) Standard conditions F: measurement performed manually in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above

Applications

OIML calibration range RS	-
OIML calibration range E1	1 kg
OIML calibration range E2	500 g - 1 kg
OIML calibration range F1	100 g - 1 kg
OIML calibration range F2	5 g - 1 kg
OIML calibration range M1	100 mg - 1 kg
OIML calibration range M2	100 mg - 1 kg
OIML calibration range M3	1 g - 1 kg
ASTM E617 calibration range Class 000	-
ASTM E617 calibration range Class 00	1 kg
ASTM E617 calibration range Class 0	300 g - 1 kg
ASTM E617 calibration range Class 1	200 g - 1 kg
ASTM E617 calibration range Class 2	100 g - 1 kg
ASTM E617 calibration range Class 3	30 g - 1 kg
ASTM E617 calibration range Class 4	10 g - 1 kg
ASTM E617 calibration range Class 5	1 g - 1 kg
ASTM E617 calibration range Class 6	10 mg - 1 kg
ASTM E617 calibration range Class 7	10 mg - 1 kg

Optional Accessories

External calibration weight	1 kg E2 YCW552-02
Climate module, uncalibrated, for all MCM models	YCM20MC
Calibration of a YCM20MC climate module with DAkkS calibration certificate	YCM20DAkkS
Climate module with DAkkS calibration certificate for all MCM models	YCM20MC-DAkkS
Optional secondary draft shield	YDS24C
Weighing table	YWT03

Germany

Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Straße 20
37079 Göttingen
Phone +49 551 308 0

USA

Sartorius Corporation
565 Johnson Avenue
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906

 For further information, visit
www.sartorius.com