SARTURIUS

Cubis® MCM106

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,...B, A 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, AB₁...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
- For display and evaluation, complete electronics and power supply separated from the weighing system to prevent heat from affecting the results



- 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
- Graphical level indicator for interactive user guidance during levelling
- Easy logging of reference weight data
- Continuous weighing range display: any weight between
 O 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	111 g
Application range	0 – 111 g
Readability	1 μg
Repeatability, optimal 1)	1 μg
Repeatability, standard E ²)	2 μg
Repeatability, E 1/10 load 2)	1 μg
Repeatability standard, F ³)	5 μg
Electronic weighing range and tare range	61 g
Substitution weights	50 g
Linearity	8 µg
Eccentric load deviation	1μg mm
Stabilization time	5 s
Cycle time, ABBA in s	90 s

Basic Equipment	
Interfaces	RS232C USB LAN
isoCAL	\checkmark
Draft shield	\checkmark
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% 4h
Power supply	100-240 V AC/50-60 Hz
Power consumption	<35 VA

Dimensions	
Weighing pan diameter	50 mm
Sample size (D × H)	50 × 120 mm
Weigh cell (W × D × H)	222 × 431 × 301 mm
Electronic unit (W × D × H)	239 × 320 × 56 mm
Net weight	16 kg
Gross weight	31.3 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:

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 mannually 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	5 g - 100 g
OIML calibration range E1	50 mg-100 g
OIML calibration range E2	1 mg - 100 g
OIML calibration range F1	1 mg - 100 g
OIML calibration range F2	1 mg - 100 g
OIML calibration range M1	1 mg - 100 g
OIML calibration range M2	100 mg - 100 g
OIML calibration range M3	1 g - 100 g
ASTM E617 calibration range Class 000	1 g - 100 g
ASTM E617 calibration range Class 00	1 g - 100 g
ASTM E617 calibration range Class 0	0.05 mg - 100 g
ASTM E617 calibration range Class 1	0.1 mg - 100 g
ASTM E617 calibration range Class 2	0.2 mg - 100 g
ASTM E617 calibration range Class 3	0.3 mg - 100 g
ASTM E617 calibration range Class 4	0.5 mg - 100 g
ASTM E617 calibration range Class 5	0.5 mg - 100 g
ASTM E617 calibration range Class 6	0.5 mg - 100 g
ASTM E617 calibration range Class 7	10 mg - 100 g

Optional Accessories	
External calibration weight	20 g E2 YCW422-02
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

For further information, visit www.sartorius.com

USA

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