

Cubis[®] MCM60K2

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
- External sensor-equipped climate module 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
- 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 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	64 kg
Application range	0 – 64 kg
Readability	10 mg
Repeatability, optimal ¹⁾	6 mg
Repeatability, standard E ²⁾	10 mg
Repeatability standard, F ³⁾	25 mg
Electronic weighing range and tare range	64 kg
Linearity	50 mg
Eccentric load deviation	3.5 mg mm
Stabilization time	5 s
Cycle time, ABBA in s	120 s

Basic Equipment

Interfaces	RS232C USB LAN
Application programs	Basic weighing, mass unit conversion, individual identifiers, density determination, statistics
Below-comparator weighing port	With optional accessory
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 dimensions (W × D)	400 × 300 mm
Weigh cell (W × D × H)	400 × 326 × 126 mm
Electronic unit (W × D × H)	239 × 320 × 56 mm
Net weight	15 kg
Gross weight	18.6 kg
Number of packages	1
Packaging Data	73 × 60 × 36 cm
Optimal height for setup	800 mm

Applications

OIML calibration range RS	-
OIML calibration range E1	-
OIML calibration range E2	50 kg
OIML calibration range F1	20 kg – 50 kg
OIML calibration range F2	5 kg – 50 kg
OIML calibration range M1	2 kg – 50 kg
OIML calibration range M2	500 g – 50 kg
OIML calibration range M3	200 g – 50 kg
ASTM E617 calibration range Class 000	-
ASTM E617 calibration range Class 00	-
ASTM E617 calibration range Class 0	50 kg
ASTM E617 calibration range Class 1	20 kg – 50 kg
ASTM E617 calibration range Class 2	20 kg – 50 kg
ASTM E617 calibration range Class 3	10 kg – 50 kg
ASTM E617 calibration range Class 4	5 kg – 50 kg
ASTM E617 calibration range Class 5	2 kg – 50 kg
ASTM E617 calibration range Class 6	1 kg – 50 kg
ASTM E617 calibration range Class 7	100 g – 50 kg

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

Optional Accessories

External calibration weight	50 kg E2 YCW752-02
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
Tower for climate module, for connection to MCM high-capacity models, including cable	YCM20MC-Tower
Optional draft shield	YDS05C YDS03C
Hook for below-comparator weighing	69EA0040
Lifting device for 10 kg	YAW51
Lifting device for 20 kg	YAW52
Lifting device for 50 kg	YAW53
Centermatic levelling pan	YWP03C

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