

Cubis® II MCA36MC

Manual Mass Comparator



User Benefits

- Integrated workflow control for efficient and error-free mass comparison
- Fast measurement cycles according to the ABA, ABBA or AB₁...B_nA method

Highlighted Features

- Cubis® II MCA color touch screen for fast and simple configuration of parameters and workflows
- 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
- Additional applications for density determination, statistics and individual identifiers are integrated as standard programs
- Automatic Motorized Leveling
- Easy logging of reference weight data
- Continuous weighing range display: any weight between 0 g and the maximum capacity can be displayed

Technical Specifications

Metrological Specifications

Maximum capacity	32 g
Application range	0 – 32 g
Electronic weighing range and tare range	32 g
Readability	1 µg
Repeatability, typical	1 µg
Repeatability, at nominal load (5 × ABA)	1.5 µg (20 g)
Repeatability, at low load (5 × ABA)	0.7 µg (2 g)
Linearity, tolerance	6 µg
Eccentric load deviation, typical	0.64 µg/mm (10 g)
Stabilization time, typical	3.5 s

Basic Equipment

Interfaces	USB-A, USB-B and USB-C
isoCAL	Yes
Draft shield	Large + Internal motorized draft shield
Below-comparator weighing port	Yes
PC connection cable	USB-A
Internal motorized draft shield	Yes

Ambient Conditions

Permissible operating temperature	5 – 40 °C
Recommended operating temperature	22 °C
Temperature fluctuations	0.3 °C/h 0.5 °C/12 h
Max. air current	<0.2 m/s
Humidity range	40 – 80%
Humidity fluctuations	5% 4 h
Power supply	100 – 240 V; 50 – 60 Hz
Power consumption	64.5 W

Dimensions

Weighing pan diameter	50 mm
Inner draft shield dimensions (D × H)	80 × 115 mm
Dimensions (L × W × H)	501 × 240 × 301 mm
Dimensions (L × W × H) with control unit removed	376.5 × 240 × 301 mm
Weight (approx.)	16 kg
Optimal height for setup	80 cm

Applications

OIML calibration range RS	20 g
OIML calibration range E1	50 mg – 20 g
OIML calibration range E2	1 mg – 20 g
OIML calibration range F1	1 mg – 20 g
OIML calibration range F2	1 mg – 20 g
OIML calibration range M1	1 mg – 20 g
OIML calibration range M2	100 mg – 20 g
OIML calibration range M3	1 g – 20 g
ASTM E617 calibration range Class 000	1 g – 30 g
ASTM E617 calibration range Class 00	1 g – 30 g
ASTM E617 calibration range Class 0	0.05 mg – 30 g
ASTM E617 calibration range Class 1	0.1 mg – 30 g
ASTM E617 calibration range Class 2	0.2 mg – 30 g
ASTM E617 calibration range Class 3	0.3 mg – 30 g
ASTM E617 calibration range Class 4	0.5 mg – 30 g
ASTM E617 calibration range Class 5	0.5 mg – 30 g
ASTM E617 calibration range Class 6	0.5 mg – 30 g
ASTM E617 calibration range Class 7	10 mg – 30 g

Optional Accessories

External calibration weight	20 g E2 YCW422-02
Climate module, uncalibrated	YCM20MC
Calibration of a YCM20MC climate module with DAkkS calibration certificate	YCM20DAkkS
Climate module with DAkkS calibration certificate	YCM20MC-DAkkS
Tower for climate module, for mounting YCM20MC	YCM20MC-Tower
Weighing table	YWT03

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
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**
[sartorius.com](https://www.sartorius.com)