

Cubis® II MCA66MC

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 | 61 g |
| Application range | 0 – 61 g |
| Electronic weighing range and tare range | 61 g |
| Readability | 1 µg |
| Repeatability, typical | 1 µg |
| Repeatability, at nominal load (5 × ABA) | 2 µg (50 g) |
| Repeatability, at low load (5 × ABA) | 0.7 µg (2 g) |
| Linearity, tolerance | 8 µg |
| Eccentric load deviation, typical | 0.8 µg/mm (20 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 | 50 g |
| OIML calibration range E1 | 50 mg – 50 g |
| OIML calibration range E2 | 1 mg – 50 g |
| OIML calibration range F1 | 1 mg – 50 g |
| OIML calibration range F2 | 1 mg – 50 g |
| OIML calibration range M1 | 1 mg – 50 g |
| OIML calibration range M2 | 100 mg – 50 g |
| OIML calibration range M3 | 1 g – 50 g |
| ASTM E617 calibration range Class 000 | 1 g – 50 g |
| ASTM E617 calibration range Class 00 | 1 g – 50 g |
| ASTM E617 calibration range Class 0 | 0.05 mg – 50 g |
| ASTM E617 calibration range Class 1 | 0.1 mg – 50 g |
| ASTM E617 calibration range Class 2 | 0.2 mg – 50 g |
| ASTM E617 calibration range Class 3 | 0.3 mg – 50 g |
| ASTM E617 calibration range Class 4 | 0.5 mg – 50 g |
| ASTM E617 calibration range Class 5 | 0.5 mg – 50 g |
| ASTM E617 calibration range Class 6 | 0.5 mg – 50 g |
| ASTM E617 calibration range Class 7 | 10 mg – 50 g |

Optional Accessories

| | |
|--|---------------------|
| External calibration weight | 50 g E2 YCW452-02 |
| 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)