SVISCISVS

Product Datasheet

Cubis® MCM6.7

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 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	6.1 g
Application range	0-6 g
Readability	0.1 µg
Repeatability, optimal ')	0.15 µg
Repeatability, standard E ²)	0.3 µg
Repeatability, E ¼ load ²)	0.2 µg
Repeatability standard, F ³)	0.6 µg
Electronic weighing range and tare range	6.1 g
Linearity	1 µg
Eccentric load deviation	0.25 µg mm
Stabilization time	10 s
Cycle time, ABBA in s	90 s

Dimensions	
Weighing pan diameter	16 mm
Sample size (D × H)	16 × 70 mm
Weigh cell (W × D × H)	122 × 343 × 141 mm
Electronic unit (W × D × H)	239 × 320 × 56 mm
Net weight	7.6 kg
Gross weight	11.9 kg
Number of packages	1
Packaging data 1	76 × 60 × 50 cm
Optimal height for setup	800 mm

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

5 g E2 YCW352-02
YCM20MC
YCM20DAkkS
YCM20MC-DAkkS
YDS20C
YWT03

Basic Equipment	
Interfaces	RS232C USB LAN
isoCAL	\checkmark
Draft shield	\checkmark
Application programs	Basic weighing, mass unit conver- sion, individual identifiers, density determination, statistics
Below-comparator weighing port	\checkmark
Air temperature sensor	\checkmark
Air humidity sensor	\checkmark
Air pressure sensor	\checkmark
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

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) Standards 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

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