



# Mass Metrology

Professional Equipment  
for Absolute Precision

Simplifying Progress

**SARTORIUS**



## From the Prototype Kilogram to Mass Standards

International trade requires the worldwide standardization of measurements. Mass plays an important role because the majority of commerce throughout the world is defined by the mass of substances. To make sure the same masses are used around the world, each country has a National Metrology Institute (NMI) that governs units of measurement.

Sartorius develops innovative mass comparators to the highest standards for NMIs and other metrological institutions. The last few years alone have seen the following developments with customer cooperation:

- 1 kg prototype mass comparator with an incredibly high resolution of 0.0000001 g,
- Various automatic balances,
- Fully automatic systems for determining the volumes of weights,
- Many more products at the forefront of technology.

At this point, we would like to especially thank all our partners for their outstanding cooperation in helping to make our developments exceptionally successful:

- The Bureau International des Poids et Mesures (BIPM),
- The German Physikalisch-Technische Bundesanstalt (PTB),
- The Technical University of Ilmenau (TUI), Germany.





# Products for the Determination of Mass



# Contents


<b>Cubis® MCM Manual Mass Comparators</b> .....	<b>7</b>
The Cubis® MCM Manual Mass Comparators .....	8
Cubis® MCM Performance Features .....	10
Cubis® MCM Models up to 1 kg .....	12
Cubis® MCM Models up to 10 kg .....	13
Cubis® MCM Models up to 60 kg .....	13
Automatic Buoyancy Correction .....	14
<b>Manual High-Capacity Mass Comparators</b> .....	<b>17</b>
Models up to 3,000 kg .....	18
<b>Table-Top Robotic Mass Comparators CCR10-C</b> .....	<b>22</b>
Model CCR10.7-C, range 0.1 µg / 10 g .....	25
Model CCR6.7-C, range 0.1 µg / 6 g .....	25
Model CCR10.6-C, range 1 µg / 10 g .....	25
Key Accessories .....	25
<b>Automatic Mass Comparators and Robots</b> .....	<b>26</b>
Mass Comparator CCL1007 for National Prototype Kilogram .....	28
Robot for Fully Automated Determination of Mass from 1 mg to 1 kg .....	32
<b>Metrological Equipment and Accessories</b> .....	<b>35</b>
Climate Station & Data Logger .....	36
Susceptometer & Permeability Indicator – A Complete Solution for Testing the Magnetic Properties of Weights .....	38
Volume & Density Determination .....	40
ScalesNet-M – The Individual Software Solution for Professionally Calibrating Weights ...	42
Weights and Weight Sets .....	46
<b>Detailed Technical Specifications</b> .....	<b>50</b>
Cubis® MCM Manual Mass Comparators up to 1 kg .....	52
Cubis® MCM Manual Mass Comparators 2 kg to 10 kg .....	54
Cubis® MCM Manual Mass Comparators 40 kg to 60 kg .....	55
Manual High-Capacity Mass Comparators 100 kg to 300 kg .....	56
Manual High-Capacity Mass Comparators 600 kg to 3,000 kg .....	57
Table-Top Robotic Mass Comparators .....	58
Floor-Standing Robotic Mass Comparators .....	60
Mass Comparators with Vacuum Chamber .....	61
Devices for Determination of Volume, Density, Susceptibility and Magnetism .....	62
<b>Calculations and Areas of Application</b> .....	<b>65</b>
Areas of Application in Accordance with OIML R111-1:2004 .....	66
Areas of Application in Accordance with ASTM E617 .....	68
Calculating Uncertainties .....	70





# Cubis® MCM Manual Mass Comparators

- 8 The Cubis® MCM Manual Mass Comparators
- 10 Cubis® MCM Performance Features
- 12 Cubis® MCM Models up to 1 kg
- 13 Cubis® MCM Models up to 10 kg
- 13 Cubis® MCM Models up to 60 kg
- 14 Automatic Buoyancy Correction

 Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# The Cubis® MCM Manual Mass Comparators

## Your Full-range Mass Standards Laboratory

The manual Cubis® MCM mass comparators are the first devices on the market that unite metrological weighing competence with an integrated control of workflows in line with the recommendations of the International Organization of Legal Metrology (OIML). In its international recommendations R111-1, the OIML defines metrological and technical requirements. Its primary aim is the global harmonization of mass determination. Especially the pharmaceutical industry demands that greater accuracies based on global regulations be transferred to production. As an alternative, the Cubis® MCM produces results that are all naturally also ASTM-compliant.

## Integrated Workflow Control

The integrated workflow control of the manual Cubis® MCM mass comparators minimizes the error rate during operation: During the measurement process, the device gives the user instructions about the step to perform next. This markedly keeps the factor “human” from compounding the accuracy of mass determination and makes the results more reliable. At the same time, the workflow is ergonomically optimized and gives the user a more relaxed working atmosphere.

## Integrated Climate Sensors

The sensors integrated in the mass comparator automatically log climate data like temperature, air pressure and humidity for calculating the air buoyancy correction at the site of the measurement.

The climate data can be documented via a PC. That way, it is possible at any time to check that the limits on temperature, air pressure and humidity of the respective calibration levels (E1, E2, F1 or F2) are in compliance.

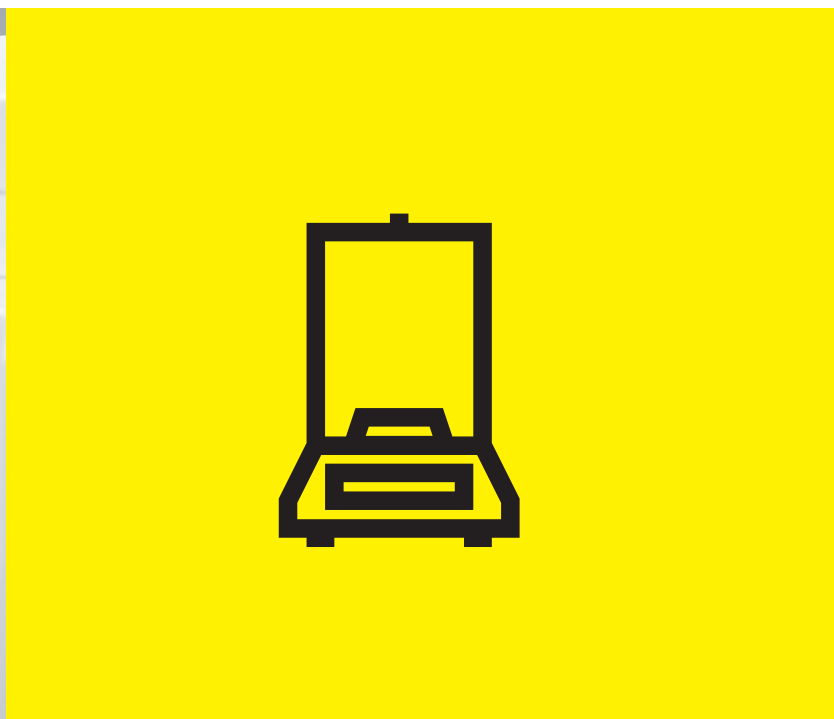
## Fastest Methodology

Compared to conventional devices, Cubis® mass comparators deliver fast-run method cycles (ABA, ABBA or AB1...BnA) for determining the conventional value of mass (also referred to as conventional mass and its combined standard uncertainty).

The manual mass comparators can seamlessly be integrated in the infrastructure of mass standards laboratory. Based on the Cubis® Q-Com communication concepts, they can be integrated in existing networks and every type of data desired can be transferred to other devices.

The Cubis® MCM mass comparators are specified under both ideal and real laboratory conditions. This ensures that they always provide you their full and reliable performance during on-site use.

With all their integrated functions and technical possibilities, the Cubis® mass comparators work like “small metrological laboratories” – but integrated in the mass comparator.







SARTORIUS

MASS COMPARATOR  
MCM 2004

Benno Gatzemeier  
24/04/2014 12:49:38  
d = 0.0001 g  
Max 2500 g  
**+ 1999.9983**  
isoCAL 0% 100%  
Current: A-B-A  
Cycle: 2 / 3  
Measuring test weight . Close Draftshield 3 sec  
Stabilization time:  
Abort

# Cubis® MCM

## Manual Mass Comparators with the Full Performance Spectrum of a Cubis® Balance

- All MCM mass comparators feature a Cubis® MSA display and control unit with color touchscreen for fast and simple configuration of parameters and workflows.
- Thanks to the continuous weighing value display, every weight value between 0 g and maximum capacity can be indicated.
- The climate sensors integrated into the draft shield log temperature, humidity and air pressure (models without a draft shield contain an external climate module with the corresponding sensors in the equipment supplied).
- Steps for calibration performed in accordance with the methods ABA, ABBA, AB1... BnA enable error-free, efficient work.
- The reference weight data and the guidance through the measurement procedure are automatically accounted for.
- The measurement uncertainty is determined in full accordance with OIML and ASTM recommendations. No external software or a climate measuring station is required for recording ambient parameters.
- Filters can be optimally adapted to the ambient conditions.



- The electronics and the power supply for display and evaluation are separate from the weighing system, in order to prevent heat from affecting the results.
- Application programs for weighing, unit conversion, individual identifiers, density determination, statistics are integrated as standard on Cubis® MCM mass comparators.
- MCM mass comparators automatically detect when the balance is skewed and provide graphic support during leveling. On the models MCM2004, MCM5004 and MCM5003, motorized leveling is also possible.
- The Cubis® MCM balances know no limits in terms of connectivity and communication. Diverse data interfaces like USB, RS-232C and Ethernet enable nearly all forms of bidirectional communication. Integration in networks or communication with external software via standardized communication protocols, SICS or Webservices is also possible.
- It is easy to store all data on an SD card and transfer them to PCs or other MCM mass comparators.
- The modern Q-App programming featured by MCM mass determination software allows you to configure the system specifically for individual customers' demands for mass determination or integration into their system. Please contact our product specialists.
- Additional draft shields are available for all models. They reduce air movement caused by air conditioning systems. Even under unfavorable ambient conditions, they keep standard deviations at a minimum.



# Models

Up to 1 kg



<b>Model 1</b>	<b>MCM6.7</b>
Maximum capacity and electrical weighing range	6.1 g
Readability	0.1 µg
Repeatability s under optimal conditions <sup>1)</sup>	0.15 µg



<b>Model 2</b>	<b>MCM36</b>	<b>MCM66</b>	<b>MCM106</b>
Maximum capacity and electrical weighing range	31 g	61 g	111 g   61 g
Readability	1 µg	1 µg	1 µg
Repeatability s under optimal conditions <sup>1)</sup>	1 µg	1 µg	1 µg



<b>Model 3</b>	<b>MCM605</b>	<b>MCM1005</b>	<b>MCM1004</b>
Maximum capacity and electrical weighing range	610 g	1,110 g   610 g	1,110 g   610 g
Readability	10 µg	10 µg	100 µg
Repeatability s under optimal conditions <sup>1)</sup>	10 µg	15 µg	50 µg



Up to 10 kg



Model 4	MCM2004	MCM5004	MCM5003
Maximum capacity and electrical weighing range	2.5 kg	5.1 kg	5.1 kg
Readability	0.1 mg	0.1 mg	1 mg
Repeatability s under optimal conditions <sup>1)</sup>	0.05 mg	0.3 mg	0.5 mg



Model 5	MCM10K3
Maximum capacity and electrical weighing range	11 kg
Readability	1 mg
Repeatability s under optimal conditions <sup>1)</sup>	0.8 mg

Up to 60 kg



Model 6	MCM40K3	MCM60K3	MCM60K2
Maximum capacity	41 kg	64 kg	64 kg
Readability	1 mg	2 mg	10 mg
Repeatability s under optimal conditions <sup>1)</sup>	2 mg	4 mg	6 mg



Model 7	MCM32002
Maximum capacity	32 kg
Readability	10 mg
Repeatability s under optimal conditions <sup>1)</sup>	10 mg

Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift  
<sup>1)</sup> Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above.

## Accessories for Cubis® MCM for Recording Temperature, Air Pressure and Humidity\*

Accessories and Service	Order Number
Climate module, without calibration certificate, for all MCM models	YCM20MC
Calibration of a climate module YCM20MC with DAkKS calibration certificate	YCM20DAkKS
Climate module with DAkKS calibration certificate for all MCM models	YCM20MC-DAkKS
Hook for below-balance weighing for models MCM40K3, MCM60K3, MCM60K2, MCM40K3-DAkKS, MCM60K3-DAkKS and MCM60K2-DAkKS	69EA0040
Tower for climate module, for mounting YCM20MC, can be ported to following models MCM10K3, MCM40K3, MCM60K3, MCM60K2, MCM10K3-DAkKS, MCM40K3-DAkKS, MCM60K3-DAkKS and MCM60K2-DAkKS incl. connection cable	YCM20MC Tower

\* Not available as an option on model MCM32002

# Automatic Buoyancy Correction

When comparing mass under atmospheric conditions, the weights are subject to buoyancy as a function of their volume and the air density; this buoyancy runs in the opposite direction of the weight force. If the material density of the object to be tested deviates from the reference weight, a buoyancy correction as well as a standardization to air density  $1.2 \text{ kg/m}^3$  must be performed to determine the conventional mass.

All models in the Cubis® MCM manual mass comparator family, except model MCM32002, are standard equipped with a module for recording air temperature, air pressure and humidity to determine air density. On comparators equipped with a draft shield, this module is located within the weighing chamber in order to record the ambient conditions that actually prevail. On devices without a draft shield, the climate module is attached to an external tower in order to record ambient conditions relevant to the mass comparison in this case as well.

Since the climate module can be plugged into the mass comparator, it is simply removed for calibration. The calibration data and related characteristic correction curves are stored in the internal memory of the climate module. That means that synchronous and metrologically traceable climate data and their measurement uncertainties are available for mass comparisons. The following standard uncertainties are specified for a calibrated climate module:

## Temperature:

$u_t = 0.15 \text{ K}$  in the range  $18^\circ\text{C}$  to  $27^\circ\text{C}$

## Pressure:

$u_p = 1 \text{ hPa}$  in the range  $800 \text{ hPa}$  to  $1,100 \text{ hPa}$

## Humidity:

$u_{hr} = 1\%$ ,  $30\% < hr < 70\%$

$u_{hr} = 2\%$ ,  $70\% < hr < 90\%$

The mass comparator calculates the air density from the measured ambient conditions and their uncertainty. When density of the test weight and reference weight and their uncertainties are known, the integrated application software "Mass Calibration" calculates the air buoyancy correction including the related uncertainty. The mass comparison results indicate the conventional mass of the calibration weight with the corresponding uncertainties.



External climate module with YCM20MC Tower



Climate module connected within the draft shield






# Manual High-Capacity Mass Comparators

18 Manual High-capacity Mass Comparators  
up to 3,000 kg



 **Find out more**  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# Manual High-Capacity Mass Comparators

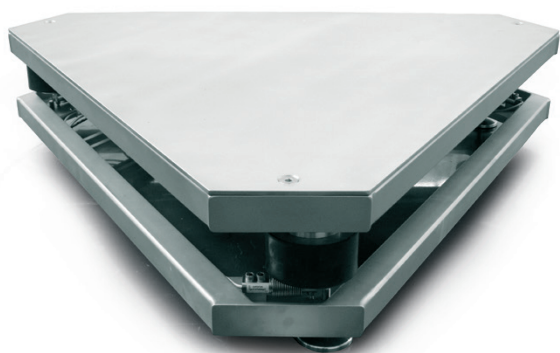
Up to 3,000 kg

All Sartorius high-capacity mass comparators are constructed of high-grade stainless steel. This design eliminates the possibility of reactions from magnetic weights having an effect on comparison weighing.

The unique sandwich construction with integrated, strain-free weigh cells guarantees excellent repeatabilities even when weights are loaded in a "heavyhanded" way.

Models of the CCS range are equipped with four high-resolution strain-test medium type load cells which are aligned with one other. Draft shields are supplied as standard for all high-capacity mass comparators.

The innovative sturdy three-point frame of the CCT models with three high-resolution strain-test medium type load cells guarantee a stable and distortion-free setup. Potential off-center loading errors are minimized due to the large distances between the weigh cells and a very high repeatability is enabled.



## CCT Range

	CCT1000K	CCT2000K
Maximum capacity	1,200 kg	2,100 kg
Readability	1 g	1 g
Repeatability (optimal) s*	2 g	5 g



## CCS Range

	CCS600K	CCS1000K	CCS3000K
Maximum capacity	605 kg	1,510 kg	3,010 kg
Readability	1 g	5 g	10 g
Repeatability (optimal) s*	2 g	5 g	10 g

## CCI Range

	CCI100K2	CCI300K
Maximum capacity	151 kg	303 kg
Readability	50 mg	1 g
Repeatability (typical) s*	200 mg	0.5 g

s\* Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift. Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an M1 mass standards laboratory.




Manual high-capacity mass comparator of the CCI range

# Table-Top Robotic Mass Comparators CCR10-C

- 25 Model CCR10.7-C, range 0.1  $\mu\text{g}$  / 10 g
- 25 Model CCR6.7-C, range 0.1  $\mu\text{g}$  / 6 g
- 25 Model CCR10.6-C, range 1  $\mu\text{g}$  / 10 g
- 25 Key Accessories



 Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)



SARTORIUS

SARTORIUS



# Table-Top Robotic Mass Comparators CCR10-C

Accurate yet compact and cost efficient solutions for mass determination and dissemination processes

The CCR-Compact Robotic Systems offer table-top sized automated solutions for mass determination or dissemination processes up to 10 g maximal capacity. With 120 magazine positions, these systems offer a walkaway solution allowing continuous processing of the measurements overnight or over the weekend. Designed for simultaneous management of reference and test weights the Sartorius dual robotic arm system with the patented multi and single weight handlers speeds up the processes with its most efficient motion sequences and handles the weights with highest precision and consistency. The climate sensors integrated into the weighing chamber additionally contribute to the reliability of the results. Automated weight handling at the range of 1 mg to 10 g not only eliminate the error rates of manual handling but significantly increase productivity which consequently leads to cost-savings long term.

## Your Benefits

- 120 magazine positions allow overnight and weekend processing
- Dual Robotic Arm Technology guarantees the most efficient motion sequences available on the market
- Table-Top design saves prestigious laboratory space
- Robustness assure best repeatability values
- Build-in Climate Sensor directly in the weighing cells to further support measurement accuracy
- 3 Models fulfill your workflow requirements

## High-throughput Allows Weekend Processing and Releases Resources

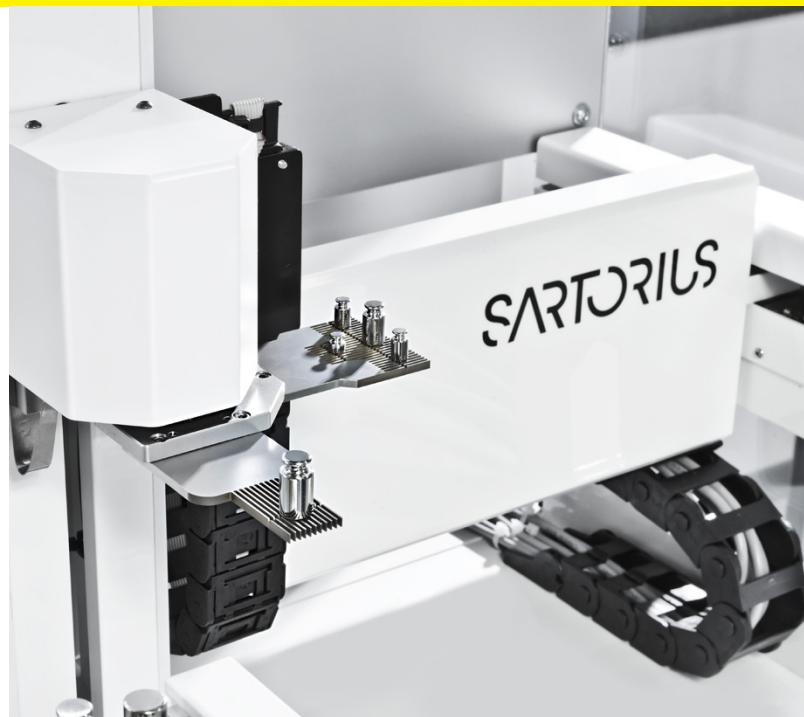
120 magazine positions enable loading of large number of weights and allow continuous processing of the measurements overnight or over the weekend. The magazine was designed to directly hold any possible weight shapes used for OIML-class and ASTM-class weights without the need of additional weight carriers or weight holders.





### Unique Robotic Arm Technology for Efficient Motion Sequences

The Sartorius dual robotic arm system with the patented multi and single weight handlers was designed for simultaneous management of reference and test weights. The two single weight handlers allow simplified and fast handling of the reference and test weight directly between the magazine and the weighing cell. The dual arm system with a multi- and single weight handler is able to collect up to 4 weights and the reference weight. The 4 weights are assembled at the collect station and measured afterwards simultaneously. Therefore, there is no need of returning the weights to the magazine or in a parking position.



### Integrated Climate Sensor

The weighing cells of all three models have a built-in climate sensor which means that the parameters for temperature, humidity and atmospheric pressure are measured at the exact location of the mass determination, therefore guaranteeing the most accurate correction for air buoyancy.



### Error Free Weight Sorting

The weight sorting plate designed specially for these systems allows an easy and error free processing of the large number of weights. Each row and column is labelled and is identically to the labeling on the weight magazine. Additionally, each row is removable which further support an intuitive and efficient weight handling.



### Variability to Adjusts to Your Workflows

Three different variants not only cover mass dissemination requirements from National Metrology Institutes and Legal Metrology Labs accredited to Class E1, but offer attractive high-throughput solutions for calibration laboratories typically accredited to E2 or F1.









# Technical Specifications

Models	CCR10.7-C	CCR10.6-C	CCR6.7C
Capacity	10 g	10 g	6 g
Readability	0.1 µg	1.0 µg	0.1 µg
Repeatability Typical	0.2 µg	0.5 µg	0.2 µg
Magazine Positions	120	120	120
Robotic Arm Technology	Dual Arm System	Dual Arm System	Dual Arm System
Weight Handlers	1 Multi & 1 Single	2 × Single	1 Multi & 1 Single
Dimensions (W × D × H mm)	1200 × 800 × 760	1200 × 800 × 760	1200 × 800 × 760

## Key Accessories

Order Code		Description
YWT12		Weighing Table made of granite, weight: 680 kg, dimensions (W × D × H mm) 800 × 1200 × 800
YWT13		Weighing Table made of granite and steel, weight: 330 kg, dimensions (W × D × H mm) 800 × 1200 × 800
YCM20MC-Tower		Climate tower monitors the influencing variables of the environment (temperature, pressure, and humidity)
YCM20DAKKS		DakKS calibration certificate for climate sensors
YAW10CCR-C		Weight sorting plate with removable trays for easy and error free processing of the large number of weights. Each row and column is labelled and is identical to the labelling on the weight magazine.
YSN03C   YSN03NC   YSN03RC		PC Software ScalesNet-M
YCW322-02		2 g Class E2 with DAkKS calibration certificate for model CCR10.7-C
YCW412-02		10 g Class E2 with DAkKS calibration certificate for model CCR10.6-C
YCW352-0		5 g Class E2 with DAkKS calibration certificate for model CCR6.7-C



Find out more  
Brochure: CCR-Compact Robotic Systems



Find out more  
White Paper: Accurate and High Throughput  
Mass Determination with Table-Top Robotic Systems




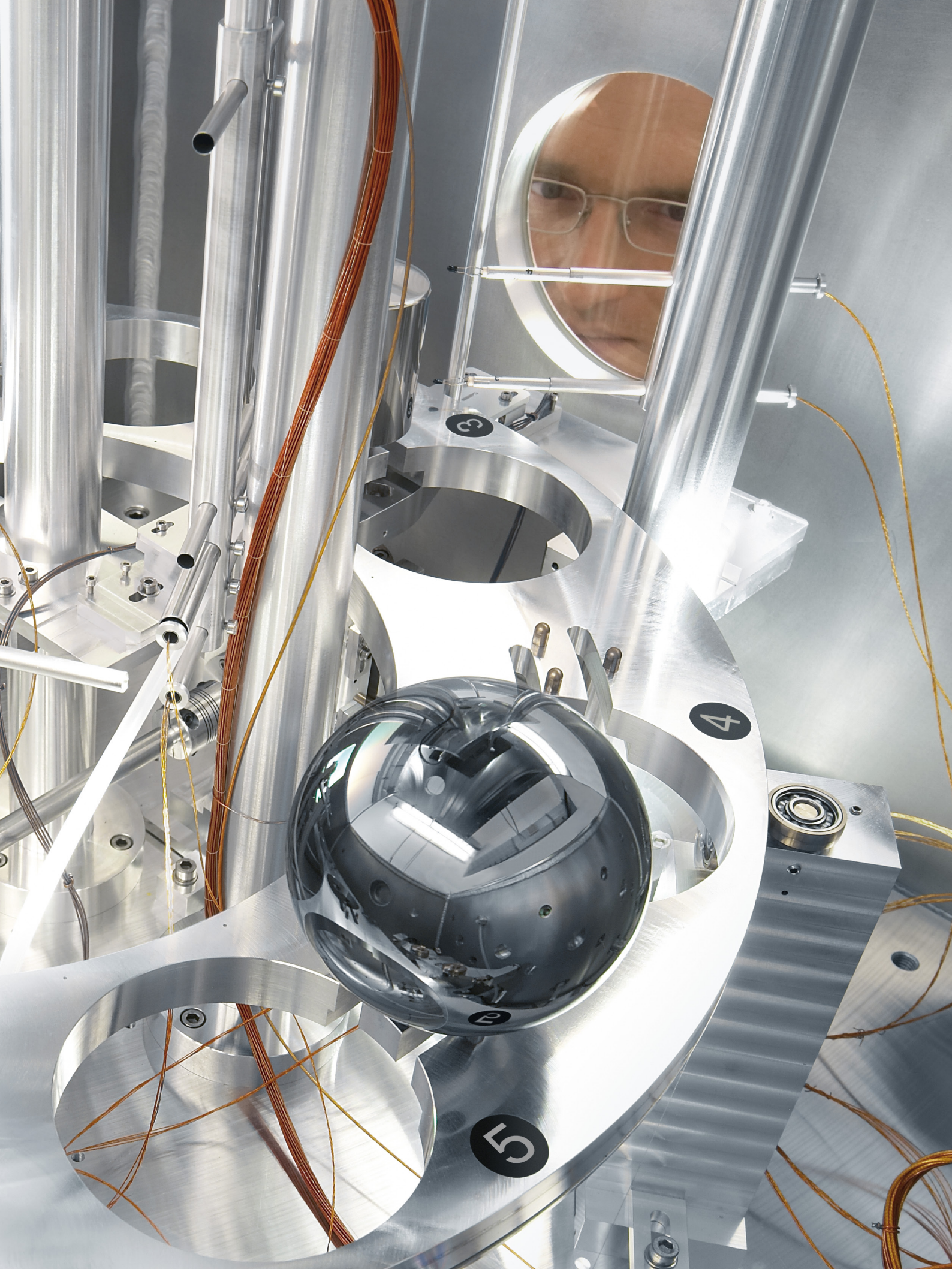
Find out more  
Datasheet: Table-Top Robotic Mass Comparators

# Automatic Mass Comparators and Robots

- 28 Mass Comparator CCL1007 for National Prototype Kilogram
- 32 Robot for Fully Automated Determination of Mass from 1 mg to 1 kg



 Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)



# Mass Comparator CCL1007 for National Prototype Kilogram

## For Mass Comparisons of Silicon Spheres and Weight Artefacts

Sartorius has mastered the core disciplines of weighing like no other company. It is also setting benchmarks in mass metrology. In cooperation with the Bureau International des Poids et Mesures and the Institute for Process Measurement and Sensor Technology of the Technical University of Ilmenau, Sartorius has developed a mass comparator – the CCL1007 – that is capable of determining differences in mass to an accuracy of 0.1 µg for weights of 1 kg – even under high-vacuum conditions.

The CCL1007 comparator is optimally protected from environmental effects by a pressure-stable vacuum chamber made of aluminum with a housing of polymethyl methacrylate (PMMA). Measurements can be performed under normal pressure conditions as well as at high vacuum up to a pressure of  $10^{-6}$  hPa. Unlike conventional stainless steel vacuum bell jars, the aluminum chamber does not cause any magnetic reactions that would affect the weighing system and weights.

The system is equipped with a fully automatic load alternator which can take up to eight weights simultaneously. Since the weights are stored on a unique threepoint frame, it is possible to use cylindrical as well as spherical weight. Thus, the device fulfills the essential prerequisites for highly accurate metrological tests for the new definition of the kilogram based on the Avogadro constant using silicon spheres.

Thanks to its automatic loading device, the weight sets can be positioned on the load alternator without having to open the entire vacuum chamber.

During loading, the center of mass of the weight is determined. A special procedure automatically corrects any de-centering of the weight which is then placed in the center of the load alternator. For contamination-free loading, the mass comparator CCL1007 can additionally be equipped with a vacuum transfer channel (YVTS01C).

The actual weigh cell is located in the upper section of the vacuum chamber. In case the reference standards on the load alternator deviate from one kilogram, the weighing pan for the substitution weights can be accessed via the upper loading hatch. Moreover, it is also possible to remove the internal one-gram calibration weight and perform calibrations separately.

All components inside the vacuum chamber are completely resistant to high vacuum. This prevents any contamination of the weights. All materials used in the vacuum chamber are solid materials with a low vapor pressure. They are free of oil and grease.

The driving forces necessary for rotating the load alternator and lifting and lowering the weights are transferred into the interior of the vacuum chamber by rotary axes and vacuum couplings. The load alternator is driven by a pulley hoist system and positioned exactly. The more than 20 vacuum flanges enable additional sensors to be introduced into the vacuum chamber.

The motors that drive the load alternator and the related control device are located outside the vacuum chamber in a control unit. Ceramic axles and spatial separation prevent any energy input from these components onto the vacuum chamber. This minimizes temporal and spatial temperature gradients in the interior of the vacuum chamber.

The system is controlled via a user-friendly operating software, which can be extended with ScalesNet-M. ScalesNet-M (YSN03C, see page 44) a management software program designed to cover all requirements and to network equipment in mass standards laboratories.

The CCL1007 places special environmental and mechanical demands on rooms, installations and operations. Specially trained personnel are also required. For that reason, our specialists are there to assist you in equipping your laboratory through setup to startup and help you troubleshoot any problems.



CCL1007 with vacuum transfer system VTS

## Areas of Application

- As a 1 kg prototype weighing system for the subdivision of the mass of the primary reference standards of National Metrology Institutes (NMIs) with national mass reference standards (kilogram prototype)
- Dissemination of the mass scale of NMIs in the range of 1 kg to 1 mg; adapter plates are required for groups of weights and weights  $\leq 200$  g
- Determination of mass for international comparison measurements (key comparisons) and calibrations for national institutes, calibration laboratories and industry in accordance with the attainable measurement uncertainties specified in the CMC tables published by the BIPM
- Experimental determination of air density by comparison weighing of special buoyancy artifacts in air and under vacuum
- Determination of the mass of 1 kg silicon spheres (also suitable within the scope of the Avogadro project for more accurate definition of the Avogadro constants and for new definition of the kilogram mass unit)
- Experimental research on the impact of cleaning procedures as well as on the effects that sorption and convection could have on mass and on the longterm stability of mass standards



# Vacuum Transfer System VTS

The vacuum transfer system (VTS) not only allows air-to-air loading on the standard model, but air-to-vacuum and air-to-protective gas. That way the measurement conditions inside the vacuum chamber remain constant while loading the load alternator with weights. Moreover, the vacuum transfer channel can be connected with special transport containers.

These containers enable weights to be stored in the lab under protective gas or a vacuum. This makes it possible to open containers inside the vacuum transfer channel under protective gas or in a vacuum and to load the load alternator automatically. That means that the weights are only ever exposed to the ambient conditions desired. The transport containers thus also enable loading of vacuum to vacuum and of protective gas to protective gas.

## Technical Specifications

	<b>CCL1007</b>
Maximum capacity	1,031 g
Readability	0.1 µg
Repeatability s*	s ≤ 0.2 µg
Repeatability in vacuum, s*	s ≤ 0.1 µg
Repeatability (typical), s*	s ≤ 0.1 µg
Linearity	≤ 1 µg
Electronic weighing range	2 g
Pressure range	10 <sup>-6</sup> – 1,000 mbar
<b>Application Ranges</b>	
OIML R111 Classes	≤ E1 1 kg
with adapter plates for groups of weights	≤ E1 1 mg ...1 kg
Silicon sphere	Ø 95 mm

s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift. Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.

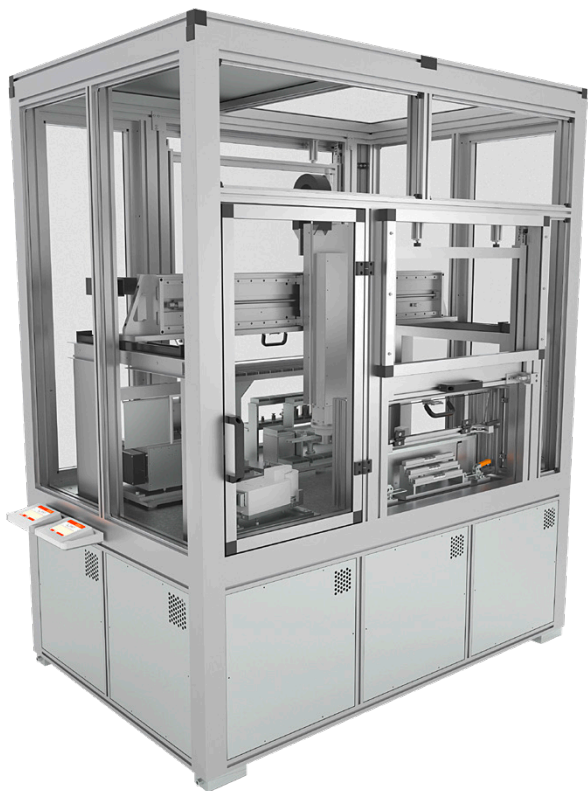
# Robot for Fully Automated Determination of Mass

From 1 mg to 1 kg

Robot systems with the latest weighing technology guarantee highest accuracy for the determination of mass. The fully automated mass comparison simplifies work in the mass laboratory. This enables complete sets of weights to be measured efficiently within the shortest time possible.

The robot systems are supplied with a user-friendly control software. An additional PC workplace is thus not needed. The CCR robot systems are equipped as standard with YCM20MC climate modules, which means that all climate parameters relevant for a mass comparison are available.

Robot systems, CCR10-1000, for determining the mass of weights (1 mg – 1 kg)



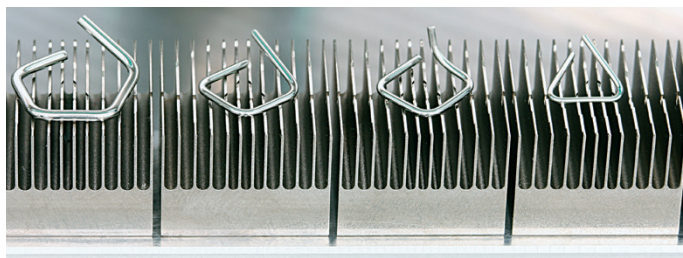
Comb-type weight grabber with 4 weights



Magazine of the CCR10-1000 (10 g – 1 kg)



Detail of the magazine on CCR10-1000 (1 mg – 10 g)





Thanks to the modular design of the robot system, the device can be adapted to customer requirements accordingly. The range of functions can also be extended at any time: The application range covers weights from 1 mg up to 10 g (CCR10) or 10 g up to 1 kg (CCR1000) and from 1 mg up to 1 kg in the maximum level of expansion (CCR10-1000). The number of magazine positions is specified by the customer: Up to 164 magazine positions can be provided for the range of 1 mg to 1 kg. Because you can already position your weights on the rotating magazines for the next measurements while running your current measurement procedure, you no longer have to sit through protracted waiting periods to acclimatize your weights.

Once the user has positioned the weights on the magazine positions and entered the weighing mode via the operating software, all mass comparisons can be performed automatically (direct comparison of weights or comparison of weight combinations). The corresponding measurement records, listing weighing and climate values, are generated and available for evaluation in a broad variety of programs. Naturally, all Sartorius robot systems can also be controlled directly by ScalesNet M. ScalesNet-M (YSN03C, see page 44) a management software program designed to cover all requirements and to network equipment in mass standards laboratories.

To allow the robot system to work quickly, the test weight and reference weight are placed on separate comb-type weight handlers. These weight handlers can accommodate a group of up to four weights. So that all weights of a weight group are placed on the load pan at the same time.

The mass comparison for E1 weights is continuously possible from 1 kg to 1 mg in one device (CCR10-1000). There is no need for time-consuming transfer of weights from device to device.

The magazine positions and comb-type weight handlers are designed so that any desired types of weight – wire or leaf weights or knob, cylindrical or disk weights – can be handled efficiently.

### Typical users are:

- National metrological institutes for presenting a mass scale from 1 kg to 1 mg
- Manufacturers of weights with maximum throughput of direct mass comparisons
- Calibration laboratories for increasing the throughput through mass comparisons even during night hours and on the weekends

## Technical Specifications

	CCR10	CCR1000	CCR10-1000
	1 mg – 10 g	10 g – 1 kg	1 mg – 1 kg
Maximum capacity	10.5 g	1,016 g	10.5 g   1,016 g
Readability	0.1 µg	1 µg	0.1 µg   1 µg
Repeatability (typical), s*	< 0.2 µg	< 2 µg	< 0.2 µg   < 2 µg
Linearity	1 µg	20 µg	1 µg   20 µg
Electronic weighing range	3.5 g	26 g	3.5 g   26 g
Magazine positions	39	23	39   23
Optional magazine positions	26 – 65	13 – 36	13 – 101


s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift. Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.





# Metrological Equipment and Accessories

- 36 Climate Station & Data Logger
- 38 Susceptometer & Permeability Indicator –  
A Complete Solution for Testing the Magnetic  
Properties of Weights
- 40 Volume and Density Determination
- 42 ScalesNet-M – The Individual Software Solution  
for Professionally Calibrating Weights
- 46 Weights and Weight Sets

 **Find out more**  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# Climate Station & Data Logger

With the YCM16C, Sartorius has developed a highly precise, unique climate station for continually recording ambient temperature data.

## Suitable for All Metrological Laboratories

The climate station YCM16C is the perfect solution for determining air density and can be used to monitor the room climate in all metrological laboratories.

## Room Monitoring

As is required for calibration laboratories, the YCM16C provides room monitoring for all of the accuracy classes given in accordance with OIML R111-1 up to class E1. The room climate data is constantly monitored and shown on the display.

## Air Buoyancy Correction

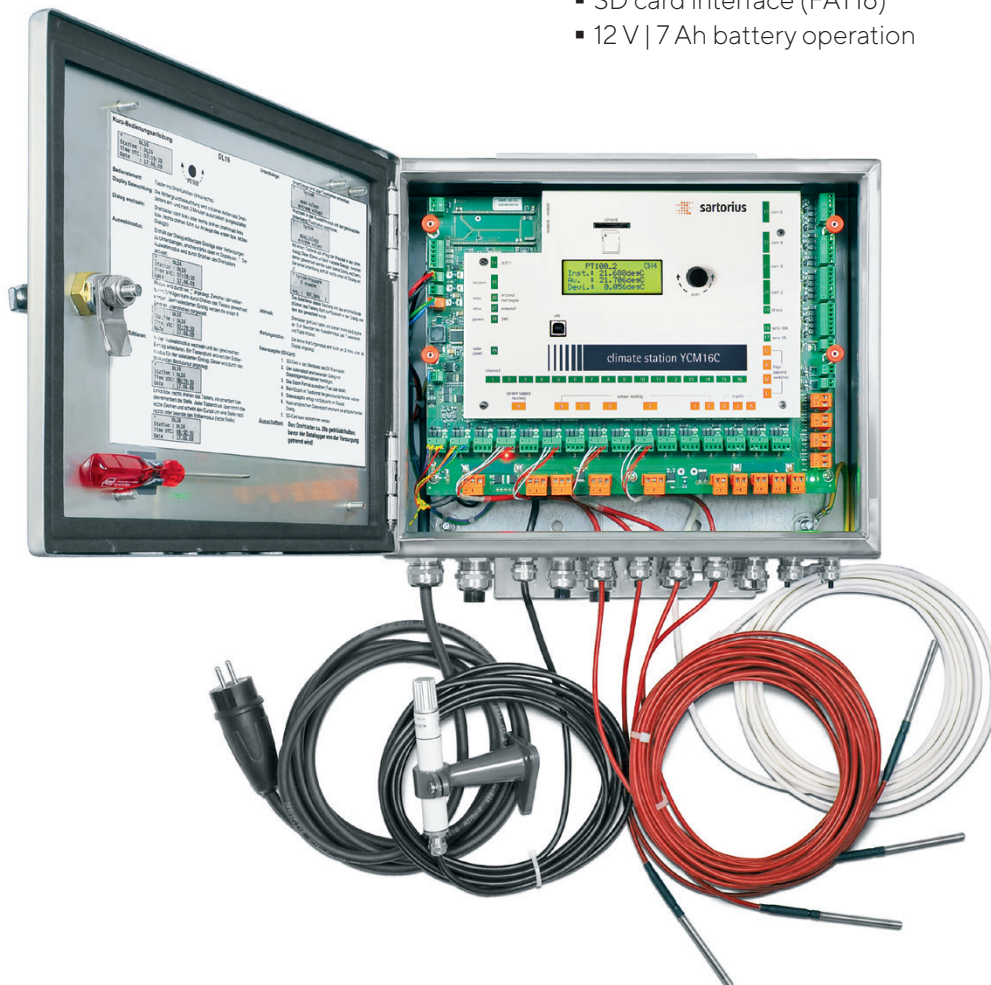
The climate station can be combined with ScalesNet-M software to make air buoyancy corrections in order to achieve an extremely high degree of accuracy when determining mass.

## Easy to Use

It can be operated directly using the touch key or from your desk using networkbased remote control. The supplied software automatically synchronizes the ambient climate data on your PC. If there is a power failure, the integrated UPS enables the climate station to keep recording reliable climate data even without a PC.

## Special Features

- Up to 16 analog and 99 digital sensors
- PC software for synchronization
- 8 MB data logger for up to 30 days of data recording
- Monitors the accuracy class in accordance with OIML R111
- Dew point calculation
- Air density calculation
- 1 integrated air pressure sensor
- 1 external humidity | temperature sensor
- WEB and FTP server
- UPS for power outages
- 4-line LCD display, 20 characters per line
- Can be operated directly using the push knob or remotely via USB | LAN | RS-232
- Firmware update via SD card
- LAN | USB | RS-485 | RS-232 interfaces
- SD card interface (FAT16)
- 12 V | 7 Ah battery operation



## Technical Specifications

	YCM16C
Temperature measuring range	-30 – +60 °C
Temperature readability	0.001°C
Humidity measuring range	0 – 100 %
Humidity readability	0.01 %
Pressure measuring range	300 – 1100 hPa
Pressure readability	0.001 hPa

## Optional Accessories

	Order number
Air temperature sensor 1/3 DIN	YCM16T
Air temperature sensor 1/10 DIN	YCM20T
Air humidity temperature sensor	YCM16H
Air pressure sensor	YCM16P
ScalesNet-M software	YSN03C



Calibrated air humidity | temperature sensor YCM16H | YCM16T



Calibrated air pressure sensor YCM16P



Air temperature sensor YCM20T

# Susceptometer & Permeability Indicator

## A Complete Solution for Testing the Magnetic Properties of Weights

### Susceptometer

It is recommended to test the magnetic properties of weights using the OIML R111-1. In fact, evidence supports the demand that both the susceptibility and magnetization of weights be accounted for.

### Susceptometer Method

There are several methods for defining the magnetic properties of weights. The susceptometer method is stipulated in OIML R111-1 for testing weights of classes E1, E2, F1, and F2. The Sartorius YSZ01C | YSZ02C susceptometers let you easily and conveniently determine the susceptibility and magnetization of weights in compliance with OIML R111-1.

This method measures the interaction between a permanent magnet and the weight to be tested using a high-resolution balance by weight force. The desired magnetic properties of the weight can be calculated on the basis of the known properties of the test magnet, the distance of the weight from the test magnet and the weight's geometry.

The Sartorius susceptometer stands out against other well-known instruments by virtue of its highly compact design and integrated components.

## Technical Specifications

Susceptometer	YSZ01C	YSZ02C
Readability	10 µg	1 µg
Application range in accordance with OIML R111	E2   F1   F2	E1   E2   F1   F2
Maximum capacity	50 kg	50 kg
$Z_0$ , distance from center of magnet to the bottom of weight in mm   field strengths $Z_0$ (A/m)	Adjustable in fixed steps: Z1 = 18   2,700 Z2 = 20   2,000 Z3 = 27   800	Z4 = 35   360 Z5 = 43   200
Turning mechanism for magnet	Easy to operate by external rotary knob, marking for N-S orientation of the magnet	
Software	Convenient application software; standard settings and easy, user-defined configuration possible; quick check function; printing of reports, exporting of results	
Data transfer protocol	HTML mode and data transfer to user-specific metrology software via CSV file format	

## Optional Accessories

Susceptometer	Order number
3 Reference magnets in a wooden case for Sartorius susceptometers YSZ01C and YSZ02C	YSZ01RMC
Susceptibility standard, as reference, in a wooden case for Sartorius susceptometers YSZ01C and YSZ02C, with PTB susceptibility certificate	YSZ01RSC



Susceptometer YSZ01C | YSZ02C

An innovative turning mechanism for adjusting the position of the test magnet and simple adjustment of the height of the loading platform also make working with the susceptometer simple and effective. This design significantly improves the repeatability of measurements. The susceptometer is available in two resolution levels. The YSZ01C features a 10- $\mu\text{g}$  resolution; the YSZ02C has a 1- $\mu\text{g}$  resolution.

### Application Software

The convenient software that comes standard with the susceptometer enables data to be automatically transferred from the unit. It moreover supports the user through all required handing steps and provides assistance with settings before and during a measurement. Menu prompts request all relevant data and guide the user through the steps necessary for performing a measurement. The results of the test are recorded and can be exported to higherlevel databases.

## Technical Specifications

Permeability Indicator	YAW61
Balance beams with a permanent magnet in the housing, including permeability reference standards [ $\mu\text{r}$ ]	1.01   1.02   1.03   1.05   1.06   1.08   1.15   1.20   1.30   1.60   1.80   1.90   5.0

### Permeability Indicator

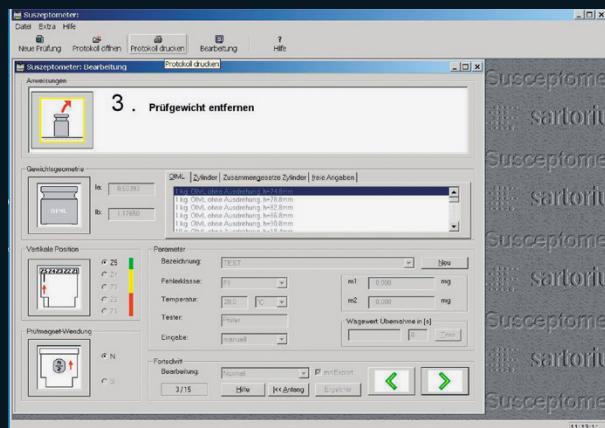
The Sartorius permeability indicator is easy to operate and is used to determine the relative permeability value of materials with small magnetic properties. The indicator operates according to the susceptometer method described in detail in OIML R111:2004 (Part 1, B.6.5). The magnetic susceptibility "x" is the result of the relationship between the relative magnetic permeability and the magnetic susceptibility ( $\mu\text{r} = 1 + x$ ).

x = magnetic susceptibility  
 $\mu\text{r}$  = relative magnetic permeability

The measuring principle of the indicator is based on the reciprocal attraction of a permanent magnet with a reference material of a known size along with an unknown material to be tested. This makes it possible to determine the relative permeability figure of the material to be tested by substituting the reference material as a limit value.



YAW61 with optional knob weight



Application software for susceptometers

### YSZ01RMC



# Volume & Density Determination

## Volume Comparators up to 1 kg

The most accurate method for determining density of solid bodies in accordance with OIML R111-1 is hydrostatic technique comparing the mass in a liquid.

Sartorius has integrated this method in its fully automatic volume comparator: With the VD1005, Sartorius supplies a system for density determination for weights in the range from 1 g up to 1 kg.

The volume comparator is equipped with two nine-position load alternators: One in a liquid bath and one in the air.

The two load alternators operate synchronously in that one position in liquid is assigned to each second position in air. By using substitution weights, this innovative dual-weight alternator design makes it possible to directly compare the mass of a single-volume reference (e.g., silicon sphere) with a variety of weights in liquid.

The load alternators and weighing pans are designed as comb-type weight grabbers so that weights and groups of weights from 1 g to 1 kg can be transferred directly from the load alternator to the suspended weighing pan. Adapter plates are not required. This has a positive effect on the accuracy of the overall measurement system.

After the data has been entered, an integrated PC with user-friendly software takes over fully automatic control of the volume comparator and evaluation of density measurement.



Volume comparator VD1005



## Application

Density determination of weights in accordance with OIML R111, class E1:  
1 g...1 kg



## Technical Specifications

	VD1005
Density uncertainty*	1 kg/m <sup>3</sup>
Volume uncertainty*	0.00015 cm <sup>3</sup>
Weight diameter	6...100 mm
Maximum sphere diameter	95 mm
<b>Comparator Technical Specifications</b>	
Maximum capacity	1,010 g
Readability	10 µg
Repeatability s*	< 40 µg
Repeatability (typical), s*	< 20 µg
Weighing range, electronic	1,010 g

\* Partial uncertainty of the volume comparator (without references and test weights)

s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift. Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.



Thermostat YVT01C

## Optional Accessories for Density Determination

	Order number
1 kg silicon sphere with PTB density certificate	YDR1000SIC
E1 set of 1 g to 1 kg weights with PTB density certificate (stackable) as a density reference and as a substitution weight set for volume comparators	YCS31-612-09



1 kg silicon sphere YDR1000SIC



E1 set of weights YCS31-612-09, stackable

# ScalesNet-M

## The Individual Software Solution for Professionally Calibrating Weights

With ScalesNet-M, Sartorius offers a customized solution for equipping a simple measurement laboratory through to the complete equipping of a national institute. Years of experience as a leading manufacturer of weights are incorporated in the development and have made ScalesNet-M into professional software for effective mass calibration.

### Individual

The modular concept of ScalesNet-M can be adapted to the highly diversified needs of our customers and supplemented at any time.

### Mass Calibration

With ScalesNet-M, mass calibration is traceable, secure and transparent at all levels. ScalesNet-M monitors and records all steps from the order arriving to the creation of calibration certificates, helping the customer reach their objective efficiently. ScalesNet-M is suitable for connection to all comparators from any manufacturer and flexible in incorporating existing climate measurement stations. At the same time, Sartorius provides a global service including initial installation, training, equipment connection and (remote) servicing and updating of ScalesNet-M.

### Automatic Monitoring

All means of testing, such as reference weights, comparators and climate sensors are automatically monitored by ScalesNet-M. Reminders are issued for all necessary calibrations and adjustments, which are then executed and documented by ScalesNet-M.

### Advantages

- Safe weight calibration
- Efficient measurement laboratory management
- Calculation and evaluation according to international guidelines such as OIML | ASTM
- Basis for an accredited measurement laboratory according to DIN EN ISO | IEC 17025
- Transparent documentation and archiving of all process-relevant data
- Automatic monitoring of test medium in laboratory
- Display of the customer and reference weight calibration history
- Incorporation of existing customer and calibration data into the ScalesNet-M database
- Individual software module compiling
- Connecting of comparators and climate measurement stations from any manufacturer
- Flawless, time-saving and automatic creation of multi-lingual, linguistically accurate calibration certificates
- Verifiable software
- Customizable service package

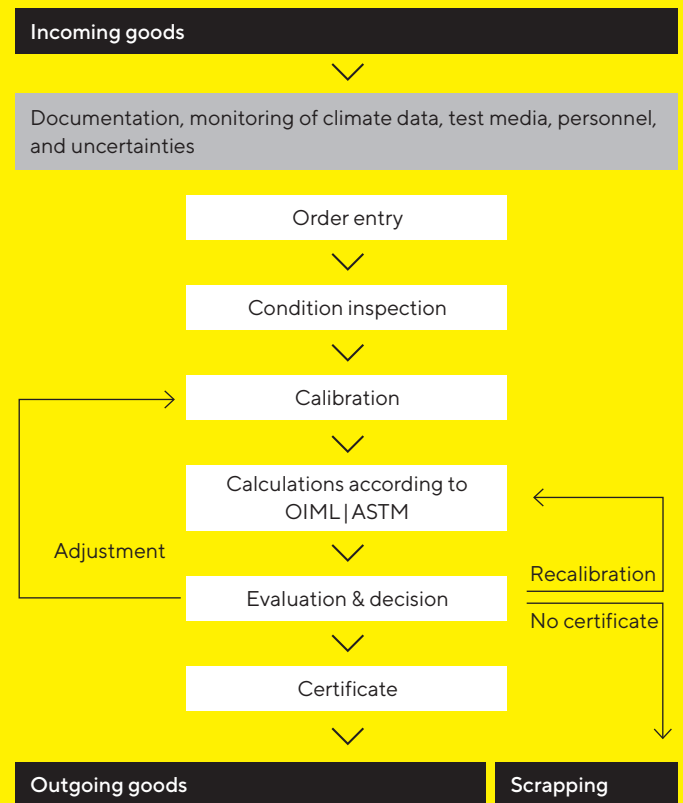


## Features

- Calibration of any class and value of weights
- Suitable for all weighing cycles in accordance with international guidelines (OIML | ASTM)
- Subdivision mass comparison for calibrating E1 weights
- Automatic evaluation of calibration results
- Rapid comparison of weights without logging
- Export function of weight values as a CSV file or JSON file
- Monitoring, recording and visual presentation of climate data
- Order preparation
- Examples integrated for weighing schemes
- Presentable history of all processes
- Cycles with and without additional weights and sensitivity weights
- All weight classes are already integrated in accordance with OIML and ASTM
- Inputting of own weight classes
- Automatic uncertainty calculation
- Continual test medium monitoring
- Automatic syntactically correct printout of calibration certificates
- DAkkS certificates and test log templates. Freely designable customer certificates in 2 languages
- SQL database structure for customer, weight and calibration data
- Administration of user rights
- Automatic generation of inventory lists
- Plausibility test when standard set and comparator are selected
- Manual input of weighing data for comparators without RS-232 connection



## Monitoring and Controlling the Calibration Process Using ScalesNet-M



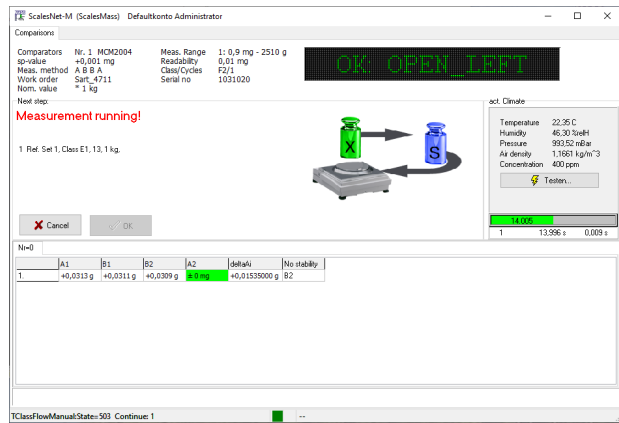
## Equipment Supplied

The YSN03C basic package contains one CD and one dongle with the following licenses:

- 1× ScalesServer, SQL database
- 1× ScalesMass, laboratory calibration
- 1× ScalesDesk, administration and user management
- 1× ScalesPrinter, printer control
- 1× ScalesPlan, data backup
- 1× License for 5× mass comparators,
- 1× License for OIML R111 F1, F2, M1, M2, M3
- ASTM 2, 3, 4, 5, 6, 7, NIST Handbooks F

## System Requirements

- PC or laptop with a resolution of at least 1024 × 768, with current Windows® version and internet access for remote maintenance
- Local administrator rights during installation for all PCs | Laptops Microsoft WORD
- PC connection cable for comparators



Software ScalesNet-M

## Reports

- DAKS calibration certificates
- Calibration report
- Inventory lists
- Device lists
- Six comparison log
- Dissemination log

## Languages

The software is available in German, English, French and Czech. Please contact us if another language is required.

## User Management

- User groups (read | write rights, administrator)
- Individual user accounts



# ScalesNet-M Module

## YSN03C – comprehensive module

Sartorius software module for highly accurate mass determination.

### Modules for supplementing the ScalesNet-M software



#### YSN03NC

##### Network module

PC license for connecting to local networks.



#### YSN03CC –

##### Module for a climate measuring station

PC license for connecting an additional climate measuring station.



#### YSN03LC

##### ScalesMass module

User license for simultaneous mass calibration on an additional PC. Only in conjunction with YSN03NC.



#### YSN03BC – Module for manual comparators

PC license for connecting up to five additional scales or manual mass comparators for simultaneous use.



#### YSN03PC

##### ScalesDesk module

Administrator license for the simultaneous use of an additional administrator on an additional PC. Only in conjunction with YSN03NC.



#### YSN03AC –

##### Module for automatic mass comparators

PC license for connecting automatic mass comparators with load alternator equipment.



#### YSN03DC

##### ScalesPrinter module

PC license for connecting an additional printer.



#### YSN03MC – Module for the mass comparison

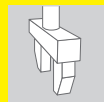
PC license for mass comparison for E1 weight classes and for calibrating weights of all accuracy classes in accordance with OIML R111 and ASTM E617.



#### YSN03EC

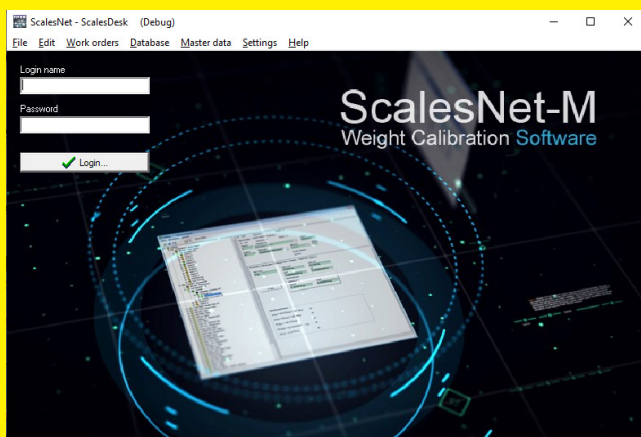
##### Class E module

PC license for calibrating weights in accordance with OIML R111 classes E and F as well as classes 0 to 4 in accordance with ASTM E617.

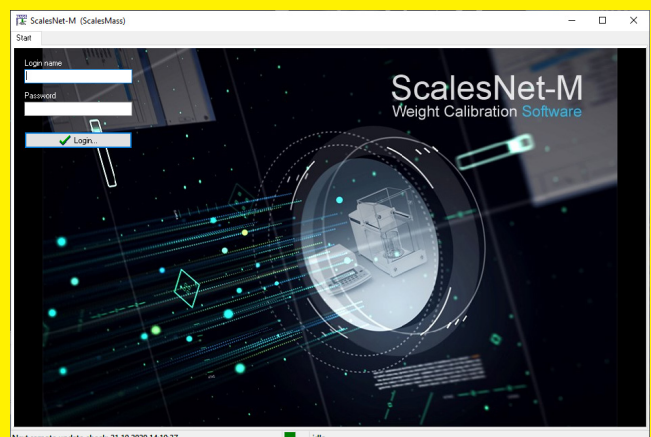


#### YSN03RC – Robot module

PC license for connecting a robot, a vacuum mass comparator (such as the CCL 1007) or a load alternator that uses the a\_control controlling software. Only in conjunction with YSN03NC.



ScalesNet-M, YSN03PC Administration Module



ScalesNet-M, YSN03LC Balance Communication Module

# Weights and Weight Sets

Regular inspection and testing of weighing instruments are a must to ensure reliable weighing results. Sartorius offers highly accurate metrological weights and weight sets with nominal mass values from 1 mg to 1,000 kg, special and test weights, as well as the accessories required for correct handling and storage of weights.

## **Unambiguousness | Traceability**

For unique classification, Sartorius weights are laser marked with a three-digit marking in accordance with the international OIML R111 recommendations. Sartorius weights are therefore uniquely identifiable.

Sartorius reference weights and sets of reference weights are the first choice for primary mass standards. They feature a first-class finish with a high gloss polished, corrosion stable surface. The quality of the reference weights exceeds the requirements of the international OIML R111 recommendation.

Sartorius test weights and sets of test weights are top-class working standards for everyday use. These weights are made of high-quality material with a polished or finely turned surface and also meet the international OIML R111 recommendation for shape, material, and markings.

They are therefore suitable for legal and general metrological applications in research and industry.



## DAkkS certificate

Sartorius reference, test, special, and custom weights are available with or without DAkkS certification.

Sartorius weights meet the requirements for traceability to the national kilogram prototype in conformance with ISO 9001:2000. They help support your quality management and quality assurance systems, and meet GLP and GMP requirements.

Details about our weights and other accessories can be found in the separate brochure or please visit our website at [www.sartorius.com](http://www.sartorius.com).

## Excerpt from Our Weight and Accessories Range

Weight sets	Order number DAkkS calibrated E1	Order number DAkkS calibrated E2
1 mg – 5 g	YCS011-351-02*	YCS011-352-02*
1 mg – 100 g	YCS011-511-02*	YCS011-512-02*
1 mg – 200 g	YCS011-521-02*	YCS011-522-02*
1 mg – 1 kg	YCS011-611-02*	YCS011-612-02*
1 mg – 5 kg	YCS011-651-02*	YCS011-652-02*
1 g – 1 kg	YCS31-611-02*	YCS31-612-02*
1 g – 5 kg	YCS31-651-02*	YCS31-652-02*
1 g – 10 kg	YCS31-711-02*	YCS31-712-02*

Glass bell jar with lower base	Order number
for 1 mg – 5 g	YAW00
for 1 mg – 200 g	YAW01
for 100 g – 1 kg	YAW02
for 2 kg – 5 kg	YAW03
for 10 kg	YAW04
for 20 kg	YAW05
for 50 kg	YAW06


\* Weights with DAkkS certificate in Sartorius's name for classes E2, F1, F2 up to 50 kg



Set of reference weights



Glass bell jar with lower base

 Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

## Accessories

Handles for lifting weights	Nominal value	Order number
Weight fork	for 500 g	YAW41
Weight fork	for 1 kg	YAW42
Weight fork	for 2 kg	YAW43
Handles for lifting weights	for 5 kg	YAW50
Handles for lifting weights	for 10 kg	YAW51
Handles for lifting weights	for 20 kg	YAW52
Handles for lifting weights	for 50 kg	YAW53
Crane with chain hoist		YLD01C
Gripper for weights with handle		YLD02C

Weighing tables	Order number
Weighing table made from synthetic stone (L x W x H) 900 x 600 x 760 mm	YWT03
Weighing table for Table-Top Robots, made of granite, weight 680 kg, dimensions (W x D x H) 800 x 1200 x 800 mm	YWT12
Weighing table for Table-Top Robots, made of granite and steel, weight 330 kg, dimensions (W x D x H) 800 x 1200 x 800 mm	YWT13

Weight Sorting Plate	Order number
Weight sorting plate designed for Table-Top Robots. The removable trays and the identical labelling to the 120 positions in the magazine ensure easy and error free processing of the large number of weights.	YAW10CCR-C

Special weighing pans	Order number
Weighing pan with set of weighing plates for weighing mass combinations for CC1000S-L	YWP04C
Centering pan for MCM40K3   MCM60K3   MCM60K2	YWP03C





Draft shield	Order number
for MCM67	YDS20C
for MCM1004   MCM2004   MCM5004   MCM5003   MCM36   MCM66   MCM106   MCM605   MCM1005	YDS24C
for MCM40K3   MCM60K3   MCM60K2   MCM10K3	YDS05C
for CCI60K2	YDS62C
for CCI100K2   CCI300K	YDS64C

Density references	Order number
1,000 g silicon sphere	YDR1000SIC
500 g silicon sphere	YDR500SIC
200 g silicon sphere	YDR200SIC
E1 set of 1 g to 1 kg weights with PTB density certificate (stackable) as a density reference and as a substitution weight set for volume comparators	YCS31-612-09

Printer	Order number
Thermo transfer and thermo direct printer for GxP printing on standard paper and self-adhesive labels	YDP30

Switch	Order number
Foot switch with T-connector	YFS01
Hand switch with T-connector	YHS02

Density determination	Order number
Below-balance weighing equipment for MCM40K3   MCM60K3   MCM60K2   CCE30002   CCE50002	69EA0040
Density kit for MCM36   MCM66   MCM106   MCM605   MCM1005	YDK01LP



Weighing table with optional accessories



Draft shield

# Detailed Technical Specifications

- 52 Cubis® MCM Mass Comparators up to 1 kg
- 54 Cubis® MCM Mass Comparators 2 kg to 10 kg
- 55 Cubis® MCM Mass Comparators 40 kg to 60 kg
- 56 Manual Mass Comparators 100 kg – 300 kg
- 57 Manual Mass Comparators 600 kg – 3,000 kg
- 58 Table-Top Robotic Mass Comparators
- 60 Floor-Standing Robotic Mass Comparators
- 61 Mass Comparator with Vacuum Chamber
- 62 Devices for Determination of Volume, Density, Susceptibility and Magnetism





# Cubis<sup>®</sup> MCM Manual Mass Comparators

Up to 1 kg



Order number with uncalibrated climate sensors	MCM6.7	MCM36	MCM66	MCM106
Order number with calibrated climate sensors with DAkkS certificate	MCM6.7-DAkkS	MCM36-DAkkS	MCM66-DAkkS	MCM106-DAkkS
Design	1	2	2	2
Maximum capacity	6.1 g	31 g	61 g	111 g
Readability	0.1 µg	1 µg	1 µg	1 µg
Application range	0 – 6 g	0 – 30 g	0 – 60 g	0 – 111 g
<b>Repeatability s</b>				
Under optimal conditions <sup>1)</sup>	0.15 µg	1 µg	1 µg	1 µg
Under standard conditions E <sup>2)</sup>	0.3 µg	1.5 µg	2 µg	2 µg
at 1/3 load <sup>2)</sup>	0.2 µg			
at 1/10 load <sup>2)</sup>		0.7 µg	0.7 µg	0.7 µg
Under standard conditions F <sup>3)</sup>	0.6 µg	4 µg	5 µg	5 µg
Electronic weighing   taring range	6.1 g	31 g	61 g	61 g
Substitution weights				50 g
Linearity	1 µg	6 µg	8 µg	8 µg
Eccentric load deviation	0.25 µg/mm	1 µg/mm	1 µg/mm	1 µg/mm
Stabilization time	10 s	3 s	3 s	5 s
<b>Standard Accessories</b>				
Data interfaces	RS-232C, USB, Ethernet, SD card (optional RS-232C, PS2, Bluetooth <sup>®</sup> )			
Draft shield	■	■	■	■
Additional application programs	Weighing, unit conversion, individual identifiers, density determination, statistics			
Below-balance weighing equipment	■	■	■	■
Climate sensors	Integrated into draft shield			
<b>Optional Accessories</b>				
Calibration weight	5 g   E2 YCW352-02	20 g   E2 YCW422-02	50 g   E2 YCW452-02	50 g   E2 YCW452-02
Climate module	YCM20MC			
Calibrated climate module	YCM20MC-DAkkS	YCM20MC	YCM20MC	YCM20MC
2nd Draft shield	YDS20C	YDS20MC-DAkkS	YDS20MC-DAkkS	YDS20MC-DAkkS
Weighing table	YWT03	YWT03	YWT03	YWT03
<b>Dimensions</b>				
Weighing pan size	Ø 16 mm	Ø 30 mm	Ø 30 mm	Ø 50 mm
Maximum object size (D × H)	16 × 70 mm	30 × 120 mm	30 × 120 mm	50 × 120 mm
Weigh cell (W × D × H)	122 × 343 × 141 mm	222 × 431 × 301 mm	222 × 431 × 301 mm	222 × 431 × 301 mm
Electronics unit (W × D × H)	239 × 320 × 56 mm	239 × 320 × 56 mm	239 × 320 × 56 mm	239 × 320 × 56 mm

Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift

<sup>1)</sup> Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>2)</sup> Standard conditions E: measured by hand in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>3)</sup> Standard conditions F: measured by hand in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above



Order number with uncalibrated climate sensors	MCM605	MCM1005	MCM1004
Order number with calibrated climate sensors with DAkkS certificate	MCM605-DAkkS	MCM1005-DAkkS	MCM1004-DAkkS
Design	3	3	3
Maximum capacity	610 g	1,110 g	1,110 g
Readability	0.01 mg	0.01 mg	0.1 mg
Application range	0 – 610 g	0 – 1,110 g	0 – 1,110 g
<b>Repeatability s</b>			
Under optimal conditions <sup>1)</sup>	10 µg	15 µg	0.05 mg
Under standard conditions E <sup>2)</sup>	20 µg	20 µg	0.07 mg
at 1/3 load <sup>2)</sup>	15 µg		
at 1/10 load <sup>2)</sup>	10 µg	15 µg	0.05 mg
Under standard conditions F <sup>3)</sup>	30 µg	50 µg	0.2 mg
Electronic weighing   taring range	610 g	610 g	610 g
Substitution weights		500 g	500 g
Linearity	100 µg	100 µg   600 g	0.1 mg   600 g
Corner load error	10 µg/mm	15 µg/mm	30 µg/mm
Stabilization time	5 s	5 s	3 s
<b>Standard Accessories</b>			
Data interfaces	RS-232C, USB, Ethernet, SD card (optional RS-232C, PS2, Bluetooth®)		
Draft shield	■	■	■
Additional application programs	Weighing, unit conversion, individual identifiers, density determination, statistics		
Below-balance weighing equipment	■	■	■
Climate sensors	Integrated into draft shield		
<b>Optional Accessories</b>			
Calibration weight	500 g   E2 YCW552-02	500 g   E2 YCW552-02	500 g   E2 YCW552-02
Climate module	YCM20MC		
Calibrated climate module	YCM20MC-DAkkS	YCM20MC	YCM20MC
2nd Draft shield	YDS24C	YDS24C-DAkkS	YDS24C-DAkkS
Weighing table	YWT03	YWT03	YWT03
<b>Dimensions</b>			
Weighing pan size	Ø 90 mm	Ø 90 mm	Ø 90 mm
Maximum object size (D × H)	135 × 140 mm	135 × 140 mm	135 × 140 mm
Weigh cell (W × D × H)	222 × 431 × 301 mm	222 × 431 × 301 mm	222 × 431 × 301 mm
Electronics unit (W × D × H)	239 × 320 × 56 mm	239 × 320 × 56 mm	239 × 320 × 56 mm

Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift

<sup>1)</sup> Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>2)</sup> Standard conditions E: measured by hand in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>3)</sup> Standard conditions F: measured by hand in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above

# Cubis<sup>®</sup> MCM Manual Mass Comparators

2 kg - 10 kg



Order number with uncalibrated climate sensors	MCM2004	MCM5004	MCM5003	MCM10K3
Order number with calibrated climate sensors with DAkkS certificate	MCM2004-DAkkS	MCM5004-DAkkS	MCM5003-DAkkS	MCM10K3-DAkkS
Design	4	4	4	5
Maximum capacity	2,500 g	5,100 g	5,100 g	11 kg
Readability	0.1 mg	0.1 mg	1 mg	1 mg
Application range	0 – 2,500 g	0 – 5,100 g	0 – 5,100 g	0 – 11 kg
<b>Repeatability s</b>				
Under optimal conditions <sup>1)</sup>	0.05 mg	0.3 mg	0.5 mg	0.8 mg
Under standard conditions E <sup>2)</sup>	0.1 mg	0.5 mg	0.8 mg	1 mg
at 1/3 load <sup>2)</sup>				
at 1/10 load <sup>2)</sup>	0.07 mg	0.3 mg	0.5 mg	0.8 mg
Under standard conditions F <sup>3)</sup>	0.3 mg	0.8 mg	1.5 mg	3 mg
Electronic weighing   taring range	2,500 g	5,100 g	5,100 g	11 kg
Linearity	1 mg	2 mg	3 mg	6 mg
Corner load error	30 µg/mm	151 µg/mm	300 µg/mm	0.5 mg/mm
Stabilization time	3 s	3 s	3 s	3 s
<b>Standard Accessories</b>				
Data interfaces	RS-232C, USB, Ethernet, SD card (optional RS-232C, PS2, Bluetooth <sup>®</sup> )			
Draft shield	■	■	■	
Additional application programs	Weighing, unit conversion, individual identifiers, density determination, statistics			
Below-balance weighing equipment	■	■	■	■
Climate sensors	Integrated into draft shield			Remote tower connected to mass comparator
<b>Optional Accessories</b>				
Calibration weight	2 kg   E2 YCW622-02	5 kg   E2 YCW652-02	5 kg   E2 YCW652-02	10 kg   E2 YCW712-02
Climate module	YCM20MC			
Calibrated climate module	YCM20MC-DAkkS	YCM20MC	YCM20MC	YCM20MC
2nd Draft shield	YDS24C	YDS24C-DAkkS	YDS24C-DAkkS	YDS24C-DAkkS
Weighing table	YWT03	YWT03	YWT03	YWT03
Lifting device for 10 kg				YAW51
<b>Dimensions</b>				
Weighing pan size (W×D)	136×136 mm	136×136 mm	136×136 mm	200×200 mm
Maximum object size (D×H)	130×200 mm	130×200 mm	130×200 mm	
Weigh cell (W×D×H)	240×276×373 mm	240×276×373 mm	240×276×373 mm	240×276×102 mm
Electronics unit (W×D×H)	239×320×56 mm	239×320×56 mm	239×320×56 mm	239×320×56 mm

Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift

<sup>1)</sup> Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>2)</sup> Standard conditions E: measured by hand in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>3)</sup> Standard conditions F: measured by hand in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above

# 40 kg - 60 kg



Order number with uncalibrated climate sensors	MCM32002	MCM40K3	MCM60K3	MCM60K2
Order number with calibrated climate sensors with DAkkS certificate	n/a	MCM40K3-DAkkS	MCM60K3-DAkkS	MCM60K2-DAkkS
Design	7	6	6	6
Maximum capacity	32 kg	41 kg	64 kg	64 kg
Readability	10 mg	1 mg	2 mg	10 mg
Application range	10 – 30 kg	0 – 41 kg	0 – 64 kg	0 – 64 kg
<b>Repeatability s</b>				
Under optimal conditions <sup>1)</sup>	10 mg	2 mg	4 mg	6 mg
Under standard conditions E <sup>2)</sup>	15 mg	3 mg	6 mg	10 mg
at 1/3 load <sup>2)</sup>	-			
at 1/10 load <sup>2)</sup>	-	2 mg	4 mg	
Under standard conditions F <sup>3)</sup>	25 mg	6 mg	10 mg	25 mg
Electronic weighing   taring range	32 kg	41 kg	64 kg	64 kg
Linearity	50 mg	20 mg	40 mg	50 mg
Corner load error	2 mg/mm	3.5 mg/mm	3.5 mg/mm	3.5 mg/mm
Stabilization time	5 s	5 s	5 s	5 s
<b>Standard Accessories</b>				
Data interfaces	RS-232C, USB, Ethernet, SD card (optional RS-232C, PS2, Bluetooth®)			
Additional application programs	Weighing, unit conversion, individual identifiers, density determination, statistics			
Below-balance weighing equipment	■	with opt. accessories 69EA0040	with opt. accessories 69EA0040	with opt. accessories 69EA0040
Climate sensors	no	Remote tower connected to mass comparator		
<b>Optional Accessories</b>				
Calibration weight	20 kg   E2 YCW722-02	20 kg   E2 YCW722-02	50 kg   E2 YCW752-02	50 kg   E2 YCW752-02
Climate module	n/a	YCM20MC		
Calibrated climate module	n/a	YCM20MC-DAkkS	YCM20MC	YCM20MC
Optional external draft shield	YDS05C   YDS03C	YDS05C   YDS03C	YDS05C   YDS03C	YDS05C   YDS03C
Lifting device for 10 kg	YAW51	YAW51	YAW51	YAW51
Lifting device for 20 kg	YAW52	YAW52	YAW52	YAW52
Lifting device for 50 kg			YAW53	YAW53
Crane with chain hoist			YLD01C	YLD01C
Gripper for weights with handle			YLD02C	YLD02C
Floor-mounted column, stainless steel				
<b>Dimensions</b>				
Weighing pan size (W × D)	Ø 230 mm	400 × 300 mm	400 × 300 mm	400 × 300 mm
Weigh cell (W × D × H)	400 × 326 × 126 mm	400 × 326 × 126 mm	400 × 326 × 126 mm	400 × 326 × 126 mm
Electronics unit (W × D × H)	239 × 320 × 56 mm	239 × 320 × 56 mm	239 × 320 × 56 mm	239 × 320 × 56 mm

Repeatability is the standard deviation “s”; it is calculated from 5 ABA cycles, after eliminating drift

<sup>1)</sup> Optimal conditions: automatic measurement without operator influence measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>2)</sup> Standard conditions E: measured by hand in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>3)</sup> Standard conditions F: measured by hand in a laboratory under at least F1 conditions, on a non-decoupled weighing stone, air conditioning and minimal drafts from above

# Manual High-Capacity Mass Comparators

100 kg - 300 kg



	CCI100K2	CCI300K
Maximum capacity	151 kg	303 kg
Readability	0.05 g	1 g
Application range	0 - 151 kg	0 - 303 kg
Repeatability s*	0.3 g	1 g
Repeatability at 1/10 load		
Repeatability (typical), s*	0.2 g	0.5 g
Electronic weighing   taring range	151 kg	303 kg
Linearity	2 g	10 g
Corner load error	25 mg/mm	50 mg/mm
Stabilization time	10 s	10 s
<b>Standard Accessories</b>		
Interfaces	2× RS-232C	2× RS-232C
Draft shield	■	
Application programs		
Below-balance weighing equipment		
PC connection cable	■	■
<b>Optional Accessories</b>		
Calibration weight	50 kg   F1 YCW753-02-DAkkS	100 kg   F1 YCW813-02
PC software ScalesNet-M	YSN03C	YSN03C
Climate station for E1	YCM16C	YCM16C
2nd Draft shield		YDS64C
Lifting device for 10 kg	YAW51	YAW51
Lifting device for 20 kg	YAW52	YAW52
Lifting device for 50 kg	YAW53	YAW53
Crane with chain hoist	YLD01C	YLD01C
Gripper for weights with handle	YLD02C	YLD02C
Floor-mounted column, stainless steel	YDH03CIS	YDH03CIS
<b>Dimensions</b>		
Weighing pan dimensions (W×D)	800×600×135 mm	800×600×135 mm
Weigh cell (W×D×H)		
Electronics unit (W×D×H)	303×195×90 mm	303×195×90 mm

s\* Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift.

Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.



## 600 kg - 3,000 kg



	CCS600K	CCT1000K	CCS1000K	CCT2000K	CCS3000K
Maximum capacity	605 kg	1,200 kg	1,510 kg	2,100 kg	3,010 kg
Readability	1 g	1 g	5 g	1 g	10 g
Application range	0 - 605 kg	0 - 1,200 kg	0 - 1,510 kg	0 - 2,100 kg	0 - 3,010 kg
Repeatability s*	2.5 g	3 g	6 g	8 g	12 g
Repeatability (typical), s*	2 g	2 g	5 g	5 g	10 g
Electronic weighing   taring range	605 kg	1,200 kg	1,510 kg	2,100 kg	3,010 kg
Linearity	30 g	25 g	200 g	30 g	500 g
Corner load error	0.2 g/mm	0.3 g/mm	0.6 g/mm	0.7 g/mm	1 g/mm
Stabilization time	20 s	30 s	20 s	30 s	20 s
<b>Standard Accessories</b>					
Interfaces	2× RS-232C	2× RS-232C	2× RS-232C	2× RS-232C	2× RS-232C
Draft shield	■	■	■	■	■
PC connection cable	■	■	■	■	■
<b>Optional Accessories</b>					
Calibration weight	500 kg   F1 YCW853-02	1000 kg   F2 YCW914-02	1000 kg   F2 YCW914-02	1000 kg   F2 YCW914-02	1000 kg   F2 YCW914-02
PC software ScalesNet-M	YSN03C	YSN03C	YSN03C	YSN03C	YSN03C
Climate station for E1	YCM16C	YCM16C	YCM16C	YCM16C	YCM16C
Floor-mounted column, stainless steel	YDH03CIS	YDH03CIS	YDH03CIS	YDH03CIS	YDH03CIS
Base for installing floor-mounted column, stainless steel	YBP03CIS	YBP03CIS	YBP03CIS	YBP03CIS	YBP03CIS
<b>Dimensions</b>					
Weighing pan dimensions (W×D)	830×1,030× 250 mm	1,510×1,370× 240 mm	830×1,030× 250 mm	1,920×1,710× 230 mm	1,000×1,250× 300 mm
Electronics unit (W×D×H)	303×195×90 mm	303×195×90 mm	303×195×90 mm	303×195×90 mm	303×195×90 mm
Net weight	150 kg	225 kg	150 kg	400 kg	300 kg
Gross weight	250 kg	338 kg	250 kg	536 kg	470 kg
Pallet	1.5×1.1×0.6 m	1.8×1.8×0.6 m	1.5×1.1×0.6 m	2.1×2.2×0.6 m	1.7×1.6×0.6 m

s\* Repeatability is the standard deviation "s"; it is calculated from 5 ABA cycles, after eliminating drift.

Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an M1 mass standards laboratory.

# Table-Top Robotic Mass Comparators

Up to 10 g



## Technical Specifications

Model	CCR10.7-C	CCR10.6-C	CCR6.7-C
Maximum capacity	10.5 g	10.1 g	6.1 g
Application range	1 mg - 10 g	1 mg - 10 g	1 mg - 6 g
Readability	0.1 µg	1 µg	0.1 µg
Repeatability typical	0.2 µg	0.5 µg	0.2 µg
Rep. under standard conditions E <sup>1</sup>	0.5 µg	0.7 µg	0.3 µg
Repeatability, at > 1 - 6 g	0.3 µg		
Repeatability, at 0 - 1 g	0.15 µg		0.15 µg
Rep. under standard conditions F <sup>2</sup>	1.5 µg	2 µg	1.5 µg
Electronic weighing range & taring range	3.5 g	10.1 g	6.1 g
Substitution weights	2 × 3.5 g	-	-
Linearity	1 µg   3.5 g	4 µg	1 µg
Off-center loading error	0.25 µg / mm	0.5 µg / mm	0.25 µg / mm
Stabilisation time	15 s	10 s	15 s

<sup>1</sup> Standard conditions E: ABA measured in a laboratory under E1 conditions, on a decoupled weighing stone, no drafts from above

<sup>2</sup> Standard conditions F: ABA measured in a laboratory under at least F1 conditions, on a non-decoupled weighing table, air conditioned and minimal drafts from above

**Important:** The stated specifications and technical data apply only under good ambient conditions. Disruptive factors at the place of installation, such as strong drafts (especially from air conditioning equipment), excessive vibrations, physical effects of the items being weighed (e.g. magnetic fields or electrostatic charges), or ambient conditions outside the allowable tolerances, may adversely affect on the specifications.

## Models

Model	CCR10.7-C	CCR10.6-C	CCR6.7-C
Type	4-axis robot	4-axis robot	4-axis robot
Robotic Arm Technology	Dual Arm System	Dual Arm System	Dual Arm System
Weight Handlers	1 Multi × 1 Single	2 × Single	1 Multi × 1 Single
Magazine positions	120	120	120
Interfaces	LAN	LAN	LAN
Draft shield	■	■	■
Enclosure	■	■	■
Laptop PC	-	-	-
PC software	■	■	■
Climate sensor integrated in the weighing chamber to measure humidity, air pressure & air temperature	■	■	■
Test certificate	Sartorius	Sartorius	Sartorius

## Ambient Conditions

Model	CCR10.7-C	CCR10.6-C	CCR6.7-C
Permissible operating temperature range	17–27°C	17–27°C	17–27°C
Recommended operating temperature	22°C	22°C	22°C
Temperature fluctuation	0.3°C/h 0.5°C/12h	0.3°C/h 0.5°C/12h	0.3°C/h 0.5°C/12h
Max. air movement	< 0.2m/s	< 0.2m/s	< 0.2m/s
Humidity range	40–60%	40–60%	40–60%
Humidity fluctuation	5%/4h	5%/4h	5%/4h
Power supply	100–240V AC/50–60Hz	100–240V AC/50–60Hz	100–240V AC/50–60Hz

## Dimensions

Model	CCR10.7-C	CCR10.6-C	CCR6.7-C
Weighing pan dimensions (W×D)	49×29 mm	49×29 mm	49×29 mm
Sample size (D×H)	18×20 mm	18×20 mm	18×20 mm
External dimensions W×D×H	1,200×800×760 mm	1,200×800×760 mm	1,200×800×760 mm
Gross weight	265 kg	265 kg	265 kg
Net weight	190 kg	190 kg	190 kg
Number of packages	1	1	1
Pallet	1400×980×1400 mm	1400×980×1400 mm	1400×980×1400 mm
Optimal height for setup	800 mm	800 mm	800 mm

## Optional Accessories

Model	CCR10.7-C	CCR10.6-C	CCR6.7-C
External Calibration weight	2g   E2 YCW322-02	10g   E2 YCW412-02	5g   E2 YCW352-02
PC Software ScalesNet-M	YSN03C + YSN03RC (+ YSN03MC optional)	YSN03C + YSN03RC	YSN03C + YSN03RC (+ YSN03MC optional)
Weight sorting plate	YAW10CCR-C	YAW10CCR-C	YAW10CCR-C
External climate Sensor	YCM20MC-Tower	YCM20MC-Tower	YCM20MC-Tower
DakkS calibration certificate for climate sensors	YCM20DAKKS	YCM20DAKKS	YCM20DAKKS
Weighing Table premium	YWT12	YWT12	YWT12
Weighing Table budget	YWT13	YWT13	YWT13

# Floor-Standing Robotic Mass Comparators

Up to 1 kg



	CCR10	CCR1000	CCR10-1000
Maximum capacity	10.5 g	1,016 g	1,021 g
Readability	0.1 µg	1 µg	0.1 µg   1 µg
Application range	1 mg – 10 g	10 g – 1 kg	1 mg – 1 kg
Repeatability s*	0.5 µg	8 µg	0.5 µg   8 µg
Repeatability at 1/2 load	0.3 µg	5 µg	0.3 µg   5 µg
Repeatability at 1/10 load	0.2 µg	3 µg	0.2 µg   3 µg
Repeatability (typical), s*	0.2 µg	2 µg	0.2 µg   2 µg
Electronic weighing   taring range	3.5 g	21 g	3.5 g   21g
Substitution weights	2×3.5 g	30   40   50, 2×100   300   400 g	2×3.5   30   40   50, 2×100   300   400 g
Linearity	1 µg   3.5 g	20 µg   21 g	1 µg   3.5 g   20 µg   21 g
Corner load error	0.25 µg / mm	0 mg	0.25 µg / mm   0 mg
Stabilization time	15 s	25 s	15 s   25 s
<b>Standard Accessories</b>			
Load alternator positions	2	2	2   2
Magazine positions	39 pos.	23 pos.	39 pos.   23 pos.
Interfaces	LAN   USB   RS-232	LAN   USB   RS-232	LAN   USB   RS-232
Centermatic	-	•	-   •
Enclosure	■	■	■
PC	■	■	■
PC software	■	■	■
Climate module	YCM20MC	YCM20MC	YCM20MC
<b>Optional Accessories</b>			
PC software ScalesNet-M	YSN03C	YSN03C	YSN03C
Calibration weight	2 g   E2 YCW322-02-DAkks	20 g   E2 YCW422-02-DAkks	2 g   E2 YCW322-02-DAkks and 20 g   E2 YCW422-02-DAkks
<b>Extensions</b>			
Upgrade kit	Y1000UPGRADE	Y10UPGRADE	
Reference magazine	Y10R:(26)	Y1000R:(13)	Y10R:(26)   Y1000R:(13)
2nd turning magazine	Y10M:(39)	Y1000M:(23)	Y10M:(39)   Y1000M:(23)
<b>Dimensions</b>			
Weighing pan dimensions (W×D)	49×29 mm	104×68 mm	49×29 mm   104×68 mm
Sample size (D×H)	18×20 mm	100×120 mm	18×20 mm   100×120 mm
Exterior measurements (W×D×H)	1,900×1,250×2,328 mm	1,900×1,250×2,328 mm	1,900×1,250×2,328 mm
Net weight	1,900 kg	1,900 kg	1,900 kg

s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift.

Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.

# Mass Comparators

With Vacuum Chamber



	CCL1007	YVP01C	YVTS01C	YVC01C
Maximum capacity	1031 g		1 kg	1 kg
Readability	0.1 µg			
Application range	1 mg – 1 kg			
Repeatability s*	0.2 µg			
Repeatability in vacuum, s*	0.1 µg			
Repeatability (typical), s*	0.1 µg			
Electronic weighing   taring range	2 g			
Linearity	1 µg			
Range sensitivity	0.2 µg – 500 mg			
Stabilization time	80 s			
Pressure range	10E-6 – 1000 mbar	10E-6 – 1000 mbar	10E-6 – 1000 mbar	10E-6 – 1000 mbar
Helium leak rate	< 10E-7 (mbar*l)/s	< 10E-7 (mbar*l)/s	< 10E-7 (mbar*l)/s	< 10E-7 (mbar*l)/s
<b>Standard Accessories</b>				
Load alternator positions	8		1	
Interfaces	LAN   USB   RS-232			
isoCAL				
Centermatic	■			
Enclosure	■			
Control unit	■		■	
PC	■			
PC software	■			
<b>Optional Accessories</b>				
Turbopump Stand	YVP01C		YVP02C	YVP02C
Vacuum Transfer System	YVTS01C			
Vacuum container	YVC01C		YVC01C	
Climate station for E1	YCM16C			
PC software ScalesNet-M	YSN03C			
Calibration weight	2 g   E2 YCW322-02			
<b>Dimensions</b>				
Sample size (D × H)	34 – 95 × 110 mm		34 – 95 × 110 mm	34 – 95 × 110 mm
Diameter range for a silicon sphere	45 – 100 mm		45 – 100 mm	45 – 100 mm
Vacuum chamber (W × D × H)	960 × 860 × 1,260 mm		1,081 × 1,044 × 1,261 mm	
Control unit (W × D × H)	600 × 800 × 2.000 mm			
Net weight	350 kg	45 kg	200 kg	

s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift.

Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.

# Devices for Determination of Volume, Density, Susceptibility and Magnetism



	VD1005	YSZ02C	YSZ01C
Maximum capacity	1,110 g	50 kg	50 kg
Readability	10 µg	1 µg	10 µg
Application range	1 g – 1 kg	2 g – 50 kg	2 g – 50 kg
Repeatability s*	40 µg	8 µg	20 µg
Repeatability (typical), s*	20 µg	5 µg	10 µg
Electronic weighing   taring range	305 g	10 g	10 g
Linearity	0.12 mg	0.05 mg	0.1 mg
Stabilization time	20 s	15 s	15 s
Uncertainty of density	1 kg/m <sup>3</sup>		
Uncertainty of volume	0.00015 cm <sup>3</sup>		
<b>Standard Accessories</b>			
Load alternator positions	2×9		
Interfaces	LAN   USB   RS-232	RS-232	RS-232
Draft shield	■	■	■
Enclosure	■		
Control unit	■		
PC	■		
PC software	■	■	■
Air temperature sensor	1		
Humidity sensor	1		
Air pressure sensor	1		
Fluid temperature sensor	2		
Test certificate	BEV	Sartorius	Sartorius
PC connection cable		■	■
Climate station	■		
<b>Optional Accessories</b>			
Calibration weight	200 g   E2 YCW522-00	10 g   E2 YCW412-00	10 g   E2 YCW412-00
Weighing table	YWT20C	YWT03	YWT03
Crane with chain hoist		YLD01C	YLD01C
Gripper for weights with handle		YLD02C	YLD02C
1 kg PTB reference susceptibility		YSZ01RSC	YSZ01RSC
Magnet calibration kit for susceptometer		YSZ01RMC	YSZ01RMC
Density reference (Si)	200   500   1,000 g		
Density reference, set of weights 1 g – 1 kg	YCS31-612-09		


s\* Repeatability is the standard deviation "s"; it is calculated from 6 ABBA cycles, after eliminating drift.

Prerequisites for accurate calculation of the standard deviation are good ambient conditions in accordance with OIML R111 for an E1 mass standards laboratory.










# Calculations and Areas of Application

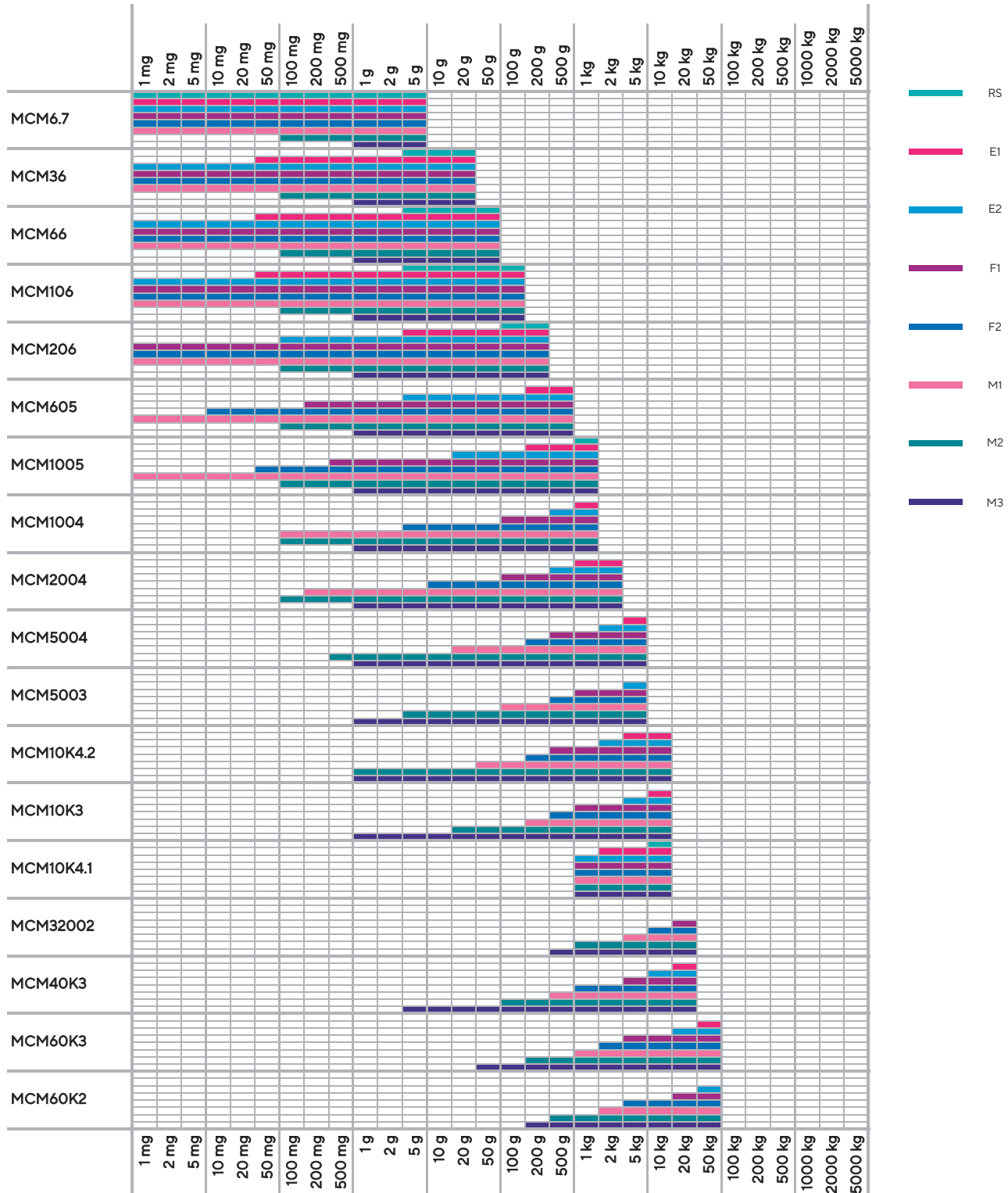
- 66 Areas of Application in Accordance with OIML R111-1:2004
- 68 Areas of Application in Accordance with ASTM E617
- 70 Calculating Uncertainties in Accordance with OIML R111

 **Find out more**  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

# Areas of Application

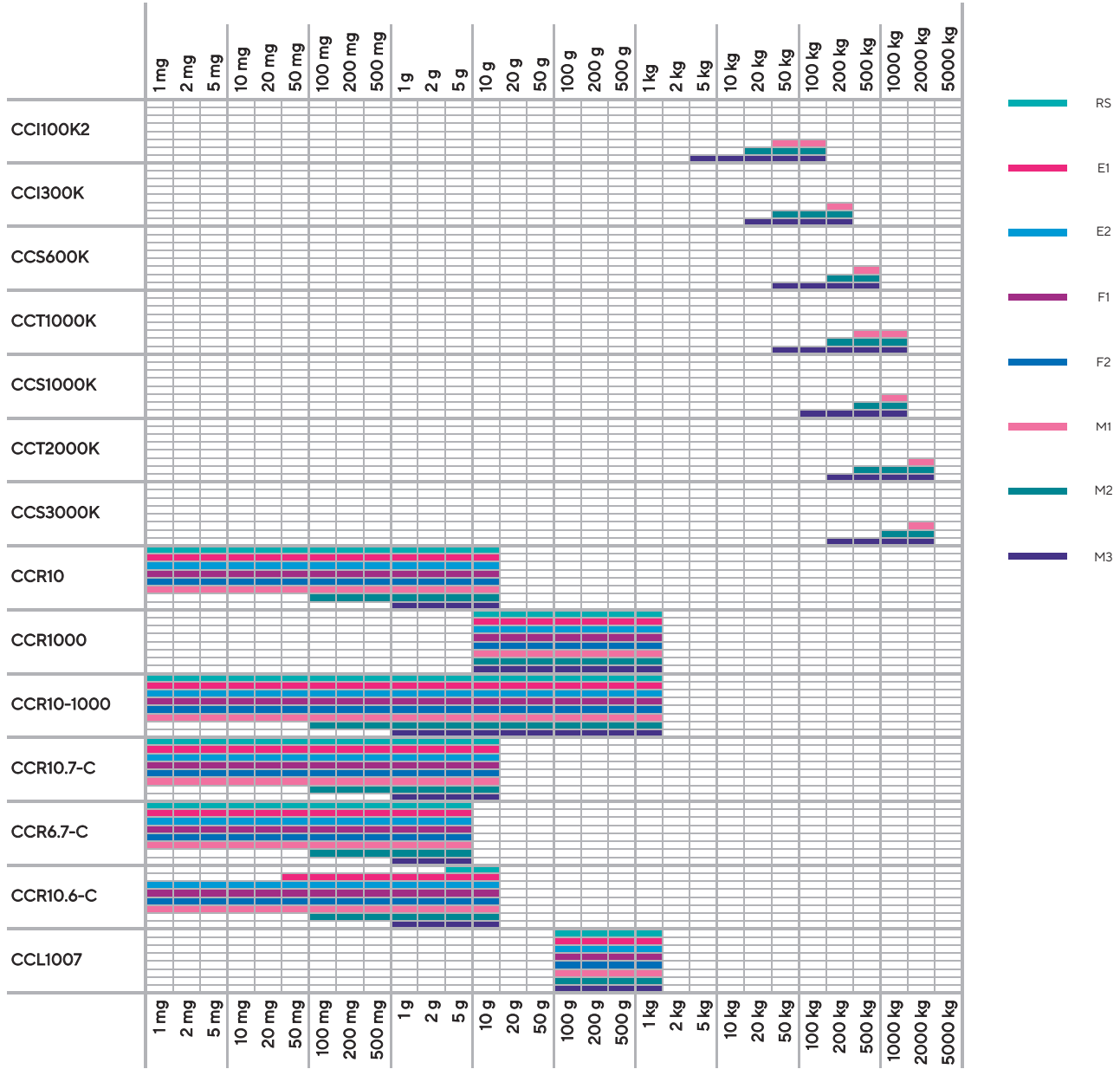
In Accordance with OIML R111-1:2004

Recommended Application Range according to OIML R111



RS: Reference Standard with 1/5 uncertainty contribution of the E1 Class tolerance limit

Recommended Application Range according to OIML R111

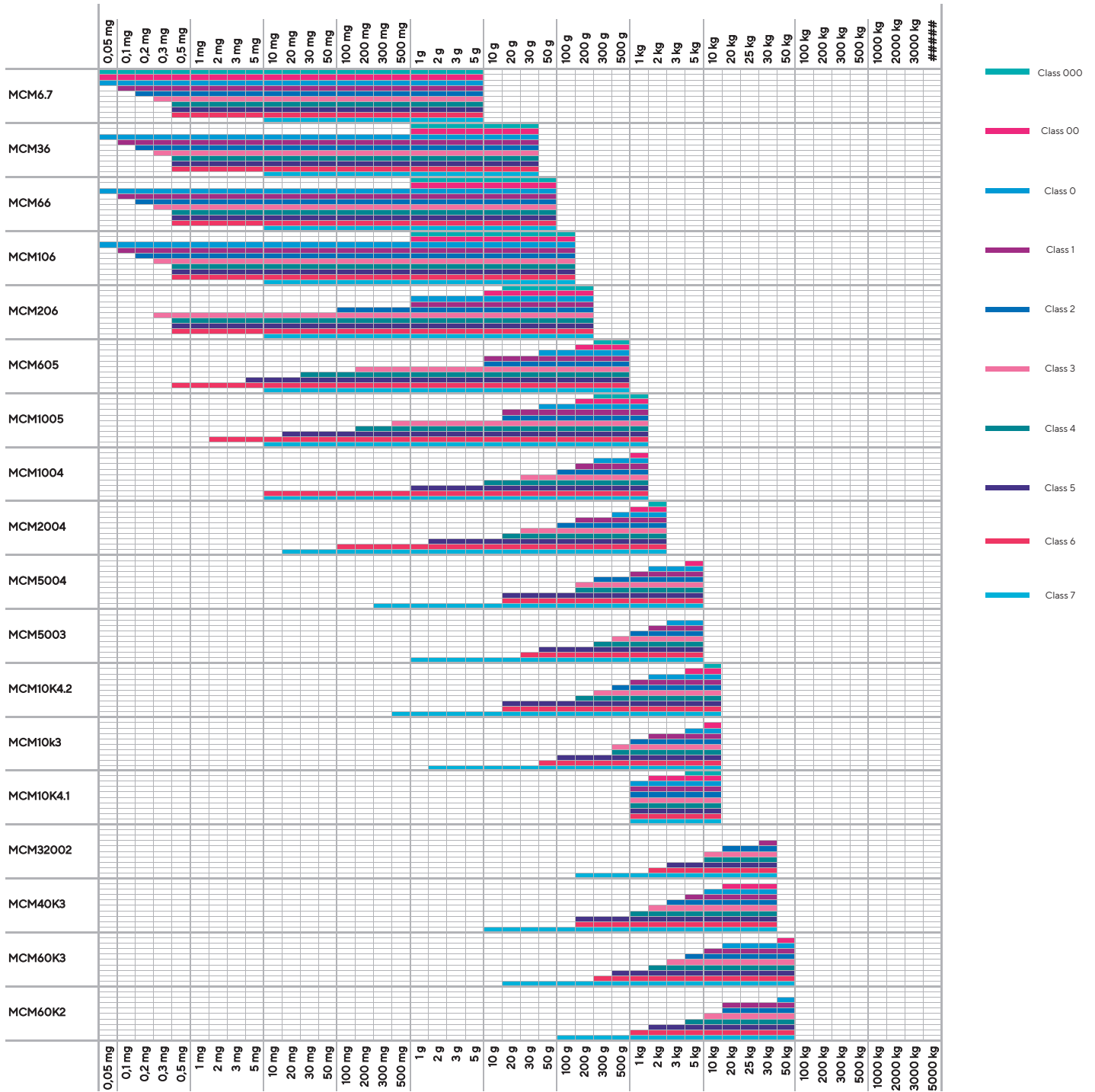


RS: Reference Standard with 1/5 uncertainty contribution of the E1 Class tolerance limit

# Areas of Application

In Accordance with ASTM E617

Recommended Application Range according to ASTM E617





# Calculating Uncertainties

## In Accordance with OIML R111

The figures "Areas of Application in Accordance with OIML R111-1 and ASTM E617-13" present an overview of the suitability of Sartorius mass comparators for mass comparisons in the various accuracy classes with a 95% confidence level ( $k = 2$ ).

The following explains the basic principles for determining these ranges.

The expanded uncertainty of the conventional mass of the test weight is produced from:

$$U(m_{ct}) = k \times u_c(m_{ct}) \quad [1, \text{Formula C.6.5-3}]$$

Each weight must fulfill this condition

$$U(m_{ct}) \leq \frac{1}{3} \text{MPE} \quad \text{cf. [1, Formula C.5.2-1, Tab. 1]}$$

(MPE: Maximum Permissible Error)

The combined uncertainty of the test weight is produced from:

$$u_c(m_{ct}) = \sqrt{u_w^2(\overline{\Delta m_c}) + u^2(m_{cr}) + u_b^2 + u_{ba}^2} \quad [1, \text{Formula C.6.5-1}],$$

from the standard uncertainty of the weighing process:

$$u_w(\overline{\Delta m_c}) = s(\Delta m_c) / \sqrt{n} \quad [1, \text{Formula C.6.1-1}],$$

from the standard uncertainty of the reference weight:

$$u(m_{cr}) = \sqrt{(U/k)^2 + u_{\text{inst.}}^2(m_{cr})} \quad [1, \text{Formula C.6.2-1}],$$

from the standard uncertainty of the air buoyancy correction with

$$\rho_{al} = \rho_0: \quad \text{cf. [1, Formula C.6.3-1], [3, Formula 34.67]}$$

$$u_b = m_{cr} \sqrt{[(\rho_r - \rho_t) / (\rho_r \times \rho_t) u(\rho_a)]^2 + (\rho_a - \rho_0)^2 (u(\rho_t)^2 / \rho_t^4 + u(\rho_r)^2 / \rho_r^4)}$$

and the combined standard uncertainty of the balance:

$$u_{ba} = \sqrt{u_s^2 + u_d^2 + u_E^2 + u_{ma}^2} \quad [1, \text{Formula C.6.4-5}].$$

This comprises the standard uncertainty due to the sensitivity of the comparator:

$$u_s = |\overline{\Delta m_c}| \sqrt{u^2(m_s) / m_s^2 + u^2(\Delta I_s) / \Delta I_s^2} \quad [1, \text{Formula C.6.4-1}, \text{ and | or simplified}]$$

$$u_s \approx |\overline{\Delta m_c}| u(m_s) / m_s \approx 5 \times 10^{-4} |\overline{\Delta m_c}| \quad \text{cf. [3, Sections 3.4.6.2-3.4.6.3],}$$

from the standard uncertainty of the display resolution:

$$u_d = d\sqrt{2}/(2\sqrt{3})$$

[1, Formula C.6.4-2],

from the standard uncertainty due to eccentricity:

$$u_E = 0$$

cf. [1, Section C.6.4.4.1]

and the standard uncertainty due to magnetism:

$$u_{ma} = 0$$

cf. [1, Section C.6.4.5]

Calculation of the air density from the climate parameters temperature, air pressure and humidity produces the standard uncertainty of the air density according to [1, Formula C.6.3-3]:

$$u(\rho_a) = \sqrt{u_T^2 + (\partial\rho_a/\partial p u_p)^2 + (\partial\rho_a/\partial t u_t)^2 + (\partial\rho_a/\partial hr u_{hr})^2}$$

with the standard uncertainty of formula:

$$u_F = 2 \times 10^{-4} \rho_a$$

cf. [1, Section E.3]

and the sensitivity coefficient:

$$\begin{aligned} \partial\rho_a/\partial p &\approx +1 \times 10^{-5} \text{Pa}^{-1} \\ \partial\rho_a/\partial t &\approx -4 \times 10^{-3} \text{K}^{-1} \\ \partial\rho_a/\partial hr &\approx -9 \times 10^{-3}, \text{ with } 0 \leq hr \leq 1 \end{aligned}$$

cf. [4, Section 2.2]

To verify the suitability of mass comparators for the various accuracy classes, uncertainty shares must be included in the uncertainty budget. It is good practice for a user to clearly apply the following:

$$u_w(\overline{\Delta m_c}) \leq \frac{4}{5} u_c(m_{ct}) \quad \text{and} \quad u(m_{cr}) = u_b = u_{ba} \leq \frac{1}{3} u_c(m_{ct})$$

If these individual requirements are observed, the combined uncertainty is always less than the required uncertainty limit.

This requirement for the standard uncertainty of the weighing process  $u_w(\overline{\Delta m_c})$  results in a similar requirement for the repeatability  $s$  of the mass comparator.

$$u_w(\overline{\Delta m_c}) = s(\Delta m_c)/\sqrt{n} \leq \frac{4}{5} u_c(m_{ct}) = \frac{2}{5} U(m_{ct}) = \frac{2}{15} \text{MPE}, \text{ i.e. } s(\Delta m_c) \leq \frac{2}{15} \text{MPE}\sqrt{n}$$

The necessary number of weighing cycles can be found in the OIML and/or ASTM for the respective accuracy class. The calculations in the diagrams for the suitability of mass comparators for the various accuracy classes are based on the respective number of ABA cycles indicated in the table.

## Example

In the following, the generation of an uncertainty budget is illustrated using the example of a calibration of a 20-kg weight of accuracy class E2 ( $n = 3$ ) on mass comparator MCM60K3. This example equally illustrates the suitability of the mass comparator for the corresponding accuracy class.

Mass comparator: MCM60K3 with calibrated internal climate module YCM20MC

Repeatability:  $s = 6 \text{ mg}$

Digital scale interval:  $d = 2 \text{ mg}$

Permissible error limit: MPE = 30 mg (20 kg, accuracy class E2)

Reference weight: Accuracy class E1, with DAkkS certificate

Conventional mass:  $m_{cr} = 20 \text{ kg} - 5 \text{ mg}$

Expanded uncertainty:  $U(m_{cr}) = 3 \text{ mg}$ ,  $k = 2$

Density:  $\rho_r = 8012.7 \text{ kg m}^{-3}$ ,  $u(\rho_r) = 1.25 \text{ kg m}^{-3}$

Test weight: Conventional mass:  $m_{ct} = 20 \text{ kg} + 15 \text{ mg}$

Density:  $\rho_t = 7950 \text{ kg m}^{-3}$ ,  $u(\rho_t) = 70 \text{ kg m}^{-3}$  [1, Tab. B.7, Stainless steel]

Ambient conditions: Measured values of the internal climate module

Temperature:  $t = 20 \text{ °C}$ ,  $u_t = 0.15 \text{ K}$

Pressure:  $p = 970 \text{ hPa}$ ,  $u_p = 1 \text{ hPa}$

Humidity:  $hr = 50 \%$ ,  $u_{hr} = 1 \%$

Air density:  $\rho_a = 1.148 \text{ kg m}^{-3}$ ,  $u(\rho_a) = 0.001 \text{ kg m}^{-3}$   
(calculated by application software on the mass comparator)



Uncertainties:

$$u_w(\overline{\Delta m_c}) = 6 \text{ mg}/\sqrt{3} = 3.46 \text{ mg}$$

$$u(m_{cr}) = 3 \text{ mg}/2 = 1.50 \text{ mg, with } u_{\text{inst.}}(m_{cr}) = 0$$

$$u_b = 1.15 \text{ mg}$$

$$u_{ba} = 0.82 \text{ mg, with}$$

$$u_s = 5 \times 10^{-4} \times 20 \text{ mg} = 0.01 \text{ mg}$$

$$u_d = 2 \text{ mg} \times \sqrt{2}/(2\sqrt{3}) = 0.82 \text{ mg}$$

$$u_E = 0 \text{ mg (already contained in } u_w)$$

$$u_{ma} = 0 \text{ mg (since weight is OIML-compliant)}$$

$$u_c(m_{ct}) = \sqrt{(3.46 \text{ mg})^2 + (1.50 \text{ mg})^2 + (1.15 \text{ mg})^2 + (0.82 \text{ mg})^2} = 4.03 \text{ mg}$$

$$U(m_{ct}) = 2 \times 4.03 \text{ mg} = \mathbf{8.06 \text{ mg}} \leq \frac{1}{3} 30 \text{ mg} = \mathbf{10 \text{ mg}}$$

The following example calculation is used to prove that the expanded uncertainty of the conventional mass of the test weight is less than 1/3 of the permissible error limit. This means the calibration undertaken with the mass comparator used is compliant with OIML R111-1.

## References


1. OIML R111-1, Weights of classes E1, E2, F1, F2, M1, M1-2, M2, M2-3 and M3, Part 1: Metrological and technical requirements, Edition 2004 (E)
2. ASTM E617-13, Standard Specification for Laboratory Weights and Precision Mass Standards, 100 Barr Harbor Drive, PO Box C700, West Conshohocken, PA 19428-2959, United States, 2013
3. Kochsiek, M., Gläser, M., "Comprehensive Mass Metrology", Wiley-VCH Verlag Berlin (2000), Sec.3.4, "Mass determination with balances" (Roman Schwartz)
4. Picard, A., Davis, R.S., Gläser, M., Fujii, K., "Revised formula for the density of moist air (CIPM-2007)", Metrologia 45 (2008) 149-155
5. GUM 1995, Guide to the expression of uncertainty in measurement, JCGM 100:2008, 2008

**Table:** Number of ABA weighing cycles

OIML class	E1	E2	F1	F2	M1	M2	M3
ASTM class	0	1 2	3 4	5	6	7	F
Cycles n	5	3	2	1	1	1	1
$s_{\max}(\Delta m_c) = \frac{2}{15} \text{MPE} \sqrt{n}$	0.30 MPE	0.23 MPE	0.19 MPE	0.13 MPE	0.13 MPE	0.13 MPE	0.13 MPE

In individual cases, it must be proven that the preset limit values of the expanded uncertainty of the conventional mass of the test weight comply.




 Find out more  
For more information, please visit  
[www.sartorius.com](http://www.sartorius.com)

**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 contacts, visit  
[www.sartorius.com](http://www.sartorius.com)