



CCR-Compact Robotic Systems

High Throughput
Mass Determination

Simplifying Progress

SARTORIUS

Table-Top Robotic Systems

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 walk-away 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.






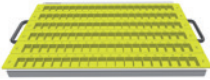
Error Free Weight Sorting

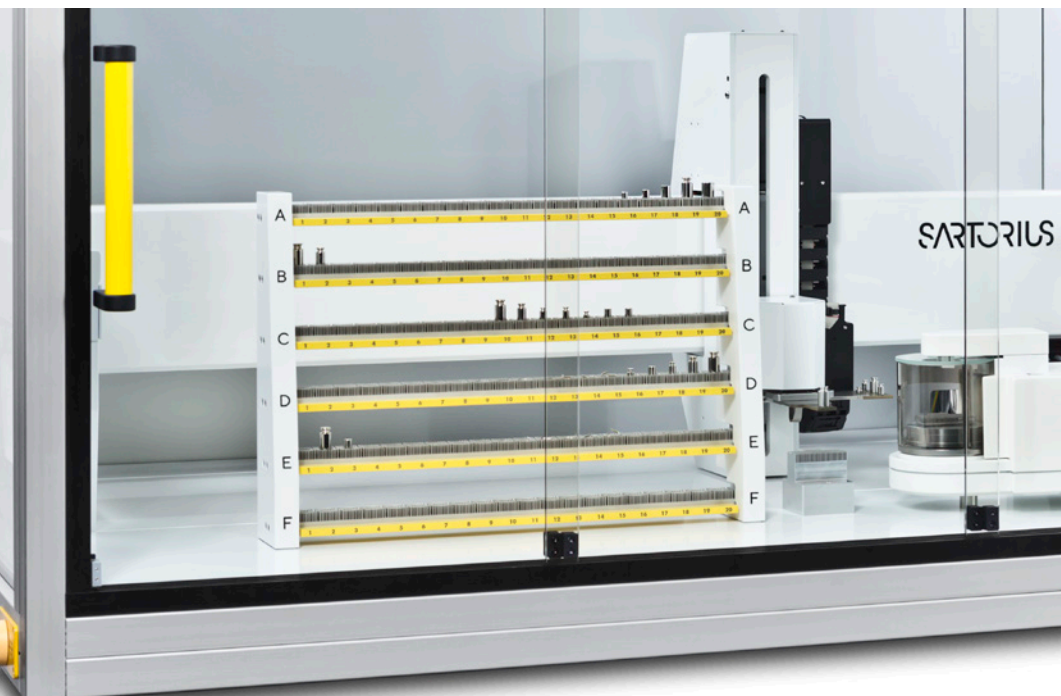
The weight sorting plate designed specially for these systems allows an easy and error free processing of the large number of weights. Each row and column is labelled and is identically to the labeling on the weight magazine. Additionally, each row is removable which further support an intuitive and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

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 x Single	1 Multi & 1 Single
Dimensions (W x D x H, mm)	1200 x 800 x 760	1200 x 800 x 760	1200 x 800 x 760

Key Accessories

Order Code		Description
YWT12		Weighing Table made of granite, weight: 680 kg, dimensions (W x D x H, mm) 800 x 1200 x 800
YWT13		Weighing Table made of granite and steel, weight: 330 kg, dimensions (W x D x H, mm) 800 x 1200 x 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



Visit our Website for More Details

- Datasheet
- White Paper

Sales and Service Contacts

For further contacts, visit
sartorius.com

Germany

Sartorius Lab Instruments
GmbH & Co. KG
Otto-Brenner-Strasse 20
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
5 Orville Drive, Suite 200
Bohemia, NY 11716
Phone +1 631 254 4249
Toll-free +1 800 635 2906