

## Moisture Trap For Cubis® II Ultra-High Resolution Balances

YCP07MC

### Compatibility:

- High-Capacity Micro Balances
- MCA/MCE36S/P,
- MCA/MCE66S/P, MCA116S
- Semi-Micro Balances  
MCA/MCE125S-3, MCA/  
MCE225S-3, MCA226S



New

# Highlights

- Fast Stabilization
- Data Integrity
- Climate Module optional

High-Capacity Micro Balances					
Product	MCA/MCE36S	MCA/MCE66S	MCA/MCE36P	MCA/MCE66P	MCA116S
Maximum Capacity	32 g	61 g	32   10.1 g	61   12 g	111 g
Readability	0.001 mg	0.001 mg	0.001   0.01 mg	0.001   0.01 mg	0.002 mg
Volume of pipetting vessel	21 mL				
Nominal volume calibration range (acc. to ISO8655:2022)*	-	From 0.5 µL to 1,000 µL		-	From 20 µL to 1,000 µL
Ingenix compatibility	Yes				

Semi-Micro Balances					
Product	MCA/MCE125S-3	MCA/MCE225S-3	MCA/MCE225P-3	MCA226S	
Maximum Capacity	120 g	220 g	120   220 g	220	
Readability	0.01 mg	0.01 mg	0.01   0.1 mg	0.005	
Volume of pipetting vessel	21 mL				
Nominal volume calibration range (acc. to ISO8655:2022)*	From 20 µL to 1,000 µL				
Ingenix compatibility	Yes				

\* without emptying the vessel during one 3x10 calibration cycle

Calibration is an important maintenance step to ensure the proper performance of your pipette during your daily work. Depending on the demands of your application a quite frequent check of pipettes in between work steps can be necessary.


With the moisture trap for Cubis® II Ultra-High Resolution Balances Sartorius provides rather convenient and comfortable accessory to convert your lab balance into a pipette calibration system.

## 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)