



Cubis Balance Sample Holders

Overview Table

Simplifying Progress



SARTORIUS

Ordering Information




Sample Holders for Cubis® II Ultra-High Resolution Balances

(Weighing modules 36S-3, 36P-3, 66S-3, 66P-3, 116S-3, 226S-3, 125S-3, 225P-3 and 225S-3)



Safe-lock Tubes

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH14-3	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Titanium alloy	Yes
YSH18-3	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Titanium alloy	No


Vial, Flask Holders

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH22-3	 Vial Holder for conical -, centrifuge-, round bottom- and test tubes up to 40 mL volume	Titanium alloy	No
YSH47-3	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Titanium alloy	Yes
YSH02-3	 Adjustable sample holder for tubes and vessels up to 40 mL	Titanium alloy	No


Filter Holders

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH35-3	 Filter Holder for up to 75 mm Ø filters	Titanium alloy	Yes
YSH30-3	 Filter Holder for up to 150 mm Ø filters	Titanium alloy	No


Stent Holders

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH12-3	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38 mm of length	Titanium alloy TA2(3.7035)	Yes

Syringe Holders

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH46-3	 Syringe Holder designed to hold analytical glass syringes typically used for in vivo applications	Titanium alloy	No


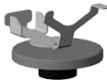
Weighing Boat Holder

Order Code	Description	Material	Compatible with inner draft shield (YDS125A or YDS125U)
YSH26-3	 Weighing Boat Holder to weigh weighing boats of up to 40 x 60 mm	Titanium alloy	No



Sample Holders for Cubis® II Semi-Micro and Analytical Balances

(Weighing modules MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S)**


Safe-lock Tubes

Order Code	Description	Material
YSH15	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Titanium alloy
YSH19	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Titanium alloy


Vial, Flask Holders

Order Code	Description	Material
YSH23	 Vial Holder for conical -, centrifuge-, round bottom- and test tubes up to 40 mL volume	Titanium alloy
YSH37	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Titanium alloy


Filter Holders

Order Code	Description	Material
YSH30	 Filter Holder for up to 150 mm Ø filters	Titanium alloy

Syringe Holders

Order Code	Description	Material
YSH46	 Syringe Holder designed to hold analytical glass syringes typically used for in vivo applications	Titanium alloy

Weighing Boat Holder


Order Code	Description	Material
YSH26	 Weighing Boat Holder to weigh weighing boats of up to 40 x 60 mm	Titanium alloy

** Please note that weighing modules 225S/P and 125S are discontinued. Corresponding sample holders are still available.




Sample Holders for Cubis® II Ultramicro- and Micro Balances

(Weighing modules MC* 2.7S; 6.6S; 3.6P; 10.6S)


Safe-lock Tubes

Order Code	Description	Material
YSH13	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Titanium alloy

Filter Holders

Order Code	Description	Material
YSH34	 Filter Holder for up to 50 mm Ø filters	Titanium alloy
YSH35	 Filter Holder for up to 75 mm Ø filters	Titanium alloy
YSH36	 Filter Holder for up to 90 mm Ø filters	Titanium alloy


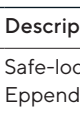
Stent Holders

Order Code	Description	Material
YSH10	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38mm of length	Titanium alloy TA2(3.7035)

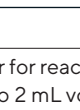
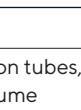
Sample Holders for Cubis® High-Capacity Micro Balances

(Weighing modules MS* 36S/P; 66S/P; 116P)**


Safe-lock Tubes

Order Code	Description	Material
YSH14	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Titanium alloy
YSH18	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Titanium alloy


Vial, Flask Holders

Order Code	Description	Material
YSH22	 Vial Holder for conical-, centrifuge-, round bottom- and test tubes up to 40 mL volume	Titanium alloy
YSH47	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Titanium alloy

Stent Holders

Order Code	Description	Material
YSH12	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38 mm of length	Titanium alloy TA2(3.7035)

Syringe Holders

Order Code	Description	Material
YSH41	Syringe Holder designed to hold analytical glass syringes typically used for in vivo applications	Titanium alloy
YSH02	 Adjustable sample holder for tubes and vessels up to 40 mL	Titanium alloy

** Please note that Cubis High-Capacity Micro Balances are discontinued. Corresponding sample holders are still available.

Germany


Sartorius Lab Instruments GmbH & Co. KG
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Corporation
North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

UK

Sartorius UK Ltd.
Longmead Business Centre
Blenheim Road
Epsom
Surrey
KT19 9QQ
Phone +44 1372 73710

 **For further contacts, visit:**
www.sartorius.com