



Cubis[®] Sample Holders

Ergonomic Sample Weighing Without Sample Loss and Headaches

Simplifying Progress

SARTORIUS

Cubis® Sample Holders: Your Advantages

While standard weighing pans are the most commonly used surfaces during the typical weighing processes, the pans reach their limit as soon as samples, vials, materials with unique form or for special applications need to be used.

Whether you want to weigh directly into safe-lock tubes, conically shaped vials, round bottomed flasks, filters, syringes or stents etc., ergonomics and accuracy of your weighing process has the ultimate importance.

Cubis® Sample Holders from Sartorius are specially designed weighing pans which enable ergonomic sample weighing in workflows where using a standard pan is difficult or impossible. The usage of Cubis Sample Holders not only simplifies your weighing process, but results in better repeatability as well.

Uncompromised Accuracy

Made of titanium, not only the highest material quality is guaranteed, but notably the repeatability is not jeopardized because of the non-magnetic properties

Convenient Usage

No more powder stuck to the weighing boat, weighing paper, etc. No more tip over. Thanks to ergonomic design weighing workflows can be executed much easier.

Great Variety of Choice

Selecting from numerous design variants you can make sure that the sample holder fits the balance and the application type you are looking for.

Customized Solutions

In the rare case that your sample holder needs are not covered in our portfolio, a fully personalized solution can be developed for you.

“The Safe-Lock Tube Holder for Cubis® II Micro Balances is easy to use and can fit the small tube directly. We can perform small amount of weighing easily and save a lot of reference standard and time as well.”

Nurul Hammizah Bt Hamidon
Industrial Biotechnology
Research Centre, SIRIM Berhad, Malaysia

“The YSH23 vial holder from Sartorius has the perfect ergonomics and enables us to perform hydrolytic resistance measurements at the highest level of precision.”

Otto Molnar
Quality Engineering
BD, Hungary



Cubis[®] Sample Holders: Great Variety of Choice



Safe-Lock Tube Holders

No more sample loss...even on a micro balance.

Unless a specific solution provided, for highly sensitive weighing workflows typically paper or weighing boats are used in the first step and then samples are transferred into safe-lock tubes. This may result in sample loss which could negatively affect the outcome. Cubis[®] Safe-Lock Tube holders are specially designed weighing pans which enable direct sample weighing into test tubes. Using these holders the risk of sample loss can be completely eliminated which ensures high measurement accuracy. Design variants for tubes between 1,5 mL and 5 mL are available and can be used in micro balances, high capacity micro-balances, semi-micro and analytical balances accommodating to your specific needs.



Vial- and Flask Holders

No more headaches.

Conically shaped tubes, round bottomed flasks are widely used in analytical-, chemical- and QC laboratories. Because of their special form it is difficult to place these type of flasks directly on a flat weighing pan. Sartorius designed special weighing pans to hold these flasks whilst not compromising the accuracy of the weighing results. Diverse types for high capacity micro balances, semi-micro and analytical balances are available to support the comfort of use.



Filter Holders

Fulfill your regulatory requirements

During exhaust gas -, air sampling or emission testing the determination of residual weight or particulate matter is a typical workflow. Sartorius offers various sizes of filter holders from 50 mm - up to 150 mm diameter for ultra-micro and micro filter balances, and semi-micro balances. Made of titanium, not only the highest material quality is guaranteed, but notably the repeatability is not compromised because of the non-magnetic properties.

Syringe Holders

Safety at first place

In areas of pharma, biopharma, cosmetics, medical devices or research facilities syringes are weighed on a daily basis. Ergonomic holding of syringes is crucial not only because of better handling but also for safety reasons. Syringe Holders from Sartorius were designed to hold standard sized analytical glass syringes (up to 2,5 mL) as well as syringes used for in vivo applications (e.g. insulin syringes) and are available for Cubis® High Capacity Micro Balances and Cubis® II Semi-Micro and Analytical Balances. Made of titanium, not only high quality is guaranteed, but better repeatability due to its non-magnetic properties.

Cubis® Custom Made Solutions in Just Five Steps

In case your sample holder needs are not covered in our portfolio, a fully personalized solution can be developed for you.



Define needs
Identify and specify your requirements



Contact us



Development & manufacturing phase
Production of your personalized solution



Consulting phase
We will work closely with you to support realizing your goals



Delivery
As soon as your request is ready



Weighing Boat Holder

Straightforward weighing

Weighing Boat Holders are easy to assemble on a Cubis® Semi-Micro and Analytical Balances, allowing direct dosing into weighing boats. Titanium material guarantees better repeatability because of its non-magnetic properties.




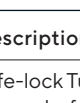


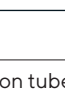
Stent Holders

Increased reproducibility and patient safety

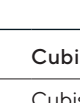

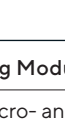

When it comes to stent weighing the standard weighing pan might not support your weighing workflows well enough. These medical products need to be weighed at the highest level of precision while handling them with caution and great delicacy which is difficult to execute using a flat weighing pan. Not only precision and carefulness play a role, but other factors as well, for ex. contact with particular materials might be avoided (e.g. contact with Aluminium). Stent Holders from Sartorius are made of Grade 2 Titanium which not only improves repeatability but is Aluminium free too. Specific holders for Cubis® ultra-micro, micro and high capacity micro balances are available for ergonomic weighing of coronary-, neurovascular- or small vascular stents up to 38 mm length. For bigger sized stents we offer custom made solutions.

Ordering Information





Safe-lock Tubes

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH13	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Cubis® II Ultramicro- and Micro Balances Models: MC* 2.7S; 6.6S; 3.6P; 10.6S
YSH14	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Cubis® High Capacity Micro Balances Models: MS* 36S/P; 66S/P; 116P
YSH15	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S
YSH18	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Cubis® High Capacity Micro Balances Models: MS* 36S/P; 66S/P; 116P
YSH19	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S



Vial, Flaks Holders

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH22	 Vial Holder for conical -, centrifuge-, round bottom- and test tubes up to 40 mL volume	Cubis® High Capacity Micro Balances Models: MS* 36S/P; 66S/P; 116P
YSH23	 Vial Holder for conical -, centrifuge-, round bottom- and test tubes up to 40 mL volume	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S
YSH37	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S
YSH47 (please check availability)	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Cubis® High Capacity Micro Balances Models: MS* 36S/P; 66S/P; 116P



Filter Holders

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH34	 Filter Holder for up to 50 mm Ø filters	Cubis® II Ultramicro- and Micro Balances Models: MC* 2.7S-F; 6.6S-F; 3.6P-F; 10.6S-F
YSH35	 Filter Holder for up to 75 mm Ø filters	Cubis® II Ultramicro- and Micro Balances Models: MC* 2.7S-F; 6.6S-F; 3.6P-F; 10.6S-F
YSH36	 Filter Holder for up to 90 mm Ø filters	Cubis® II Ultramicro- and Micro Balances Models: MC* 2.7S-F; 6.6S-F; 3.6P-F; 10.6S-F
YSH30	 Filter Holder for up to 150 mm Ø filters	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S


Stent Holders

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH10 (please check availability)	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38mm of length	Cubis® II Ultramicro- and Micro Balances Models: MC* 2.7S; 6.6S; 3.6P; 10.6
YSH12 (please check availability)	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38mm of length	Cubis® High Capacity Micro Balance Models: MS* 36S/P; 66S/P; 116P

Syringe Holders

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH46 (please check availability)	 Syringe Holder designed to hold analytical glass syringes and plastic syringes (typically used for in vivo applications)	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S
YSH41 (please check availability)	 Syringe Holder designed to hold analytical glass syringes and plastic syringes (typically used for in vivo applications)	Cubis® High Capacity Micro Balances Models: MS* 36S/P; 66S/P; 116P

Weighing Boat Holder

Order Code	Description	Cubis® Weighing Modules (Compatibility)
YSH26	 Weighing Boat Holder to weigh weighing boats of up to 40 x 60 mm	Cubis® II Semi-Micro and Analytical Balances Models: MC* 225S/P; 125S/P; 524S/P; 324S/P; 224S; 124S

Sales and Service Contacts

For further contacts, visit
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