



# Cubis® II Ultra- High Resolution Balances

## Sample Holders

Ergonomic Sample

Weighing Without Sample

Loss and Headaches

Simplifying Progress

**SARTORIUS**

# Cubis® II Ultra-High Resolution 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® Ultra-High Resolution 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 Ultra-High Resolution 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

## 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.

## Convenient Usage

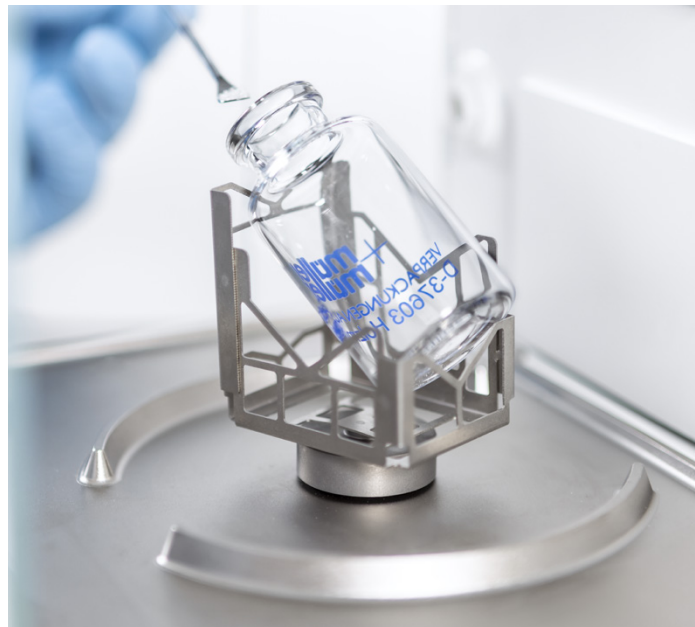
No more powder stuck to the weighing boat, weighing paper, etc. No more tip over. Thanks to ergonomic design weighing workflows can be executed much easier.

## 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.



# Ultra-High Resolution 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 Ultra-High Resolution 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 Ultra-high Resolution 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-High Resolution 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® II Ultra-High Resolution 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**



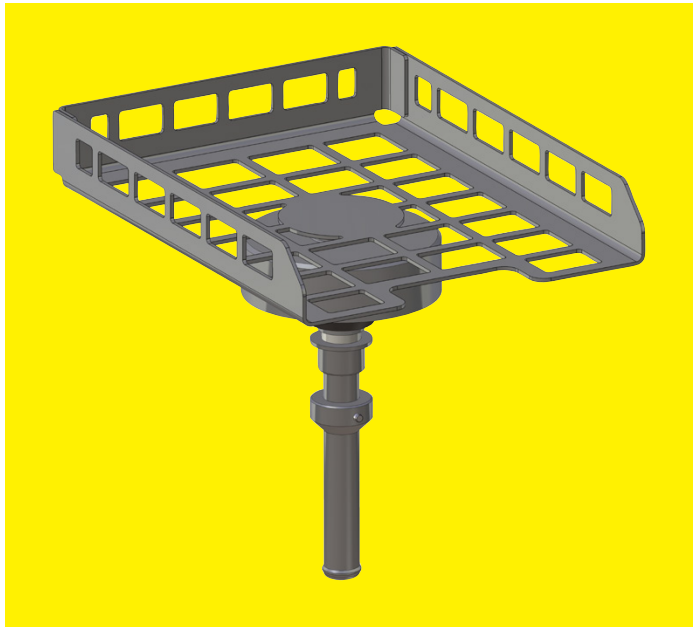
**Development & manufacturing phase**  
Production of your personalized solution



**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® II Ultra-High Resolution Balance, 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. A specific holder for Cubis® II Ultra-High Resolution Balances is 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

Sample holders for Cubis® II Ultra-High Resolution Balances  
(weighing modules 36S-3, 36P-3, 66S-3, 66P-3 and 116S-3)


## Safe-lock Tubes

Order Code	Description	Material
YSH14-3	 Safe-lock Tube Holder for reaction tubes, PCR tubes, Eppendorf tubes up to 2 mL volume	Titanium alloy
YSH18-3	 Safe-lock Tube Holder for bigger reaction- and Eppendorf tubes up to 5 mL volume	Titanium alloy


## Vial, Flaks Holders

Order Code	Description	Material
YSH22-3	 Vial Holder for conical -, centrifuge-, round bottom- and test tubes up to 40 mL volume	Titanium alloy
YSH47-3	 Holder for titration vessels, round bottom flasks and test tubes with diameter up to 50 mm	Titanium alloy
YSH02-3	 Adjustable sample holder for tubes and vessels up to 40 mL	Titanium alloy


## Filter Holders

Order Code	Description	Material
YSH35-3	 Filter Holder for up to 75 mm Ø filters	Titanium alloy
YSH30-3	 Filter Holder for up to 150 mm Ø filters	Titanium alloy


## Stent Holders

Order Code	Description	Material
YSH12-3	 Coronary Stent Holder designed to weight coronary-, neuro vascular- or small vascular stents up to 38mm of length	Titanium alloy TA2(3.7035)

## Syringe Holders

Order Code	Description	Material
YSH46-3	 Syringe Holder designed to hold analytical glass syringes typically used for in vivo applications	Titanium alloy

## Weighing Boat Holder

Order Code	Description	Material
YSH26-3	 Weighing Boat Holder to weigh weighing boats of up to 40 x 60 mm	Titanium alloy

**Germany**


Sartorius Stedim Biotech GmbH  
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 Stedim Lab Ltd.  
Sperry Way, Stonehouse Park  
GL10 3UT Stonehouse,  
Gloucestershire, UK  
Phone +44 1453 821972

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