

# Mycap<sup>®</sup> CCX Flask Stand

Stabilize Flasks During  
Cell Passaging

## Benefits

- Prevent tipping of flasks during cell passaging
- No product contact
- Durable materials for multiple uses



## Product Information

Aseptic cell passaging with Mycap<sup>®</sup> CCX allows media and culture to be transferred aseptically by peristaltic pump. Flasks can easily tip from handling or the weight of tubing and fittings. The Mycap<sup>®</sup> CCX Flask Stand stabilizes the flask on the balance during the transfer to enhance accuracy and simplify handling.

## Applications

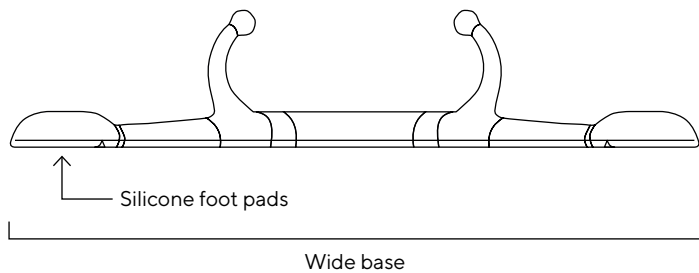
The Mycap® CCX Flask Stand is a reusable accessory that sits atop the balance pan and stabilizes Mycap® CCX flasks.

### Design Principle

Mycap® CCX have integral tubing so transfer of media and culture into a flask can be done aseptically outside a biosafety cabinet. Flasks become unstable when the center of gravity moves outside the base of the flask. The center of gravity is mobilized during handling of tubing causing the flask to become unstable and slide or tip.

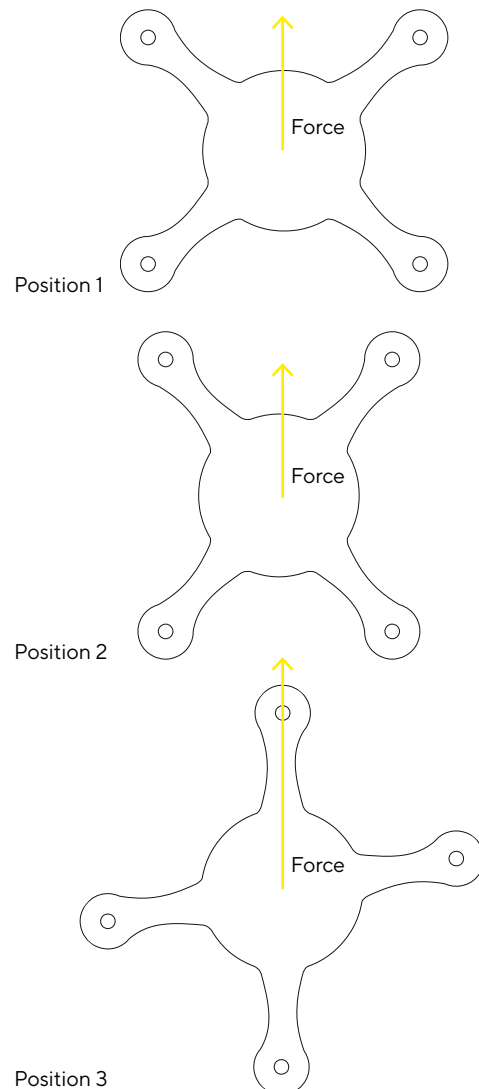
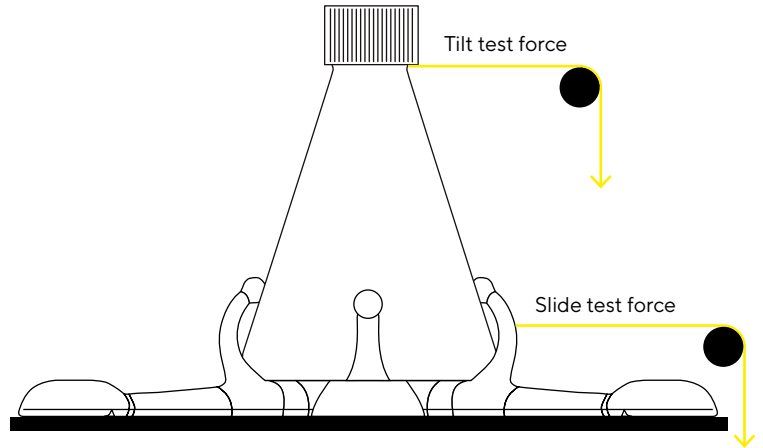
### The Mycap® CCX Flask Stand restabilizes the flask by two mechanisms:

- The stand creates a significantly wider base thus increasing the stability.
- The four feet of the stand have silicone pads. The pads are slightly concave to create suction so the stand resists movement.



## Performance

Sartorius conducted a Slide Test and Tilt Test to quantify the performance of the flask stand. Force was applied from three different directions. Three replicates were performed on flasks held in the Mycap® CCX Flask Stand and not in the stand, as a control.



## Materials of Construction

Mycap® CCX Flask Stand is a 3D printed multi-use accessory. The Flask Stand is not product contact and is constructed of:

- Platinum-Cured Silicone
- MPU 100 (Multi-purpose polyurethane)

The Mycap® CCX Flask Stand may be wiped down with 70% IPA or water. Routinely wipe down feet to restore tackiness of silicone pads.

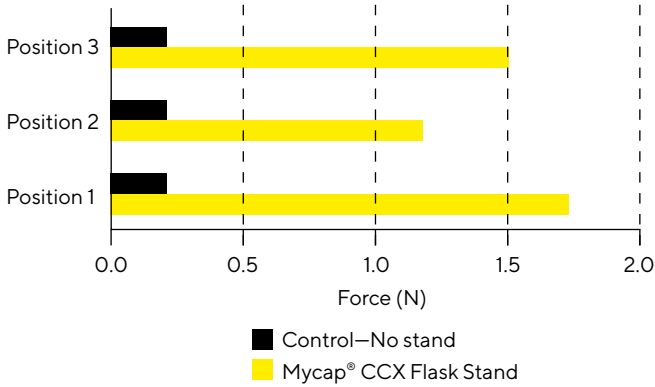
The Mycap® CCX Flask Stand cannot be autoclaved.

Recommended operating temperatures:  
15–25 °C (60–75 °F)

### Results—Slide Test

The Mycap® CCX Flask Stand increases the force required to slide a flask by an average of 704% from the control.

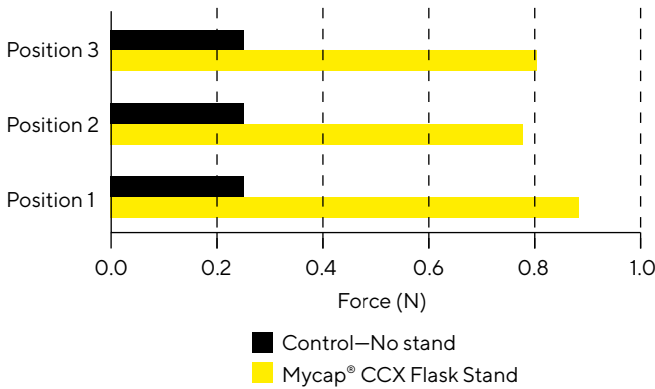
#### Slide Test



### Results—Tilt Test

The Mycap® CCX Flask Stand increases the force required to tilt a flask by an average of 331% from the control.

#### Tilt Test

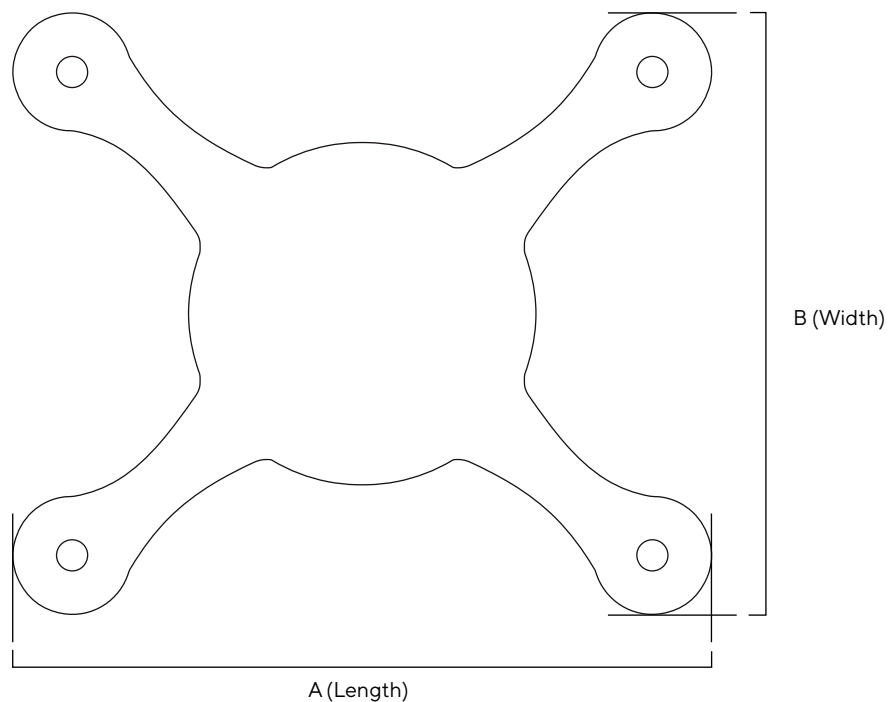


The Mycap® CCX Flask Stand meets the requirements to stabilize the flask during use.



# Ordering Information and Dimensions

Order Code	Description	A (Length)	B (Width)
MCX0250STDCRN	Mycap® CCX Flask Stand; 250 mL Corning Erlenmeyer Flask	7.73" (196.3 mm)	6.66" (199.2 mm)
MCX0500STDCRN	Mycap® CCX Flask Stand; 500 mL Corning Erlenmeyer Flask	7.73" (196.3 mm)	6.66" (199.2 mm)
MCX1000STDCRN	Mycap® CCX Flask Stand; 1,000 mL Corning Erlenmeyer Flask	7.73" (196.3 mm)	6.66" (199.2 mm)



## Germany

Sartorius Stedim Biotech GmbH  
August-Spindler-Strasse 11  
37079 Goettingen  
Phone +49 551 308 0

## USA

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

 For more information, visit

[www.sartorius.com](http://www.sartorius.com)