



sartorius



8

in one  
go!

HPLC Sample Preparation with the  
Claristep<sup>®</sup> Filtration System

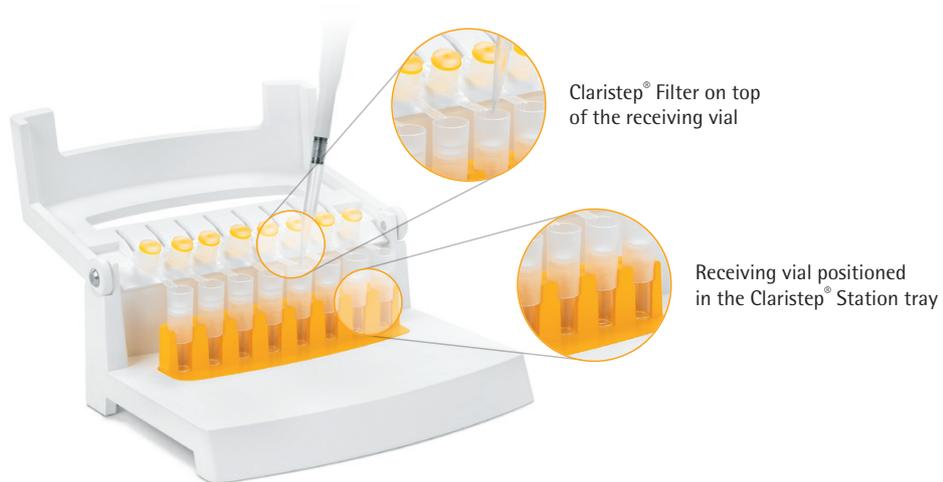
# Claristep<sup>®</sup> Filtration System



What makes the Claristep<sup>®</sup> Filtration System so easy to use?  
The Power of Simplicity!

Do you process dozens of samples each day? This syringeless solution will help you reduce time, effort and waste – and minimize hand strain.

- **Less time**  
Filter 8 samples simultaneously – without needing any power supply or a vacuum | pressure source.
- **Less stress**  
Simply place the filters on your vials, gently close the station and press on the station lid to filter – that's it!
- **Less waste**  
Filter up to 600 µL sample volumes without the use of syringes before you perform analysis, such as HPLC.



The Most Ergonomic Clarification  
How does it work?



Close the station lid. The grooves align the caps automatically, securely sealing every single Claristep<sup>®</sup> Filter unit for the most convenient processing.



Press down on the station lid so that the left and right corners touch the base plate. Hold the lid in place for 3 seconds to ensure all sample liquid is filtered through.

What if...

...you need more than 600 µL sample volumes?

Is this simply because you need a minimum filling height in your receiving vial?  
In this case, you can use vials with an insert or thick-walled vials to raise the liquid level.  
This also makes Claristep® suitable for low-volume samples.



### Claristep® Components and Ordering Information

EFD <sup>1</sup>	Membrane	Housing	Pore Size	Sterile	Qty.   Pkg.	Order No.
9.7 mm	RC	PP	0.2 µm	No	96	17C07FT---96
9.7 mm	RC	PP	0.2 µm	No	480	17C07FT--480
9.7 mm	RC	PP	0.45 µm	No	96	17C06FT---96
9.7 mm	RC	PP	0.45 µm	No	480	17C06FT--480
Claristep® Station, complete					1	17C--M8
Claristep® Single Tray					1	17C--S1

<sup>1</sup> Effective Filtration Diameter  
RC = Regenerated Cellulose

2.



Apply slight uniform pressure with your hand to start sample clarification. You will feel a certain resistance while liquid is pressed through each membrane.

4.



Claristep® Filter units press liquid through each membrane by an air pocket that forms over each filter unit when the station lid is closed. This air pocket is released when you stop holding down the lid – you will feel it in your fingertips!



## Contact

Sartorius Lab Instruments  
GmbH & Co. KG  
Otto-Brenner-Strasse 20  
37079 Goettingen

Phone +49.551.308.0  
Fax +49.551.308.3289



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

