



Venting Needle

for Ampoules, Collapsible Bags and Vials

Sterile Venting



Typical performance data

Filter area	2.0 cm ²
Diameter housing	25 mm
Diameter needle	1.60 mm
Total length	75 mm
Needle length	40 mm
Air Flow at 100 mbar	1.3 L/min
Water penetration point	>3.0 bar
Bubble Point (IPA)	> 1.0 bar
Sterility	Gamma sterile single packed
Order #	16596-HNK

Typical applications

- Venting of glass vials with rubber septum tops
- Venting of rigid plastic vials
- Use in sterility testing of ampoules or collapsible bags (during the process of rinsing or media filling)
- Venting for vial-to-vial aseptic transfer of large volume parenterals

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.91.3586102
Japan +81.3.3740.5407

Specifications subject to change without notice. Printed and copyrighted by Sartorius Stedim Biotech GmbH
W - G
Publication No.: SL-2025-e09103
Order No.: 85030-522-27
Ver. 10 | 2009