

User Guide

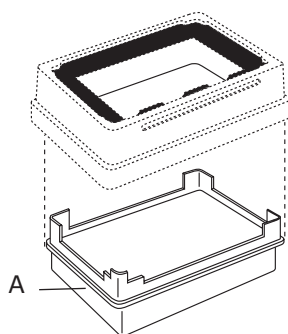
Collar Holder

Accessory for the MultiScreen[®]_{HTS} Vacuum Manifold

MSVMHTS0H

Collar Holder

The collar holder accessory (A) is recommended for use on most robotic systems and creates a stable position for placement of the manifold collar during assembly/disassembly routines.



Specifications

Materials of construction

Collar holder Acetal plastic

Collar holder dimensions length × width × height

Footprint	12.7 × 8.6 cm (5.0 × 3.4 in.)
Overall	14.2 × 10.2 × 5.1 cm (5.6 × 4.0 × 2.0 in.)

Product Ordering

Products can be purchased at SigmaAldrich.com.

Description	Cat. No.	Qty/Pk
Collar Holder	MSVMHTS0H	1
MultiScreen [®] _{HTS} Vacuum Manifold	MSVMHTS00	1

Notice

We provide information and advice to our customers on application technologies and regulatory matters to the best of our knowledge and ability, but without obligation or liability. Existing laws and regulations are to be observed in all cases by our customers. This also applies in respect to any rights of third parties. Our information and advice do not relieve our customers of their own responsibility for checking the suitability of our products for the envisaged purpose.

The information in this document is subject to change without notice and should not be construed as a commitment by the manufacturing or selling entity, or an affiliate. We assume no responsibility for any errors that may appear in this document.

Technical Assistance

Visit the tech service page on our web site at SigmaAldrich.com/TechService.

Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at SigmaAldrich.com/Terms.

Contact Information

For the location of the office nearest you, go to SigmaAldrich.com/Offices.