

User Guide

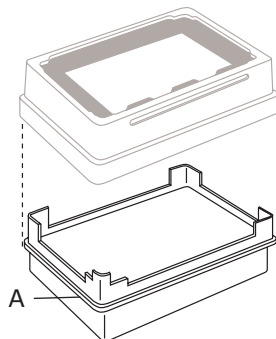
Collar Holder

Accessory for the MultiScreen®_{HTS} Vacuum Manifold

MSVMHTS0H

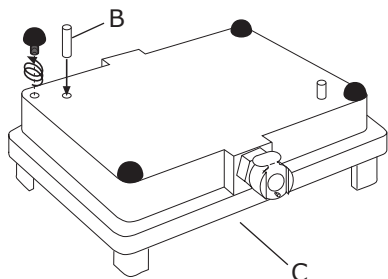
Collar Holder

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. It comes with detent pins that can be used to adapt the MultiScreen®_{HTS} Vacuum Manifold base (see next section).



Adapting the MultiScreen®_{HTS} Vacuum Manifold Base for Beckman® Systems

Detent pins (B) are supplied with the collar holder for use with Beckman® robotic systems. After removing the rubber feet from the manifold base (C), place the pins in the holes on the underside of the base to hold it in place on the robotic deck.



Specifications

Materials of construction

Collar holder	Acetal plastic
Detent pins	Stainless steel

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 Information

This section lists the catalogue numbers for the collar holder and the MultiScreen®_{HTS} Vacuum Manifold. See the Technical Assistance section for contact information. You can purchase these products on-line at SigmaAldrich.com.

Description	Catalogue 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.

Merck, Millipore, Multiscreen, and Sigma-Aldrich are trademarks of Merck KGaA, Darmstadt, Germany or its affiliates. All other trademarks are the property of their respective owners. Detailed information on trademarks is available via publicly accessible resources.
© 2004-2024 Merck KGaA, Darmstadt, Germany and/or its affiliates. All Rights Reserved.

The life science business of Merck operates
as MilliporeSigma in the U.S. and Canada.

P36558ug Rev 09/24

