

## User Guide

# Deep Well Collar

Accessory for the MultiScreen®<sub>HTS</sub> Vacuum Manifold

### MSVMHTS0D

The deep well collar is designed to accommodate the use of 2 mL deep well collection plates with MultiScreen® Deep Well Solvinert™ filter plates on the MultiScreen®<sub>HTS</sub> Vacuum Manifold. For collection plates handling other volumes, contact Technical Service.

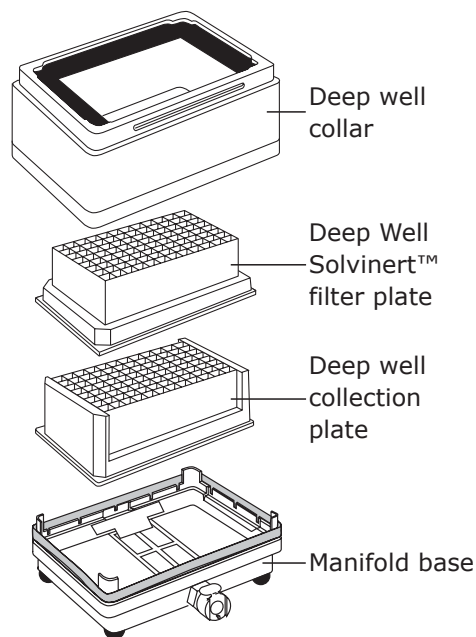
## Manifold Configurations for Use with the Deep Well Collar

The figures below show how to set up the MultiScreen®<sub>HTS</sub> Vacuum Manifold for both filtrate collection (DirectStack™ mode) and filter-to-waste applications.

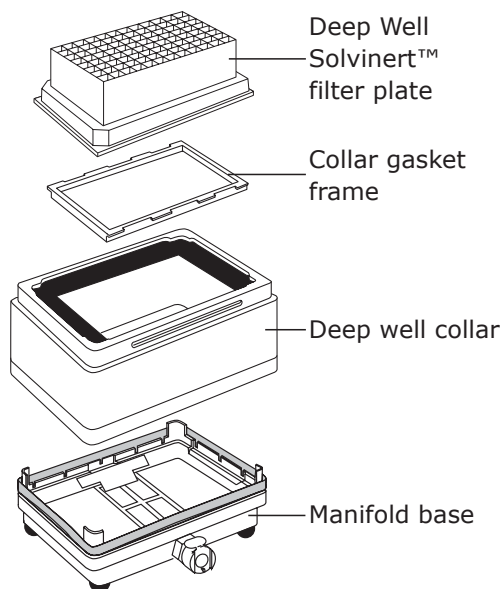
**Note:** Remove the collar gasket frame when running in Direct Stack™ mode.

Ensure that the collar gasket frame is in place for filter-to-waste mode.

### DirectStack™ mode



### Filter-to-waste mode



## Specifications

Chemical compatibility for the Deep Well Collar can be found in publication number PR4772EN, available at [SigmaAldrich.com](https://www.sigmaaldrich.com) (enter the publication number in the search box). The compatibility is the same as for the Standard Collar.

---

### Materials of construction

Deep well collar	Nylon
Gasket	Silicone

---

Deep well collar dimensions	length × width × height
	15.0 cm × 10.9 cm × 6.9 cm
	(5.9 in. × 4.3 in. × 2.7 in.)

---

Assembled height of manifold with deep well collar in place	8.6 cm (3.4 in.), excluding plate and feet
-------------------------------------------------------------	--------------------------------------------

EU/device pressure rating temperature	≤ 2.15
---------------------------------------	--------

Flow rate temperature	21 °C
-----------------------	-------

## Product Ordering Information

This section lists the catalogue numbers for the Deep Well Collar and associated products for use with the MultiScreen<sup>®</sup><sub>HTS</sub> Vacuum Manifold. See the Technical Assistance section for contact information. You can purchase these products on-line at [SigmaAldrich.com](https://www.sigmaaldrich.com).

---

Description	Cat. No.	Qty/Pk
Deep Well with Collar	MSVMHTS0D	1
MultiScreen <sup>®</sup> <sub>HTS</sub> Vacuum Manifold	MSVMHTS00	1
MultiScreen <sup>®</sup> Deep Well Solvinert <sup>™</sup> Plate, 0.45 µm, phobic, non-sterile	MDRLN0410	10
MultiScreen <sup>®</sup> Deep Well Solvinert <sup>™</sup> Plate, 0.45 µm, phobic, non-sterile	MDRPN0410	10
MultiScreen <sup>®</sup> Deep Well Solvinert <sup>™</sup> Plate, 0.45 µm, phobic with prefilter, non-sterile	MDRPNP410	10
MultiScreen <sup>®</sup> Deep Well Collection Plate, 96-well, non-sterile	MDCPNM50	50

---

## 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](https://SigmaAldrich.com/TechService).

## Terms and Conditions of Sale

Warranty, use restrictions, and other conditions of sale may be found at [SigmaAldrich.com/Terms](https://SigmaAldrich.com/Terms).

## Contact Information

For the location of the office nearest you, go to [SigmaAldrich.com/Offices](https://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.  
© 2005-2024 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.

P36537ug Rev 09/24

