Vacuum manifold

Optimized for the vacuum filtration of multi-well plates. The manifold is constructed of a chemical-resistant polymer and is designed to facilitate the filtration of liquid through a multi-well filter plate.

Ordering Information

Order Online »

ADI-987-500 1ea

Manuals, SDS & CofA

View Online »



Handling & Storage

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Application Notes For multiplexing.

Last modified: May 29, 2024



info-

eu@enzolifesciences.com