

Operating Instructions | Betriebsanleitung | Mode d'emploi |
Instrucciones de manejo | Manuale d'uso

16842

Combisart Manifold | Combisart Absaugleiste |
Rampe Combisart | Colector Combisart |
Collettore Combisart

© 2022
Sartorius Lab Instruments GmbH & Co. KG
Otto-Brenner-Str. 20
37079 Goettingen, Germany

JW | Publication No.: SL-6158-p221004

85032-535-37



SARTORIUS

Dear Customer,

thanks for deciding for a Combisart manifold with sterile vented filter station(s). Releasing the vacuum under sterile conditions after filtration, prevents secondary contamination on the underside of the membrane filter.

The graphic below shows the four different positions of the tap:

Valve
Position
Open
6 o'clock

Valve Position
Vent | Close
9 o'clock

Valve
Position
Finish
3 o'clock

Valve
Position
Autoclave
12 o'clock



EN

For Filtration

The full vacuum draws the sample through the membrane filter. The vent filter is "off-line."

After Filtration

The vacuum between the valve and membrane filter is released under sterile conditions, without any particles being drawn into the system. Secondary contamination of the bottom of the filter is ruled out entirely.

After the Filtration Run

The residual vacuum between the pump and valve is released via the sterilizing-grade filter.

For Autoclaving

For reliable sterilization, the steam flows freely through all openings.

Valve
Position
Open
6 o'clock



Valve Position
Vent | Close
9 o'clock



Valve
Position
Finish
3 o'clock



Valve
Position
Autoclave
12 o'clock



DE

Zur Filtration

Das volle Vakuum zieht die Probe durch das Membranfilter. Das Belüftungsfilter ist abgetrennt.

Nach der Filtration

Das Vakuum zwischen Hahn und Membranfilter wird steril und partikelfrei entspannt. Sekundärkontamination der Filterunterseite wird sicher ausgeschlossen.

Nach der Filtrationskampagne

Das Restvakuum zwischen Pumpe und Hahn wird über das Sterilfilter entspannt.

Zum Autoklavieren

Zur sicheren Sterilisation strömt der Dampf ungehindert durch alle Öffnungen.

FR

Pour filtrer

Le vide complet attire l'échantillon à travers la membrane filtrante. L'aération est fermée.

Après la filtration

Le vide entre le robinet et la membrane filtrante est libérée de manière stérile et sans particules. Des contaminations secondaires de la face inférieure du filtre sont exclues de manière certaine.

Après le cycle de filtration

Le vide restant entre la pompe et le robinet est supprimé par l'intermédiaire du filtre stérile.

Pour autoclaver

La vapeur passe librement à travers toutes les ouvertures pour garantir une parfaite stérilisation.

Valve
Position
Open
6 o'clock



Valve Position
Vent | Close
9 o'clock



Valve
Position
Finish
3 o'clock



Valve
Position
Autoclave
12 o'clock



ES

Para la filtración

El vacío total hace pasar la muestra por el filtro de membrana. El filtro del venteo está separado.

Después de la filtración

El vacío entre la llave y el filtro de membrana se expande esterilizado y libre de partículas. La contaminación secundaria de la cara inferior del filtro está descartada.

Después de la sesión de filtración

El vacío restante entre la bomba y la llave se expande sobre el filtro esterilizado.

Para el autoclavado
Para la esterilización segura, el vapor pasa sin dificultades a través de todas las aberturas.

IT

Per la filtrazione

Il sotto vuoto aspira il campione attraverso il filtro a membrana. Il filtro di aerazione è staccato.

Dopo la filtrazione

Il sotto vuoto tra il rubinetto e la membrana viene ridotto in modo sterile e privo di particelle. È esclusa del tutto la contaminazione secondaria del lato inferiore della membrana.

Dopo le operazioni di filtrazione

Il sotto vuoto restante tra la pompa ed il rubinetto viene ridotto attraverso il filtro sterile.

Per la sterilizzazione in autoclave

Per una sicura sterilizzazione il vapore scorre liberamente attraverso tutte le aperture.