

Whatman™ 850-DS 8-Channel Filter Plate

Product Information Sheet

Introduction

Important

Read these instructions carefully before using the products.

Intended use

The products are intended for research use only, and shall not be used in any clinical or *in vitro* procedures for diagnostic purposes.

Safety

For use and handling of the products in a safe way, refer to the Safety Data Sheets.

Description

The Whatman disposable filter plates are designed for use exclusively with an Agilent™ 850-DS Dissolution Sampling Station equipped with the optional filter module.

Load the filter plates in the inlet stack (see Figure 1 below) of the 850-DS and program the method parameters accordingly for filter plate usage.



Fig 2. Inlet face of the filter plates



NOTICE

Do not re-use filters. Blockage could occur which could result in internal leakage.



Fig 1. Inlet stack of 850-DS Dissolution Sampling Station

In order to prevent jamming, make sure the orientation of the filter plate corresponds to the inlet stack and that the plates are seated properly (lie flat, with no gaps between the plates, outer edges aligned) prior to starting the test.

One corner is notched (see Figure 2) to make sure the filter plates are loaded correctly. When loaded correctly, the printing on the filters is facing up. Do NOT force the filters into position – when properly aligned the filters slide gently into position.

Technical data: 850-DS 8-Channel Filter Plate

Weight:	35 g (max)
Filtration area/ filter:	4.2 cm ²
Housing:	Polypropylene
Fluid retention:	Full housing 0.35 ml
With air purge:	<0.2 ml
Flow direction:	Flow should enter from the inlet

Order information

850-DS Filter Plates are available with different membranes and glass fiber media.

Visit [cytiva.com](http://www.cytiva.com) for more information. A useful guide for validating your choice of filters is available from Agilent. This document is available at the following link: http://www.agilent.com/lifesciences/filter_validation.

Product information

Part number	Membrane & pore Size (µm)	Product material
7707-3300	0.7 GMF	Glass microfiber
7707-3900	1.0 GMF	Glass microfiber
7707-3100	0.45 NYLON	Nylon
7707-3500	0.2 NYLON	Nylon
7707-3200	0.45 PES	polyethersulfone

Part number	Membrane & pore Size (µm)	Product material
7707-3600	0.2 PES	polyethersulfone
7707-3000	0.45 PTFE	polytetrafluoroethylene
7707-3400	0.2 PTFE	polytetrafluoroethylene
7707-3700	0.2 PVDF	polyvinylidene difluoride
7707-3800	0.45 PVDF	polyvinylidene difluoride
7707-3700	0.2 PVDF	polyvinylidene difluoride
7707-3800	0.45 PVDF	polyvinylidene difluoride

Dissolution filter selection is dependent on the specific method and the drug formulation under test. The volumetric accuracy of the 850-DS is dependent on the membrane type, the pore size and the cannula filter (Full Flow Filter) being used. The accuracy is also dependent on the drug product and its concentration, the excipient load as well as the media. Other factors that can impact the volume accuracy are the prime volume, pumping speed and dwell time (all of which can be controlled by the 850-DS). Due to high pressures encountered with the use of sub-micron filters, the accuracy specifications for the 850-DS may not be guaranteed for all filter types and drug products.

Chemical resistance summary

The data below is to be used as a guide only.

Testing prior to application is recommended.

Solvent	PVDF	GMF	NYL	PES	PTFE
Acetic acid, 5%	R	R	R	R	R
Acetic acid, glacial	R	L	LR	R	R
Acetone	NR	R	R	NR	R
Acetonitrile	R	R	NR	R	R
Ammonia, 6N	LR	R	R	R	R
Amyl acetate	LR	R	NR	R	R
Amyl alcohol	R	R	R	R	R
Benzene ¹	R	R	LR	R	R
Benzyl alcohol ¹	R	R	LR	NR	R
Boric acid	R	R	LR	-	R
Butyl alcohol	R	R	R	R	R
Butyl chloride ¹	R	R	NR	-	R
Carbon tetrachloride ¹	R	R	LR	NR	R
Chloroform ¹	R	R	NR	NR	R
Chlorobenzene ¹	R	R	NR	NR	R
Citric Acid	R	R	LR	R	R
Cresol ¹	NR	R	NR	NR	R
Cyclohexanone	R	R	NR	NR	R

Solvent	PVDF	GMF	NYL	PES	PTFE
Cyclohexane ¹	R	R	NR	NR	R
Diethyl acetamide	NR	R	R	-	R
Dimethyl formamide	NR	R	R	NR	R
Dioxane	LR	R	R	LR	R
DMSO	LR	R	R	NR	R
Ethanol	R	R	R	R	R
Ethers ¹	LR	R	R	R	R
Ethyl acetate	NR	R	R	NR	R
Ethylene glycol	R	R	R	R	R
Formaldehyde ¹	R	R	R	R	R
Formic acid	R	R	NR	R	R
Freon TF ¹	R	R	NR	R	R
Hexane	R	R	R	R	R
Hydrochloric acid, conc ¹	R	R	NR	R	R
Hydrofluoric acid ¹	R	NR	NR	-	R
Isobutyl alcohol	R	R	R	-	R
Isopropyl alcohol	R	R	R	-	R
Methanol	R	R	R	R	R
Methyl ethyl ketone	NR	R	R	NR	R
Methylene chloride ¹	R	R	NR	NR	R
Nitric acid, conc ¹	R	R	NR	NR	R
Nitric acid, 6N ¹	R	R	NR	LR	R
Nitrobenzene ¹	R	R	LR	NR	R
Pentane ¹	R	R	R	R	R
Perchloro ethylene ¹	R	R	LR	NR	R
Phenol 0.5%	R	R	NR	NR	R
Pyridine	NR	R	LR	NR	R
Sodium hydroxide, 6N	NR	NR	LR	R	R
Sulfuric acid, conc ¹	NR	R	NR	NR	R
Tetrahydrofuran	R	R	R	NR	R
Toluene ¹	R	R	LR	NR	R
Trichloroethane ¹	R	R	LR	NR	R
Trichloroethylene ¹	R	R	NR	NR	R
Water	R	R	R	R	R
Xylene ¹	R	R	LR	LR	R

¹ Short term resistance of housing.

R = Resistant; LR = Limited resistant; NR = Not resistant.

Technical support

For questions related to the use of these filter plates contact Cytiva. Cytiva are happy to assist in resolving problems you may encounter with the 850-DS, the filter changer or the filter plates. To contact the Cytiva technical support group, find the appropriate telephone number for your country at cytiva.com.

cytiva.com

Cytiva and the Drop Logo are trademarks of Global Life Sciences IP Holdco LLC or an affiliate.

Whatman is a trademark of Global Life Sciences Solutions USA LLC or an affiliate doing business as Cytiva.

Agilent is a trademark of Agilent Technologies, Inc.

All other third-party trademarks are the property of their respective owners.

© 2020–2022 Cytiva

For local office contact information, visit cytiva.com/contact

29086372 AI V:6 02/2022

