



Revised: September 19, 2018

# **Product Information**

# Mini Syringe Filters and Mini Syringe Filtration Kit

# Contents

Component	22025 Set of 5 filters	22026 Kit for 5 samples
Millex®-GV 4 mm, 0.22 um filter (individually wrapped)	5 filters	5 filters
Sterile 1 cc syringe (individually wrapped)	Not included	5 syringes
Sterile 2 mL self-standing collection vial with O-ring screw cap	Not included	5 vials

# Storage and Handling

Store at room temperature.

#### **Product Description**

Mini Syringe Filters are 4 mm diameter, 0.22 um pore size sterile filters for filter sterilization of 1 mL or less of aqueous solution. With only 10 uL hold-up volume, they can be used to filter as little as 100 uL, and are suitable for filtering solutions of our Azide-Free CF® Dye Annexin V Conjugates for live cell imaging.

#### Filter specifications

- Millex®-GV hydrophilic 0.22 um pore size PVDF membrane
- · Ultra-low protein binding
- 4 mm diameter, 19.7 mm height
- · Female Luer-Lok® inlet
- Male Luer slip outlet, stepped
- Maximum inlet pressure: 14 bar (200 psi)
- Filter 100 uL to 1 mL with 10 uL hold-up volume
- Ethylene oxide sterilized

# Instructions

Caution: When using syringe filters, we recommend always wearing eye protection or a face shield, or performing filtration behind a hood screen. Applying excessive force to syringe filters can lead to filter failure, causing liquid to spray out.

- Peel back the end of a 1 cc syringe wrapper. Holding the syringe through the wrapper, remove the plunger and place it on a clean surface. Leave the barrel inside the wrapper for now.
- Place the sterile collection vial in a rack. Remove the cap and place it on a clean surface with the open end up.
- Peel back the wrapper on the filter blister pack from the side near the wide (female) end of the filter.
- Remove the syringe barrel from its wrapper, taking care to touch only the wide end near the top.
- While grasping the filter through its wrapper, insert the syringe tip into the female end of the filter and firmly attach with a twist.
- Remove the filter blister pack, taking care not to touch the filter or lower end of the syringe barrel.
- Place the filter and syringe inside the receiving vial so that the syringe stands up
- 8. Carefully pipette the solution into the open barrel of the syringe.
- Replace the plunger in the syringe, and slowly filter the antibody in a dropwise fashion into the collection vial. Do not use maximum force to push the plunger when using a 1 cc syringe, which this will exceed the maximum filter input pressure and may cause the filter to fail.
- 10. Discard the syringe and filter and cap the vial. Do not reuse the filters.

#### Chemical compatibility

The Millex®-GV PVDF membrane is intended for filtration of aqueous solutions. The filter manufacturer reports that the membrane is compatible with the following chemicals. It is NOT recommended for use with dimethylsulfoxide (DMSO) or dimethylformamide (DMF).

Acetic acid (glacial) Acetic acid (5%)	Carbon tetrachloride Ethanol (ethyl alcohol)	Isopropyl acetate Methyl alcohol
Aliphatic ethers	Ethylene glycol	Mineral spirits
Ammonium hydroxide (6N)	Formaldehyde	Paraldehyde
Amyl acetate	Freon® solvent, TF or PCA	Pentane
Amyl alcohol	Glycerol (glycerine)	Pet base oils
Benzyl alcohol (1%)	Hexane	Petroleum ether
Boric acid	Hydroflouric acid	Phenol (0.5%)
Brine (sea water)	Hydrogen peroxide (3%)	Pyridine
Butyl alcohol	Isobutyl alcohol	

#### Related Products

Catalog number	Product	
29009R-5ug	Annexin V, CF®405M Conjugate, Azide-Free and Lyophilized	
29005R-5ug	Annexin V, CF®488A Conjugate, Azide-Free and Lyophilized	
29010R-5ug	Annexin V, CF®568 Conjugate, Azide-Free and Lyophilized	
29011R-5ug	Annexin V, CF®594 Conjugate, Azide-Free and Lyophilized	
29014R-5ug	Annexin V, CF®640R Conjugate, Azide-Free and Lyophilized	
29007	Annexin V, CF®680 Conjugate, Azide-Free and Lyophilized	
29070	Annexin V, CF®680R Conjugate, Azide-Free and Lyophilized	
29082	Annexin V, CF®700 Conjugate, Azide-Free and Lyophilized	
29006	Annexin V, CF®750 Conjugate, Azide-Free and Lyophilized	
29046	Annexin V, CF®770 Conjugate, Azide-Free and Lyophilized	
29047	Annexin V, CF®790 Conjugate, Azide-Free and Lyophilized	
29078	Annexin V, CF®800 Conjugate, Azide-Free and Lyophilized	
10403	NucView® 488 Caspase-3 Substrate for live cells, 1 mM in PBS	
10408	NucView® 530 Caspase-3 Substrate for live cells, 1 mM in PBS	
10407	NucView® 405 Caspase-3 Substrate for live cells, 1 mM in PBS	
40061	RedDot™2 Far-Red Nuclear Stain (for dead or fixed cells)	
40083	NucSpot® 470 Green Nuclear Stain (for dead or fixed cells)	
40017	Propidium lodide, 1 mg/mL in water	
40084	7-AAD Solution, 1 mg/mL	
40050	Ethidium Homodimer III	
10126	Aquaphile™ Coelenterazine Native, lyophilized solid	
10127	Aquaphile™ Coelenterazine h, lyophilized solid	
22003	Mini-Cell Scrapers	
22023	Paraformaldehyde, 4% in PBS, Ready-to-Use Fixative	
22020	10X Phosphate-Buffered Saline (PBS)	

Please visit our website at <a href="www.biotium.com">www.biotium.com</a> for information on our life science research products, including fluorescent organelle stains for live cells, Azide-Free Annexin V, dead cell stains, NucView® Caspase Substrates, Mix-n-Stain™ CF® Dye Antibody Labeling Kits, and more.

Freon is a registered trademark of The Chemours Company. Millex is a registered trademark of Merck KGaA. Luer-Lok is a registered trademark of Becton Dickinson.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.