



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Mini Syringe Filters and Mini Syringe Filtration Kit

These sterile 0.22 μ m filters are useful for filtering 100 μ L to 1 mL of aqueous solutions with only 10 μ L hold-up volume.



Product attributes

Storage Conditions

Store at room temperature

Product Description

Mini Syringe Filters are 0.22 μ m sterile filters for filter sterilization of 100 μ L to 1 mL of aqueous solution with only 10 μ L hold-up volume. For convenience, we offer these filters in a pack of 5 or in a kit for filtering 5 samples.

Mini Syringe Filters are ideal for filtering solutions of our [Azide-Free CF® Dye Annexin V Conjugates](#) and other solutions for live cell imaging. These filters are only suitable for aqueous solutions, and not compatible with DMSO (see the product information sheet for more information on chemical compatibility).

Millex is a registered trademark of Merck KGaA. Luer-Lok is a registered trademark of Becton Dickinson

This datasheet was generated on January 8, 2026 at 09:51:58 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/mini-syringe-filters-and-filtration-kit/>