



Catalog #	Description
12012577	ProFlow Cell Filter, nonsterile, 30 µm, disposable, cup fitting, 100 pieces
12012582	ProFlow Cell Filter, nonsterile, 50 µm, disposable, cup fitting, 100 pieces
12012575	ProFlow Cell Filter, nonsterile, 70 µm, disposable cup fitting, 100 pieces
12012576	ProFlow Cell Filter, sterile, 30 µm, disposable cup fitting, 100 pieces
12012574	ProFlow Cell Filter, sterile, 50 µm, disposable cup fitting, 100 pieces
12012573	ProFlow Cell Filter, sterile, 70 µm, disposable cup fitting, 100 pieces

For research use only.

Intended Use

ProFlow Cell Filters are disposable cup-fitting filters used to remove aggregates. Researchers needing uniform single-cell suspensions can use ProFlow Cell Filters to rapidly prepare sample suspensions from tissue, fluids, plant material, nuclei, etc.

Using the ProFlow Cell Filters

Each ProFlow Cell Filter has a porous membrane with effective surface area of 80 mm², enabling large quantities of cell suspensions to be filtered. The cup-fitting ProFlow Filters fit 5 ml tubes (12 x 75 mm) and are ideal for quantities of less than 5 ml. The microporous polyethylene funnel section is designed to pass the biological material to be retrieved while helping remove debris by trapping waste particles less than 1.5 µm in size.

Note: Sterile ProFlow Cell Filters are EtO sterilized. Use them when sterility is needed. ProFlow Cell Filters should not be sterilized by the user as that may damage them.

Do not use ProFlow Filters more than once, as subsequent specimens could be contaminated.

1. For best results, wet the filter membrane with a drop of the cell or nuclei suspension medium. Sticky cell types may adhere to the filter membrane if it is left dry.
2. Avoid touching the membrane, allow only the liquid to come into contact with the membrane.
3. Ensure used ProFlow Filters are properly disposed of. Micro-residues of biological materials remain in the plastics and membrane.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc.

All trademarks used herein are the property of their respective owner.