



Glowing products for science

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

Ultrafiltration Vials

Ultrafiltration devices for simple spin-column purification of proteins or other molecules by size separation with two molecular weight cut-off options: MWCO=10 kD (10K) and MWCO=3 kD (3K).



Product attributes

MWCO	10K, 3K
------	---------

Product Description

These centrifugal ultrafiltration vials are designed for simple spin-column purification of proteins or other molecules by size separation. There are two molecular weight cut-off options: MWCO=10 kD (10K) and MWCO=3 kD (3K).

- Concentrate proteins or other molecules
- Remove small molecules (ie, Tris, glycine, glycerol)
- Remove free dye or biotin from antibodies after conjugation
- 10 kD or 3 kD Molecular Weight cut-offs
- 500 uL maximum sample volume
- Typically 80-90% protein recovery
- Pack of 5 columns

The ultrafiltration membrane allows molecules with molecular weight smaller than the cut-off to pass through into the filtrate collection tube, while molecules larger than the cut-off are retained on top of the membrane.

Ultrafiltration can be used to rapidly and easily remove small molecules such as Tris, glycerol and glycine from antibody solutions, or to concentrate antibodies prior to conjugation. They also can be used to remove free dye or free biotin from antibodies after conjugation.

For ultrafiltration of proteins, it is recommended to use a MWCO 3- to 6-times smaller than the protein to be retained. Ultrafiltration vials with MWCO=10K are recommended for ultrafiltration of antibodies and other proteins larger than 30 kDa. Vials with MWCO=3K are recommended for ultrafiltration of proteins larger than 10 kDa.

Ultrafiltration is not effective for fractionating proteins (separating one protein from another protein based on size) unless there is at least a 10-fold difference in the protein molecular weights.

This datasheet was generated on January 13, 2026 at 04:07:14 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/ultrafiltration-vial/>