

ReadiUse™ 10KD Spin Filter

 Catalog number: 60502
 Unit size: 5

Component	Storage	Amount
ReadiUse™ 10KD Spin Filter	Room temperature (10-25 °C)	5 units

OVERVIEW

ReadiUse™ 10KD Spin Filter is a disposable ultrafiltration device for the concentration of biological samples such as antibodies. Maximum initial sample volume is ~500 µL. It can be used in either a swing bucket or fixed angle rotors accepting 2.0 mL centrifuge tube at around 10000 g. The dead stop volume is ~20-25 µL. The vertical membrane design and thin channel filtration chamber minimizes membrane fouling and provides high speed concentration, even laden solutions. It is specifically designed with low internal surface and membrane area to achieve superior recovery from dilute solutions.

AT A GLANCE
Overview

1. Polypropylene housing and polyethersulfone (PES) membrane
2. MWCO: 10K
3. Sample Volume: 25 µL-500 µL
4. High sample recovery
5. Application: Protein concentration, desalting, buffer exchange, and deproteinization of biological samples
6. Maximum Centrifugal Force: 14,000 g
7. Recommended Centrifugal Force: 5,000 - 10,000 g

SAMPLE EXPERIMENTAL PROTOCOL
Concentrate volume

Add protein sample to the filter, centrifuge at 8,000 - 10,000 g.

Note it takes about 10 - 15 mins to concentrate from 500 µL to ~50 µL.

Desalting

Desalting, buffer exchange or diafiltration are important methods for removing salts or solvents in solutions containing biomolecules. The removal of salts or the exchange of buffers can be accomplished in the ReadiUse™ 10KD Spin Filter by concentrating the sample, then reconstituting the concentrate to the original sample volume with any desired solvent. The process of "washing out" can be repeated until the concentration of the contaminating microsolute has been sufficiently reduced.

Deproteinization of biological samples

Add cell lysates or tissue homogenates to the filter, centrifuge at ~10,000 g for 10 - 15min, collect the elution.

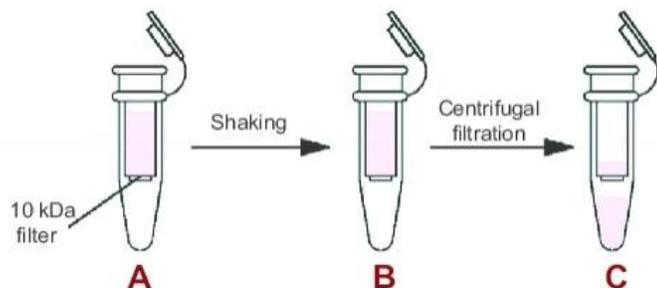
EXAMPLE DATA ANALYSIS AND FIGURES


Figure 1. ReadiUse™ 10KD Spin Filters are disposable ultrafiltration devices for the concentration of biological samples such as antibodies. Maximum initial sample volume is ~500 µL. They can be used in either a swing bucket or fixed angle rotors accepting 2.0 ml centrifuge tube at around 10000 g.

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.