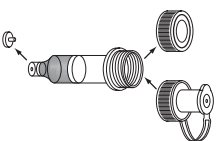
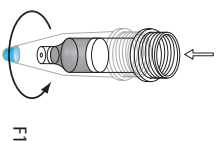


1 Dilute and filter sample

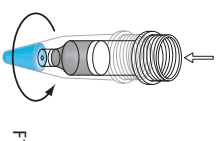
Dilute 14–16-µL human serum sample to 200 µL with Buffer A. Consult cartridge certificate for true sample capacity. Filter through 0.22-µm spin filter.

2 Prepare spin cartridge

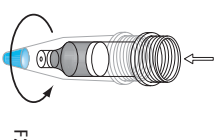
Remove cartridge cap and plug, attach Luer-lock adapter to cartridge, draw 4 mL of Buffer A into syringe and dispense through cartridge via Luer-Lock, remove excess Buffer A from top of resin bed with transfer pipette.

3 Apply sample

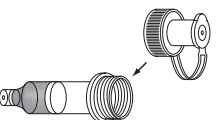
Remove Luer-Lock adapter and add 200-µL diluted serum sample. Cap cartridge loosely or leave open. Place in 1.5-mL collection tube labeled "Flow-through fraction 1" (F1). Centrifuge 1.5 min at 100 × g.

4 Wash and collect flow-through F1

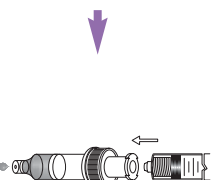
Add 400-µL Buffer A. Centrifuge 2.5 min at 100 × g. Collect in F1 tube.

5 Wash and collect flow-through F2

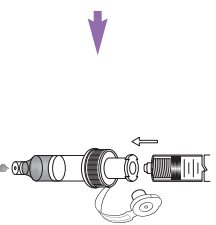
Place spin cartridge in new collection tube labeled "Flow-through fraction 2" (F2). Add 400-µL Buffer A. Centrifuge 2.5 min at 100 × g. Collect in F2 tube.

6 Prepare for elution

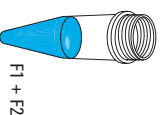
Remove spin cartridge from F2 tube and attach Luer-Lock adapter tightly to top of cartridge.

7 Elute bound fraction

Fill 5-mL Luer-Lock plastic syringe with 2-mL Buffer B and attach to Luer-Lock adapter. Slowly push Buffer B through cartridge to elute bound proteins into new collection tube. Save eluent with targeted high-abundant proteins for analysis or discard.

8 Re-Equilibration

Fill new 5-mL plastic syringe with 4-mL Buffer A and attach to Luer-Lock adapter. Slowly push Buffer A through cartridge to re-equilibrate the cartridge for the next sample or store wetted with Buffer A (at 4 °C). Recap both ends for storage.

9 Analyze F1 + F2

Fractions F1 and F2 can be analyzed individually or combined. Concentrate and analyze these fractions containing low-abundant proteins.



Agilent Technologies

For more detailed instructions or information on accessories refer to the Agilent High Capacity Multiple Affinity Removal Spin Cartridge Instruction Guide

Visit www.agilent.com/chem/bioreagents for more products that will help make your proteomics research more efficient:

- Multiple Affinity Removal LC columns for automated serum/plasma processing
- Lys Tag 4H reagent for improving MALDI-MS analysis of peptides and proteins
- Tryptic digestion kit for digesting gel-isolated proteins
- C18 Cleanup Pipette Tips and Spin Tubes for purifying proteomics samples
- Spin filters and concentrators for cleaning, concentrating, and buffer-exchanging samples
- MALDI-MS matrices and peptide/protein standards for your proteomics studies
- mRP-C18 column for extremely high recovery protein fractionation or desalting



For More Information

For more information on our products and services, visit our Web site at www.agilent.com/chem/bioreagents.

Or call your local authorized Agilent distributor.

Agilent shall not be liable for errors contained herein or for incidental or consequential damages in connection with the furnishing, performance, or use of this material.

Information, descriptions, and specifications in this publication are subject to change without notice.

© Agilent Technologies, Inc. 2005

Printed in the USA
March 29, 2005
5188-5931EN



For Human Serum

Agilent High Capacity Multiple Affinity Removal Spin Cartridge

Part Number 5188-5341

Quick Reference Guide

