

ProteoMiner™ Sequential Elution Large-Capacity Kit Protocol



Downstream Application Compatibility

The ProteoMiner sequential elution kit is available for the ProteinChip® surface enhanced laser desorption/ionization (SELDI) system or other downstream protein separation analysis methods.

This kit is NOT compatible with 2-D gel electrophoresis.

Sample Preparation

This kit is optimized for plasma and serum samples; best results are obtained with ≥ 50 mg protein load. Samples should be free of precipitate. If needed, centrifuge the samples at 10,000 x g for 10 min to clarify.

Perform the following steps to complete the protocol:

Procedure	Add*	Rotate/ Vortex	Centrifuge** 1,000 x g	Repeat Step
Remove top and bottom caps. Centrifuge column for 30–60 sec to remove storage material. Discard flowthrough.	—	—	30–60 sec	—
Add 600 μ l wash buffer to column. Rotate column for 5 min. Centrifuge for 30–60 sec. Discard flowthrough and repeat entire step.	600 μ l wash buffer	5 min	30–60 sec	x1
Add 1 ml sample to column. Rotate for 2 hr at room temperature. Centrifuge for 30–60 sec. Discard flowthrough and repeat centrifugation.	1 ml sample	2 hr	30–60 sec	—
Add 600 μ l wash buffer to column. Rotate column for 5 min. Centrifuge for 30–60 sec. Discard flowthrough and repeat entire step three times.	600 μ l wash buffer	5 min	30–60 sec	x3
Add 600 μ l DI water on all sides of the column. Centrifuge for 30–60 sec. Discard flowthrough.	600 μ l DI water	—	30–60 sec	—
Add 100 μ l elution reagent 1 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	100 μ l elution reagent 1	10 min	30–60 sec	x2
Add 100 μ l elution reagent 2 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	100 μ l elution reagent 2	10 min	30–60 sec	x2
Add 100 μ l elution reagent 3 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	100 μ l elution reagent 3	10 min	30–60 sec	x2
Add 100 μ l elution reagent 4 to column. Vortex gently for 5 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	100 μ l elution reagent 4	5 min	30–60 sec	x2

* Remove top cap and assure bottom cap is secure on column prior to each addition step. Replace top cap prior to rotating (inverting).

** Remove top and bottom caps before centrifugation. Perform centrifugation steps at 1,000 x g. Refer to manual for detailed protocol.

*** Eluates may be pooled or collected individually; if collected individually, additional collection tubes will be required.

BIO-RAD

ProteoMiner™ Sequential Elution Small-Capacity Kit Protocol



Downstream Application Compatibility

The ProteoMiner sequential elution kit is available for the ProteinChip® surface enhanced laser desorption/ionization (SELDI) system or other downstream protein separation analysis methods.

This kit is NOT compatible with 2-D gel electrophoresis.

Sample Preparation

This kit is optimized for plasma and serum samples; best results are obtained with ≥ 10 mg protein load. Samples should be free of precipitate. If needed, centrifuge the samples at 10,000 x g for 10 min to clarify.

Perform the following steps to complete the protocol:

Procedure	Add*	Rotate/ Vortex	Centrifuge** 1,000 x g	Repeat Step
Remove top and bottom caps. Centrifuge column for 30–60 sec to remove storage material. Discard flowthrough.	—	—	30–60 sec	—
Add 200 μ l wash buffer to column. Rotate column for 5 min. Centrifuge for 30–60 sec. Discard flowthrough and repeat entire step.	200 μ l wash buffer	5 min	30–60 sec	x1
Add 200 μ l sample to column. Rotate for 2 hr at room temperature. Centrifuge for 30–60 sec. Discard flowthrough.	200 μ l sample	2 hr	30–60 sec	—
Add 200 μ l wash buffer to column. Rotate column for 5 min. Centrifuge for 30–60 sec. Discard flowthrough. Repeat entire step two times	200 μ l wash buffer	5 min	30–60 sec	x2
Add 200 μ l DI water on all sides of the column. Centrifuge for 30–60 sec. Discard flowthrough.	200 μ l wash buffer	—	30–60 sec	—
Add 20 μ l elution reagent 1 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	20 μ l elution reagent 1	10 min	30–60 sec	x2
Add 20 μ l elution reagent 2 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	20 μ l elution reagent 2	10 min	30–60 sec	x2
Add 20 μ l elution reagent 3 to column. Vortex gently for 10 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	20 μ l elution reagent 3	10 min	30–60 sec	x2
Add 20 μ l elution reagent 4 to column. Vortex gently for 5 min. Place into a clean tube and centrifuge for 30–60 sec. Repeat entire step twice.***	20 μ l elution reagent 4	5 min	30–60 sec	x2

* Remove top cap and assure bottom cap is secure on column prior to each addition step. Replace top cap prior to rotating (inverting).

** Remove top and bottom caps before centrifugation. Perform centrifugation steps at 1,000 x g. Refer to manual for detailed protocol.

*** Eluates may be pooled or collected individually; if collected individually, additional collection tubes will be required.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 4BIORAD **Australia** 61 02 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400
Canada 905 364 3435 **China** 86 21 6426 0808 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65
Germany 089 318 84 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 455 8800 **India** 91 124 4029300 **Israel** 03 963 6050
Italy 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666 **New Zealand** 0508 805 500
Norway 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04 **Singapore** 65 6415 3188 **South Africa** 27 861 246 723
Spain 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55 **Taiwan** 886 2 2578 7189 **United Kingdom** 020 8328 2000