

## Example Protocol: Antibody–Antigen Capture Using Protein G

**Note:** This protocol is intended as a guideline; you may need to optimize the method for your particular sample type and experimental design.

1. Apply 2  $\mu$ l of 0.5 mg/ml protein G and 2  $\mu$ l 50 mM NaHCO<sub>3</sub> buffer, pH 9.2 to spot. Incubate the array for 1 hour at RT in a humid chamber.
2. Remove protein G and wash with 5  $\mu$ l 1x PBS for 2 minutes on shaker.
3. Remove buffer and apply 2  $\mu$ l 100  $\mu$ M BSA and 2  $\mu$ l of 1x PBS, pH 7.2 for 1 hour at RT to block any remaining active sites.
4. Remove BSA and add 5  $\mu$ l of 1x PBS. Shake for 2 minutes on a shaker and then remove buffer from spot. Repeat wash step.
5. Remove the buffer and apply 2  $\mu$ l of 0.1 mg/ml antibody and 2  $\mu$ l 50 mM NaHCO<sub>3</sub> buffer, pH 9.2 to spot. Incubate the array for 2 hours at RT in a humid chamber.
6. Remove antibody and wash with 5  $\mu$ l 1x PBS for 2 minutes on shaker.
7. Remove buffer. Prepare the antigen in 50% human serum diluted in a 0.1% Triton X-100 in 1x PBS solution, pH 7.2. Apply 2  $\mu$ l of antigen and incubate for 2 hours at RT.
8. Remove antigen and add 5  $\mu$ l of 1x PBS and wash for 30 seconds on shaker.
9. Remove buffer and add 5  $\mu$ l urea CHAPS buffer (50 mM Tris + 1 M urea + 0.1% CHAPS + 0.5 M NaCl, pH 7.2); wash buffer for 30 seconds on shaker.
10. Remove buffer and add 5  $\mu$ l of 1x PBS, pH 7.2; wash for 2 minutes on shaker.
11. Quickly bulk-rinse the array with 5 mM HEPES, pH 7.2 in a 15 ml tube for 10 seconds.
12. Allow the array surface to air-dry (5 minutes).
13. Add 1  $\mu$ l of 50% saturated SPA EAM solution to each spot. Allow to air-dry. Repeat addition of SPA solution and allow to dry.
14. Analyze the array using the ProteinChip SELDI system.

## Ordering Information

Catalog #	Description
C55-30082	<b>ProteinChip RS100 Arrays</b> , A–H format, 12
C30-00002	<b>ProteinChip SPA Energy Absorbing Molecules (EAMs)</b> , 5 mg/vial, 20

Triton is a trademark of Union Carbide.

The SELDI process is covered by US patents 5,719,060, 5,894,063, 6,020,208, 6,027,942, 6,124,137, 6,225,047, 6,528,320, 6,579,719, and 6,734,022. Additional US and foreign patents are pending.

# ProteinChip® RS100 Array (Reactive Surface)

## Instruction Manual

Catalog #C55-30082

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://www.bio-rad.com) **USA** 800 4BIORAD  
**Australia** 02 9914 2800 **Austria** 01 877 89 01  
**Belgium** 09 385 55 11 **Brazil** 55 21 3237 9400  
**Canada** 905 712 2771 **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 4029300/5013478 **Israel** 03 963 6050  
**Italy** 39 02 216091 **Japan** 03 5811 6270 **Korea** 82 2 3473 4460  
**Mexico** 55 5200 05 20 **The Netherlands** 0318 540666  
**New Zealand** 64 9415 2280 **Norway** 23 38 41 30  
**Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700  
**Russia** 7 095 721 14 04 **Singapore** 65 6415 3188  
**South Africa** 27 0861 246 723 **Spain** 34 91 590 5200  
**Sweden** 08 555 12700 **Switzerland** 061 717 95 55  
**Taiwan** 886 2 2578 7189/2578 7241  
**United Kingdom** 020 8328 2000

For technical support,  
call your local Bio-Rad office, or  
in the US, call **1-800-4BIORAD**  
**(1-800-424-6723)**.

**BIO-RAD**

## Uses

- Covalent immobilization of biomolecules for the subsequent capture of proteins from complex biological samples

## How It Works

The ProteinChip RS100 array is preactivated with carbonyl diimidazole chemistry that reacts with amine groups on the surface of biomolecules to form stable covalent linkages. In turn, these immobilized biomolecules capture proteins from biological samples through specific, noncovalent interactions.

## Packaging and Storage

ProteinChip RS100 arrays in sealed tubes are vacuum-packaged and protected from moisture. Store the arrays between  $-20^{\circ}\text{C}$  and  $-50^{\circ}\text{C}$  in the original packaging. Open the array tube just prior to use. Once opened, the arrays have limited stability.

## Technical Considerations

- Optimization of assay parameters will be required for each specific application
- Proteins primarily couple through amine groups
- For coupling, use pure biomolecules without carrier proteins whenever possible
- Generally, couple biomolecules at pH 7.2–9.2, using phosphate buffered saline (PBS) or sodium bicarbonate buffers
- Caution: During coupling, avoid buffers containing free amines (e.g., glycine, Tris), free thiols (e.g., dithiothreitol,  $\beta$ -mercaptoethanol), and azide
- Avoid physical contact with spot surface and surrounding coating
- Array design allows sample containment of up to 5  $\mu\text{l}$  for on-spot application and up to 180  $\mu\text{l}$  for bioprocessor application
- Include denaturants, salts, and chaotropic agents in binding and wash buffers, as required to modify binding stringency and reduce nonspecific binding
- Prepare energy absorbing molecule (EAM) solution according to instructions in EAM product insert

## Recommended Buffers

### Coupling Buffers

- PBS or sodium bicarbonate, pH 7.2–9.2
- Avoid buffers containing free amines, free thiols, or azide

### Blocking Buffers

- Ethanolamine (0.5 M), pH 8
- Tris-HCl or glycine (0.1–0.5 M), pH 8
- Bovine serum albumin (BSA) (100  $\mu\text{M}$  in 1x PBS), pH 7.2

### Washing Buffers

- Post-blocking, use buffers and additives more stringent than binding conditions. Repeated washing and pH cycling may be necessary
- Post-binding, use buffers containing 0–2.0 M salt, 0–2.0 M denaturant (urea), and 0.1–0.5% detergent (Triton X-100), prepared in PBS, pH 7–7.5 or 50 mM Tris
- Include nonionic detergent, salts, and chaotropic agents as required
- A final rinse (10–20 seconds) with 5 mM HEPES is often required to remove salts and detergents from washing buffers

### Binding Buffers

- PBS, pH 7–7.5 or buffer of choice
- Include nonionic detergent (e.g., 0.1–0.5% Triton X-100 or CHAPS) as needed
- Include salt (0.15–1.0 M), other modifiers (e.g., ethylene glycol), and carrier protein (e.g., 1% BSA), if necessary

## Example Protocol: Antibody–Antigen Capture Assay

**Note:** This protocol is intended as a guideline; you may need to optimize the method for your particular sample type and experimental design.

1. Apply 2  $\mu\text{l}$  of 0.25 mg/ml antibody and 2  $\mu\text{l}$  50 mM  $\text{NaHCO}_3$  buffer, pH 9.2.
2. Incubate the array for 1 hour at room temperature (RT) in a humid chamber.
3. Remove antibody and apply 2  $\mu\text{l}$  100  $\mu\text{M}$  BSA and 2  $\mu\text{l}$  of 1x PBS, pH 7.2 for 1 hour at RT to block any remaining active sites.
4. Remove the blocking solution and add 5  $\mu\text{l}$  of 1x PBS. Shake for 2 minutes on a shaker and then remove buffer from spots. Repeat wash step.
5. Prepare the antigen in 50% human serum diluted in 0.1% Triton X-100 in 1x PBS solution, pH 7.2. Apply 2  $\mu\text{l}$  of antigen and incubate for 2 hours at RT.
6. Remove antigen and add 5  $\mu\text{l}$  of 1x PBS and wash for 30 seconds on a shaker.
7. Remove buffer and add 5  $\mu\text{l}$  urea CHAPS buffer (50 mM Tris + 1 M urea + 0.1% CHAPS + 0.5 M NaCl, pH 7.2); wash buffer for 30 seconds on shaker.
8. Remove buffer and add 5  $\mu\text{l}$  of 1x PBS, pH 7.2; wash for 2 minutes on shaker.
9. Quickly bulk-rinse the array with 5 mM HEPES, pH 7.2 in a 15 ml tube for 10 seconds.
10. Allow the array surface to air-dry (5 minutes).
11. Add 1  $\mu\text{l}$  of 50% saturated sinapinic acid (SPA) EAM (catalog #C30-00002) solution to each spot. Allow to air-dry. Repeat addition of SPA solution and allow to dry.
12. Analyze the array using the ProteinChip SELDI system.