

The SELDI process is covered by US patents 5,719,060, 6,225,047, 6,579,719, 6,818,411, and other issued patents and pending applications in the US and other jurisdictions.

ProteinChip® Gold Array

Instruction Manual

Catalog #C55-30033

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site 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

Application

The ProteinChip gold array can be used to conduct conventional matrix-assisted laser desorption/ionization (MALDI) experiments with the ProteinChip SELDI system.

Principle

The gold surface provides a nonreactive surface that can be used to evaluate purified proteins.

Storage

Store the array at room temperature.

Sample Preparation

1. Mix analyte with energy absorbing molecule (EAM) solution.
2. Apply the solution to the ProteinChip gold array and air-dry for approximately 10 minutes.

ProteinChip Gold Array Preparation

1. Wash the ProteinChip array with acetone before use.
2. Rock-wash (bulk-wash) the array in a 0.01N HCl solution for five minutes. Discard the used wash solution.
3. Rock-wash (bulk-wash) in a 50:50 mixture of 0.01N HCl/methanol (HPLC grade) for five minutes. Discard solution.
4. Rock-wash (bulk-wash) in methanol (HPLC grade) for five minutes. Discard the used wash solution and air-dry the array for 10 minutes.

ProteinChip gold arrays can be cleaned and reused; however, performance may deteriorate with use.

Ordering Information

Catalog #	Description
C55-30033	ProteinChip Gold Array , A–H format
C30-00001	ProteinChip CHCA Energy Absorbing Molecules (EAMs) , 5 mg/vial, 20
C30-00002	ProteinChip SPA Energy Absorbing Molecules (EAMs) , 5 mg/vial, 20
C30-00003	ProteinChip EAM-1 Energy Absorbing Molecules (EAMs) , 5 mg/vial, 20