

Ordering Information

Catalog #	Description
K20-00004	ProteinChip CM High-Stringency Buffer , 200 ml
K20-00003	ProteinChip CM Low-Stringency Buffer , 200 ml
C57-30075	ProteinChip CM10 Arrays , A-H format, 12
C50-30011	ProteinChip Cassette-Compatible Bioprocessor , includes ProteinChip array forceps, cassette hold-down frame, 12 blank ProteinChip arrays
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

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



**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 905 712 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
Israel 03 963 6050 **Italy** 39 02 216091
Japan 03 5811 6270 **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

ProteinChip® CM High-Stringency Buffer

Instruction Manual

Catalog #K20-00004

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

Uses

- Protein profiling and biomarker discovery
- Rapid protein analysis to determine purity, mass confirmation, or both

Introduction

ProteinChip CM high-stringency buffer is designed for use with the ProteinChip CM10 array. The ProteinChip CM10 array incorporates a carboxylate chemistry (negatively charged) that acts as a weak cation exchanger. The carboxymethyl (CM) surface binds proteins that are positively charged at a given pH. To control selectivity, the pH of the binding buffer is increased or decreased, depending on the need. By using the higher pH ProteinChip CM high-stringency buffer, an overall net negative charge is imparted on some of the proteins, resulting in fewer proteins binding, i.e., higher stringency. Using a low-stringency buffer imparts an overall net positive charge on a greater number of proteins within the sample, resulting in more proteins binding, i.e., lower stringency.

Storage

Store buffer at 2–8°C.

Buffer Composition

50 mM HEPES, pH 7.0, antimicrobial preservatives, 200 ml — Amount is sufficient to run 12 ProteinChip arrays in a ProteinChip cassette-compatible bioprocessor (catalog #C50-30011), using binding and elution buffer volumes as outlined in the suggested protocol below.

Suggested Protocol

1. Assemble ProteinChip arrays in the ProteinChip bioprocessor.
2. Add 150 µl of the ProteinChip CM high-stringency buffer to each well. Shake for 5 minutes at room temperature.
3. Remove buffer from wells.
4. Repeat steps 2–3 for a total of two washes.
5. Add 90 µl of ProteinChip CM high-stringency buffer to each well.

6. Add 10 µl of sample to each well. Shake for 30 minutes at room temperature.
7. Remove samples from wells.
8. Wash each well with 150 µl of ProteinChip CM high-stringency buffer for 5 minutes, with agitation. Repeat twice for a total of three buffer washes.
9. Remove wash buffer from wells and rinse each well with deionized water.
10. Drain wells and remove arrays from the ProteinChip bioprocessor.
11. Allow arrays to air-dry.
12. Apply 1.0 µl of energy absorbing molecule (EAM) solution per spot. Two applications of EAM solution can be used in order to increase the intensities of high-mass proteins. Allow arrays to air-dry.
13. Analyze arrays using the ProteinChip SELDI reader.