



SELDI TECHNOLOGY

ProteinChip® System Starter Kit

- Video tutorial
- Software basics
- External mass calibration
- Skill level testing

Hit the Ground Running With the ProteinChip System Starter Kit

Introduction

The new ProteinChip system starter kit has been redesigned to deliver a step-by-step focused introduction to ProteinChip SELDI technology. The kit is a great training tool that contains everything needed to generate protein profiles from serum and *E. coli* samples. Using the included analysis tools, determine the reproducibility and optimal peak count of your protein profiles. Learn what to expect in terms of reproducibility by comparing your results to results from the included premade ProteinChip arrays. Improve your results by employing valuable tips and techniques outlined in the manual.

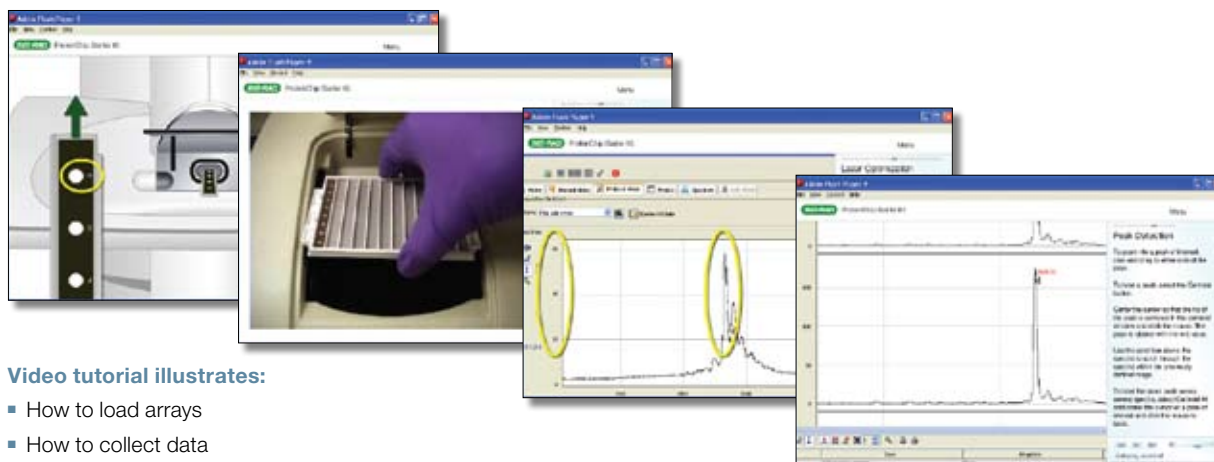
Learn the Basics Fast With a Video Tutorial

The video tutorial included with the starter kit illustrates the basic technology of the ProteinChip SELDI system. It starts with a visual description of the anatomy of a ProteinChip array and ends with instructions for performing mass calibrations. Watch the video to become familiar with the concepts, then perform the same tasks using the ProteinChip peptide standard array and instructions included with the kit. The video is also available online at www.bio-rad.com/proteinchip/starterkit/.



BIO-RAD

Everything You Need to Get Started



Video tutorial illustrates:

- How to load arrays
- How to collect data
- How to perform mass calibrations

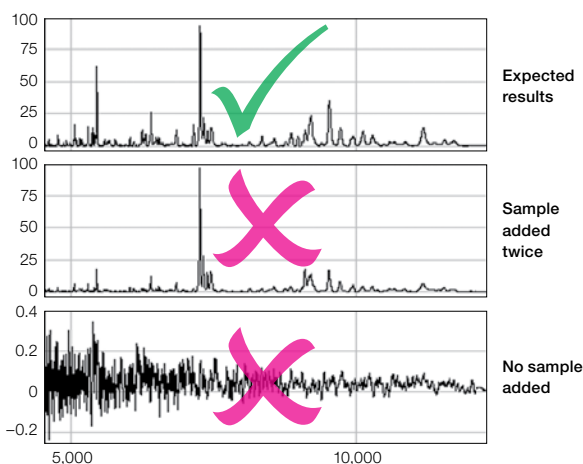
Exercise Guidelines for the ProteinChip System Starter Kit

The starter kit provides all the necessary arrays, reagents, and instructions to perform two exercises, up to three times. Measure your success with feedback on reproducibility with each practice run.

First prepare a ProteinChip NP20 array using serum and *E. coli* samples, then read your array and compare it to the premade NP20 array provided in the kit. The starter kit includes analysis tools that provide you with technical information, such as peak count and signal intensities, and pass or fail results. The manual provides useful information and suggestions on how to improve your technique so that you can obtain better results and a passing score.

Next, generate protein profiles of human serum using the ProteinChip CM10 array. Apply the serum sample to the array, then wash off unbound sample components using ProteinChip CM low-stringency buffer. Apply matrix, read the array, and use the analysis tools included in the kit according to the instructions. As with the NP20 array, measure and monitor your technical improvement until you obtain adequate reproducibility.

With one starter kit, you can prepare ProteinChip arrays while learning the best methods and obtaining technical feedback.



Spectra examples, as included in the manual, help to identify possible technical errors that can cause problems. Data illustrate the results of an error made in the number of additions of sample to the CM10 array during the binding step.

Ordering Information

Catalog #	Description
C70-00068	ProteinChip System Starter Kit , includes 3 each of CM10 and NP20 ProteinChip arrays, 1 each of CM10 and NP20 ProteinChip premade arrays, 1 ProteinChip peptide standard array, 2 vials of ProteinChip SPA matrix, ProteinChip CM low-stringency buffer, acetonitrile, human serum, <i>E. coli</i> , 1% TFA, CD-ROM with video tutorial and protocols, instructions
C70-00070*	ProteinChip Peptide Mass Calibration Kit , includes 1 ProteinChip peptide standard array, sufficient for up to 160 calibrations, CD-ROM with video tutorial and protocol, instructions

* The ProteinChip peptide standard array has a 12-week shelf life after opening.

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 905 364 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