



PROTEIN INTERACTION ANALYSIS

ProteOn™ Protocol Development Kits

- Designed for both new and experienced users of the ProteOn XPR36 system
- Gain familiarity with experimental design and instrument operation
- Use as a benchmark for developing new protocols

Explore the Power of the ProteOn 6 x 6 Interaction Array

Benefits

The ProteOn protocol development kits provide the reagents and sensor chips needed to perform and analyze a complete experiment using the ProteOn XPR36 protein interaction array system. New users will gain familiarity with instrument setup and operation, experimental design, and data analysis. Experienced users will find the kits valuable for system benchmarking and as positive control reagents for protocol development. Each kit contains sufficient materials for a complete exercise in the operation of the ProteOn XPR36 system.

Three Protocol Development Kits

One-shot Kinetics™ kit — The interaction between the cytokine IL-2 and an IL-2 antibody demonstrates a detailed kinetic analysis in a single injection cycle without regeneration (One-shot Kinetics). The kit also demonstrates a useful method for controlling ligand immobilization levels during protocol development and optimization.



Multiple protein interaction kit — The interaction between TEM1 β -lactamase and b-lactamase inhibitor protein (BLIP) demonstrates the power of the ProteOn XPR36 system to produce a detailed kinetic analysis of multiple simultaneous interactions and to map protein interfaces.



Protein-small molecule kit — The interaction between carbonic anhydrase (CA) II and 4-carboxybenzenesulfonamide (CBS) demonstrates the capability of the ProteOn XPR36 system to detect low molecular weight analytes. The kit also provides an additional benchmark of optimal system performance.



BIO-RAD

ProteOn One-shot Kinetics Kit

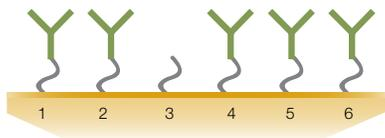
The ProteOn One-shot Kinetics kit demonstrates a complete, detailed kinetic analysis of a protein-protein interaction in a single injection cycle.

This kit also demonstrates the utility of the ProteOn XPR36 system for protocol optimization. Ligand is immobilized to different levels in five parallel channels of a sensor chip, followed by a single injection cycle of an analyte concentration series in the six orthogonal channels (Figures 1 and 2). The third ligand channel is used as a reference channel. A set of sensorgrams providing detailed kinetic analysis data is produced from each interaction channel (Figure 3). Optimal immobilization conditions may thus be rapidly and efficiently determined using the parallel processing capability of the ProteOn XPR36 system. (For details, request bulletin 3172.)

The ProteOn One-shot Kinetics kit includes:

- IL-2/IL-2 antibody pair, with sufficient material to immobilize IL-2 antibody in up to 5 channels and to analyze a complete concentration series of IL-2 across each channel
- GLC sensor chip
- Amine coupling kit
- Acetate buffer, 10 mM, pH 4.5
- Instruction manual

Step 1.



Step 2.

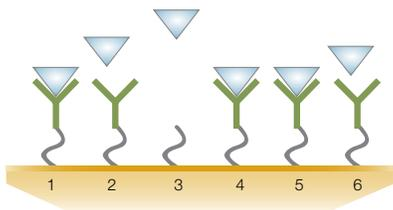


Fig. 1. Workflow of the ProteOn one-shot kinetics kit. Step 1, immobilize IL-2 antibody on chip surface. Step 2, inject IL-2 antigen and monitor interaction with immobilized IL-2 antibody. (Channel 3 is a reference channel.)

Step 1. Immobilize IL-2 antibody.

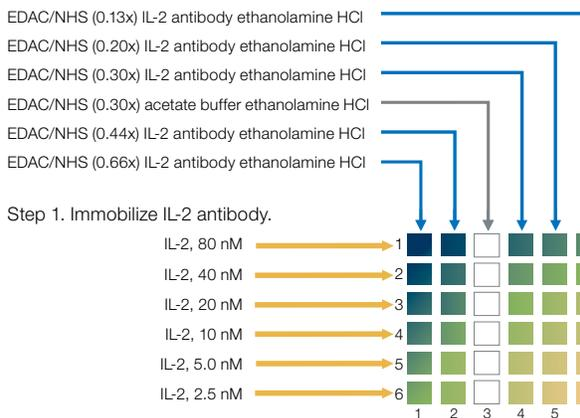


Fig. 2. Protocol optimization using the ProteOn One-shot Kinetics kit. Acetate buffer was at 10 mM, pH 4.5.

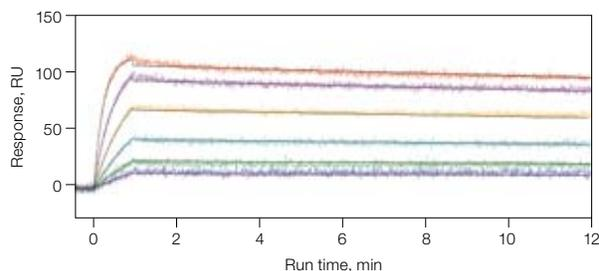


Fig. 3. Detailed kinetic analysis of the IL-2/IL-2 antibody interaction. Results obtained using the ProteOn One-shot Kinetics kit (—, 80 nM; —, 40 nM; —, 20 nM; —, 10 nM; —, 5 nM; —, 2.5 nM).

ProteOn Multiple Protein Interaction Kit

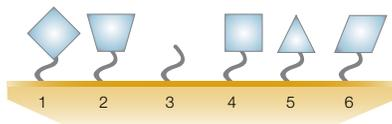
The ProteOn multiple protein interaction kit demonstrates the ProteOn XPR36 system's capability to analyze multiple protein-protein interactions simultaneously.

The kit contains five different mutant proteins of TEM1 and BLIP. Detailed kinetic data are obtained on the interaction of BLIP with each of the five TEM1 mutant proteins in a single injection cycle. The five TEM1 mutant proteins are immobilized in five of the six parallel channels on the sensor chip, and a BLIP concentration series is then injected into the six orthogonal channels (Figures 4 and 5). The third ligand channel is used as a reference channel. Five sets of six sensorgrams are simultaneously produced for rapid analysis and comparison of the binding kinetics of each TEM1 mutant protein with BLIP (Figure 6). (For details, request bulletins 5358 and 5368.)

The ProteOn multiple protein interaction kit includes:

- TEM1/BLIP protein set, with sufficient material to immobilize 5 TEM1 mutant proteins in 5 channels and to perform a detailed kinetic analysis of the interaction of each of the mutant proteins with BLIP
- GLC sensor chip
- Amine coupling kit
- Acetate buffer, 10 mM, pH 4.0
- Instruction manual

Step 1.



Step 2.

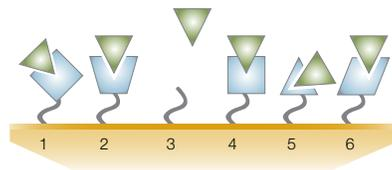
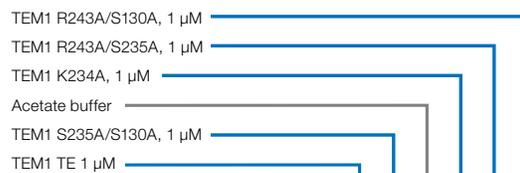


Fig. 4. Workflow of the ProteOn multiple protein interaction kit.

Step 1, immobilize five TEM1 mutant proteins on five channels. Step 2, inject BLIP analyte proteins and monitor interaction with immobilized TEM1 proteins. (Channel 3 is a reference channel.)

Step 1. Immobilize five TEM1 mutant proteins.



Step 2. Inject BLIP analyte protein panel.

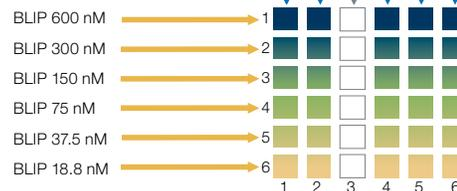


Fig. 5. One-shot Kinetic analysis of the TEM1/BLIP interaction.

Acetate buffer was at 10 mM, pH 4.0.

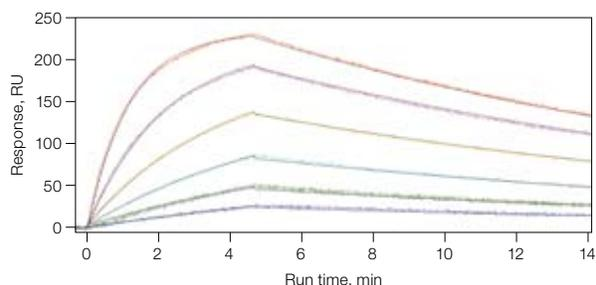


Fig. 6. Detailed kinetic analysis of BLIP protein interacting with the TEM1 mutant protein TEM1 K234A. Shown is a set of six sensorgrams from vertical channel 4 (—, 600 nM; —, 300 nM; —, 150 nM; —, 75 nM; —, 37.5 nM; —, 18.8 nM) obtained in a single injection cycle using the ProteOn multiple protein interaction kit.

ProteOn Protein-Small Molecule Kit

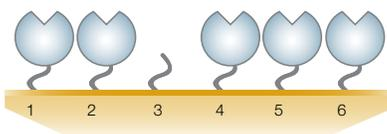
The ProteOn protein-small molecule kit demonstrates the capability of the ProteOn XPR36 system to accurately measure and resolve the kinetics of a small molecule interaction with an immobilized protein.

The ProteOn protein-small molecule kit contains the enzyme inhibitor pair CA II and CBS (Figures 7 and 8). The affinity and kinetics of the interaction between CA II and CBS have been well characterized by SPR technology (Figure 9). The kit also supplies a reagent set designed for assessing the capability of the ProteOn XPR36 system to measure small molecule interactions and for establishing a benchmark for system performance.

The ProteOn protein-small molecule kit includes:

- CA II/CBS pair, with sufficient material to immobilize up to 5 channels with CA II and to analyze a complete concentration series of CBS
- GLC sensor chip
- Amine coupling kit
- Acetate buffer, 10 mM, pH 5.0
- Instruction manual

Step 1.



Step 2.

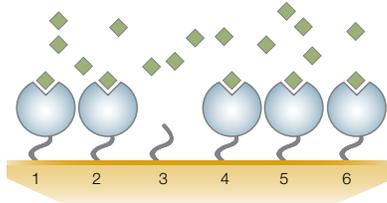


Fig. 7. Workflow of the ProteOn protein-small molecule kit.

Step 1, immobilize CA II on the chip surface. Step 2, inject CBS analyte (MW 201), and monitor interaction with CA II (31 kD). (Channel 3 is a reference channel.)

Step 1. Immobilize CA II.

EDAC/NHS (1.0x) CA II, 0.125 mg/ml ethanolamine HCl
 EDAC/NHS (1.0x) CA II, 0.125 mg/ml ethanolamine HCl
 EDAC/NHS (1.0x) CA II, 0.125 mg/ml ethanolamine HCl
 EDAC/NHS (1.0x) acetate buffer ethanolamine HCl
 EDAC/NHS (0.1x) CA II, 0.125 mg/ml ethanolamine HCl
 EDAC/NHS (0.1x) CA II, 0.125 mg/ml ethanolamine HCl

Step 2. Inject CBS analyte panel.

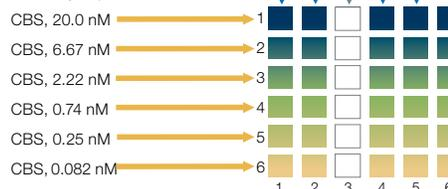


Fig. 8. ProteOn protein-small molecule kit representative protocol. Acetate buffer was at 10 mM, pH 5.0.

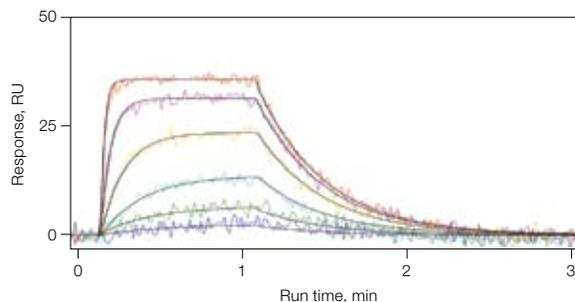


Fig. 9. Detailed kinetic analysis of the CA II/CBS interaction. Results obtained using the ProteOn protein-small molecule kit (—, 20 nM; —, 6.07 nM; —, 2.22 nM; —, 0.74 nM; —, 0.25 nM; —, 0.082 nM).

Ordering Information

Catalog # Description

ProteOn Protocol Development Kits

176-1010	ProteOn One-shot Kinetics Kit , includes IL-2/IL-2 antibody pair, GLC sensor chip, amine coupling kit, 50 ml acetate buffer, pH 4.5, instructions
176-1020	ProteOn Multiple Protein Interaction Kit , includes TEM/BLIP protein set, GLC sensor chip, amine coupling kit, 50 ml acetate buffer, pH 4.0, instructions
176-1030	ProteOn Protein-Small Molecule Kit , includes carbonic anhydrase II/CBS pair, GLM sensor chip, amine coupling kit, 50 ml acetate buffer, pH 5.0, instructions

To obtain bulletins 3172, 5358, and 5368, contact your local Bio-Rad sales representative or download them from www.bio-rad.com.

BIO-RAD

**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com USA 800 424 6723 Australia 61 2 9914 2800 Austria 01 877 89 01 Belgium 09 385 55 11 Brazil 55 31 3689 6600
 Canada 905 364 3435 China 86 20 8732 2339 Czech Republic 420 241 430 532 Denmark 44 52 10 00 Finland 09 804 22 00 France 01 47 95 69 65
 Germany 089 31 884 0 Greece 30 210 777 4396 Hong Kong 852 2789 3300 Hungary 36 1 459 6100 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