

Versatile Separation Capabilities of the PROTEAN® i12™ IEF System

ChengJun Sun, Adriana Harbers, Tom Berkelman, and Katy McGirr,
Life Science Group, Bio-Rad Laboratories Inc., 6000 James Watson Dr.,
Hercules, CA 94547

Introduction

Two-dimensional (2-D) gel electrophoresis is a powerful and widely used analytical tool for resolving and separating complex biological samples (Garfin 2003). The technique employs two sequential electrophoretic separation steps. The first step or dimension is isoelectric focusing (IEF), a separation based on protein charge. The second dimension is SDS-PAGE, a separation based on protein size. Currently, the first dimension is most commonly performed on IPG (immobilized pH gradient) strips, as this provides the best reproducibility and ease of use (Görg et al. 2004). IPG strips are available from Bio-Rad in different lengths (7, 11, 17, 18, and 24 cm) and each of them is available in a variety of pH ranges including broad ranges (pH 3–10 and pH 3–10 nonlinear), narrow ranges (pH 3–6, pH 4–7, pH 5–8, and pH 7–10), and micro ranges (pH 3.9–5.1, pH 4.7–5.9, pH 5.5–6.7, and pH 6.3–8.3). Broad-range IPG strips are best for an initial view, but narrow- and micro-range IPG strips offer higher resolution separations and often allow the visualization of many proteins that cannot be distinguished on broad-range IPG strips. In many cases, it is desirable to analyze the same sample on multiple pH ranges.

IPG-IEF is conducted at high voltage (typically several thousand volts) and the current is typically limited at 50 μ A per IPG strip. Due to these unique running conditions, IPG-IEF is conducted with specialized equipment. Currently available instruments for IPG-IEF are designed to run several IPG strips simultaneously and use a single power supply to

apply voltage across all of the IPG strips. The current limit specified is an average current per IPG strip. If IPG strips of differing conductivity are run on the same instrument, they will experience different current-voltage profiles in an uncontrolled manner. It is therefore recommended that one use only samples of similar composition in the same run. Since individual IPG strip pH ranges have differing intrinsic conductivities, it is also recommended that one restrict a single run to IPG strips of the same pH range.

Prefractionation of samples by liquid-phase IEF in the MicroRotor™ cell is employed prior to 2-D electrophoresis to reduce sample complexity, allow increased protein loading, and facilitate detection of less abundant proteins (Davidsson et al. 2002, Rachinsky et al. 2007). However, this strategy generates fractions of differing conductivities that must be run on different pH-range IPG strips. Analysis of these fractions using current equipment must be done on multiple instruments or sequentially in a time-consuming manner.

Bio-Rad's PROTEAN i12 IEF system represents an improvement over traditional IEF systems by offering individual lane power control. The individual lane control feature of the PROTEAN i12 IEF system allows each IPG strip to be run independently, making it possible to process different samples and different types of IPG strips in a single run. This feature greatly shortens the time needed for protocol optimization for a specific sample. In this study, we demonstrate the versatility and flexibility of the PROTEAN i12 IEF system using soybean protein samples prefractionated on the MicroRotor cell.

Methods

Protein Extraction

Soybeans were hydrated overnight in aerated water and homogenized with a mortar and pestle in a solution composed of 7 M urea, 2 M thiourea, 4% (w/v) CHAPS, and 40 mM DTT. For each gram of tissue, 9 ml of solution was used. The homogenized sample was centrifuged at 14,000 g for 15 min and the supernatant was stored at -80°C until use.

MicroRotorfor Cell Fractionation

Soybean protein samples (1.26 mg) were diluted in 7 M urea, 2 M thiourea, and 0.2% biolyte 3–10 to a final volume of 3 ml and 2.5 ml were loaded onto the MicroRotorfor cell (Bio-Rad Laboratories, Inc.) with 0.1 M H_3PO_4 at the anode and 0.1 M NaOH at the cathode. The run was conducted at 1 W constant power until the voltage stabilized. Total fractionation time was 50 minutes. Ten fractions were harvested at the end of the run. After measuring the pH of each fraction using pH paper, the samples were used directly for the following 2-D experiments.

2-D Electrophoresis

The first dimension was conducted using 11 cm ReadyStrip™ IPG strips pH 3–6, pH 5–8, and pH 7–10 depending on the pH of the samples. Fifteen to 40 μl of each fraction were diluted in a rehydration solution consisting of 7 M urea, 2 M thiourea, 2% (w/v) CHAPS, 0.2% (w/v) Bio-Lyte® ampholyte pH 3–10 (or ReadyStrip 7–10 buffer for pH 7–10 IPG strips), 50 mM DTT, and 0.001% (w/v) bromophenol blue dye to a final volume of 185 μl and loaded onto the strips via passive rehydration. The strips were rehydrated overnight and focused with a program of 250 V for 30 min followed by 8,000 V for 35,000 Vh under a current limit of 50 μA .

Following isoelectric focusing, the IPG strips were equilibrated for 15 min in equilibration buffer I composed of 6 M urea, 50 mM Tris-HCl, pH 8.8, 2% (w/v) SDS, 1% (w/v) DTT, 30% (w/v) glycerol, and 0.001% bromophenol blue. This was followed by another 15 min in equilibration buffer II composed of 6 M urea, 50 mM Tris-HCl, pH 8.8, 2% (w/v) SDS, 2.5% (w/v) iodoacetamide, 30% (w/v) glycerol, and 0.001% bromophenol blue. Second-dimension separations for 11 cm IPG strips were performed on Criterion™ 8–16% Tris-HCl gels. Equilibrated IPG strips were applied to the second-dimension gels and run using Tris-Glycine-SDS buffer at 200 V for 55 min.

Gel Staining and Imaging

After running, all gels were washed three times for 5 min with deionized water and stained with Bio-Safe™ Coomassie stain (Bio-Rad) for 1 hr. After destaining with water, the gels were scanned on a GS-800™ calibrated densitometer using standard settings (Bio-Rad).

Results

MicroRotorfor fractionation

Adjacent fractions collected from the MicroRotorfor cell differed from each other by 0.5 to 1.0 pH units, as judged using pH paper, and covered most of the pH range from pH 3 to pH 10 (Table 1). Narrow-range IPG strips for subsequent 2-D analysis were chosen based on sample pH (Table 1). The acidic, neutral to near-neutral, and basic fractions were focused on pH 3–6, pH 5–8, and pH 7–10 ReadyStrip IPG strips, respectively.

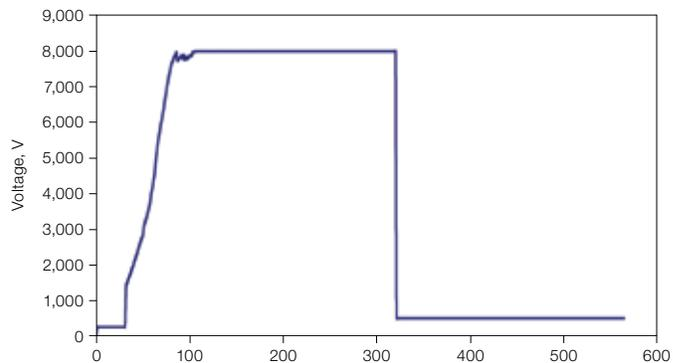
Table 1. MicroRotorfor fraction pH and corresponding IPG strips used.

Fraction #	pH	pH range of IPG strips
1	3.0	3–6
2	4.0	3–6
3	4.5	3–6
4	5.5	5–8
5	6.0	5–8
6	7.0	5–8
7	7.5	5–8
8	8.5	7–10
9	9.0	7–10
10	>9.0	7–10

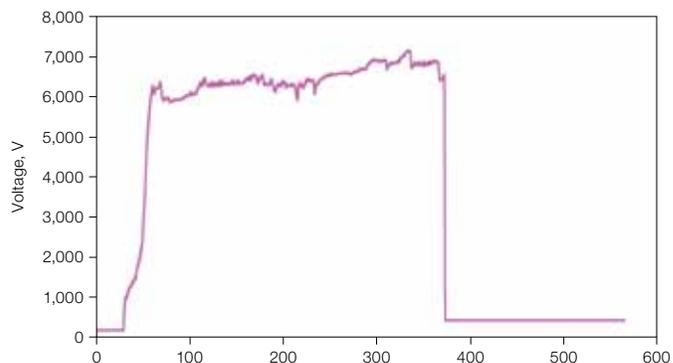
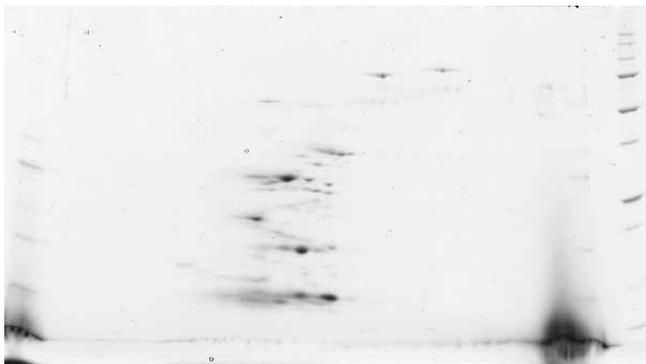
PROTEAN i12 IEF System Performance with Different Samples and Different IPG Strips in a Single Run

To test the flexibility of the PROTEAN i12 IEF system, we tested its performance by simultaneously running all the fractions from a MicroRotorfor separation. As shown in Table 1, each fraction has a distinct pH. Protein composition and resistance are therefore expected to differ from fraction to fraction. Indeed, each sample gave a distinctive voltage profile when focused with a 50 μA current limit. In a traditional focusing system, the current through each IPG strip would have been uncontrolled, resulting in unpredictable results. On the PROTEAN i12 IEF system, with individual power control for each lane, there is no effect of the other concurrently run IPG strips on any of the individual runs. As shown in Figure 1, samples ranging from very acidic to very basic pH loaded on IPG strips with different pH gradients can all be focused in one run. In addition, the gel images show that different amounts of proteins were loaded on each strip. Therefore, the PROTEAN i12 IEF system can accommodate differences in sample pH, composition, and load, as well as different IPG pH gradients, at the same time.

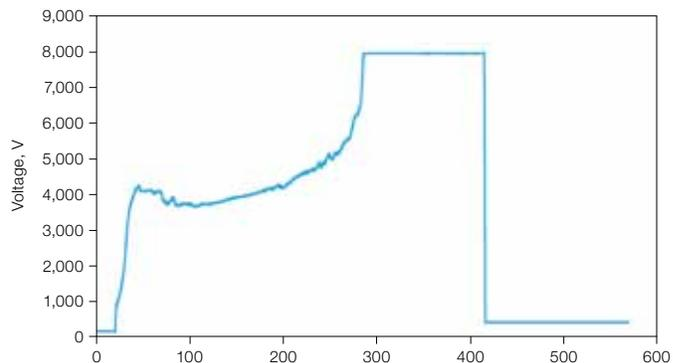
Fraction #1 run on pH 3-6 strip.



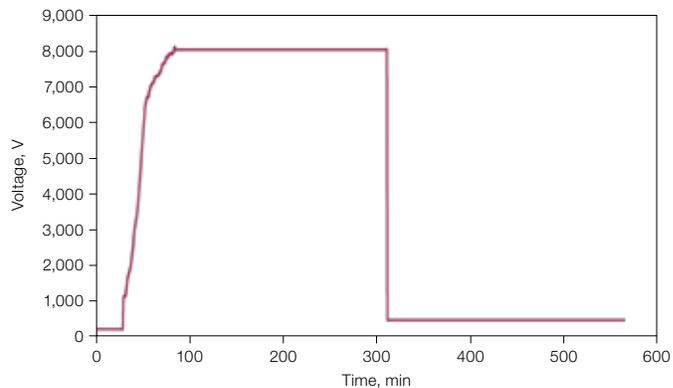
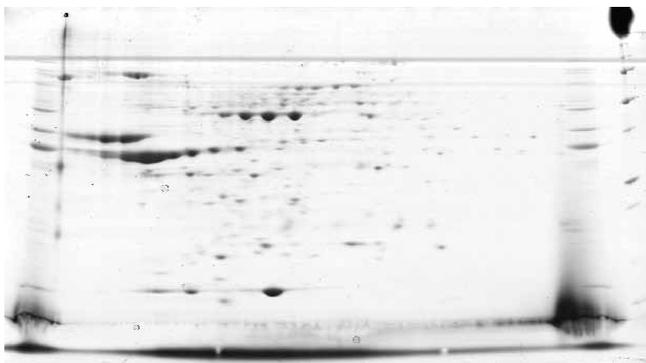
Fraction #2 run on pH 3-6 strip.



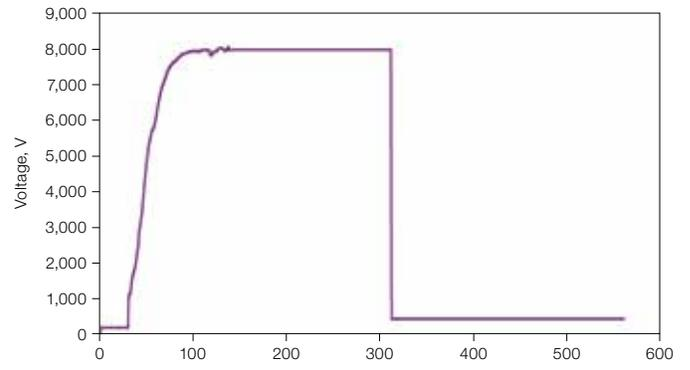
Fraction #3 run on pH 3-6 strip.



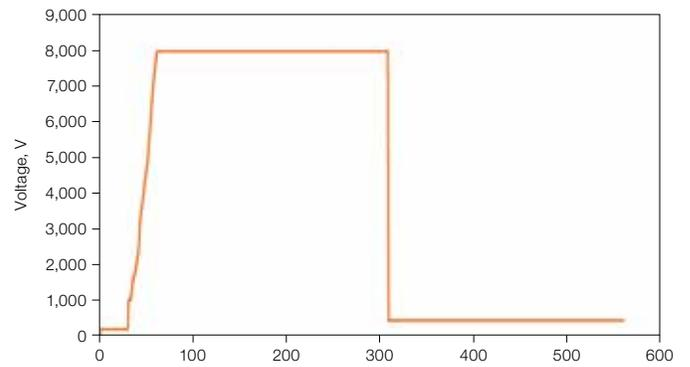
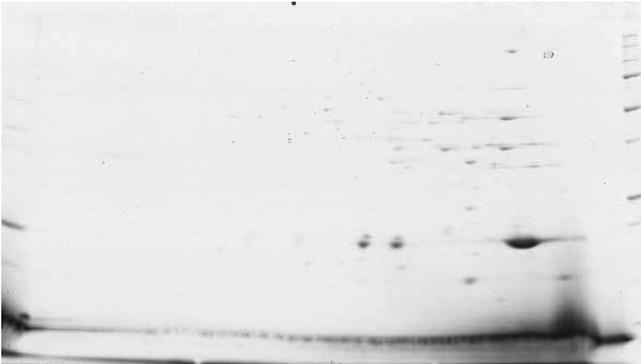
Fraction #4 run on pH 5-8 strip.



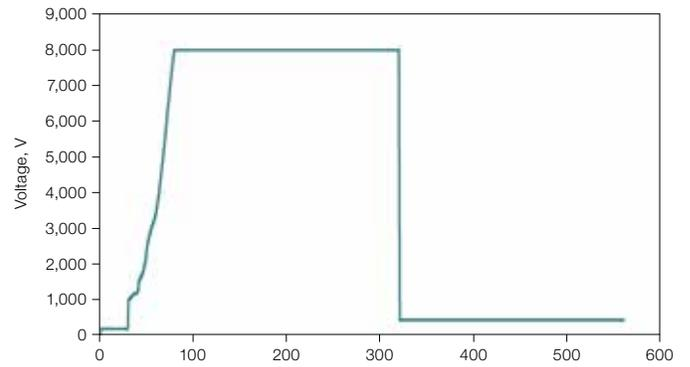
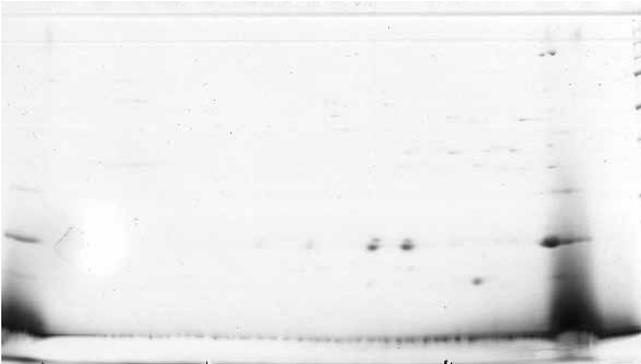
Fraction #5 run on pH 5-8 strip.



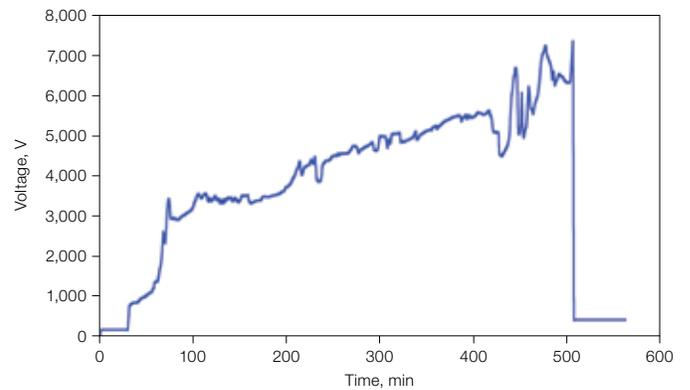
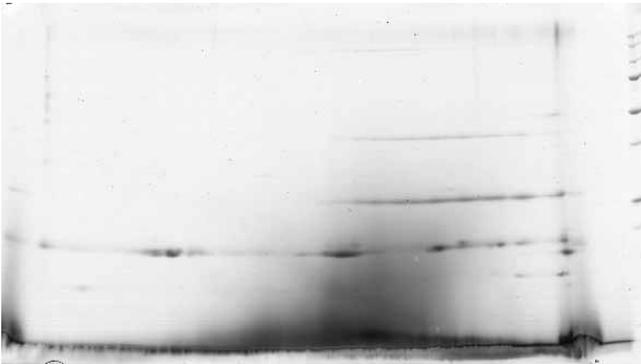
Fraction #6 run on pH 5-8 strip.



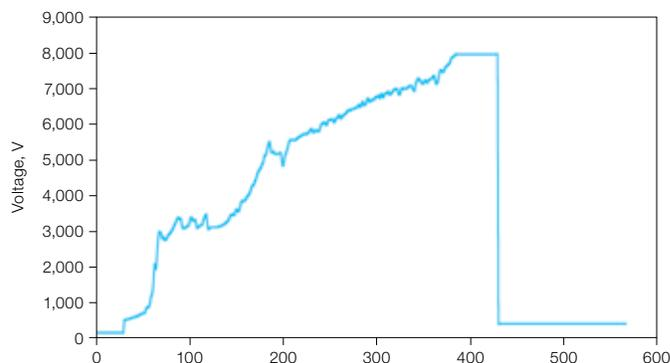
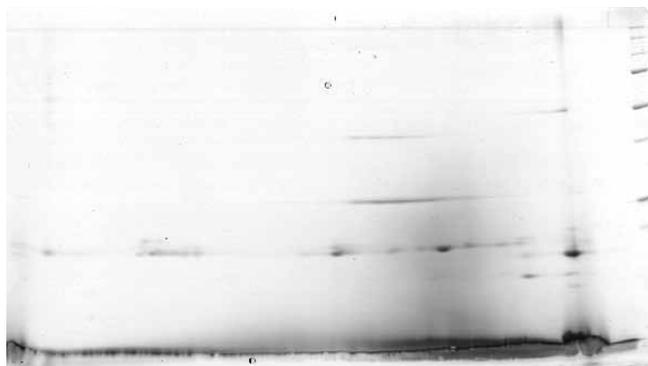
Fraction #7 run on pH 5-8 strip.



Fraction #8 run on pH 7-10 strip.



Fraction #9 run on pH 7–10 strip.





**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 21 6169 8500 **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 **Malaysia** 60 3 2117 5260 **Mexico** 52 555 488 7670
The Netherlands 0318 540666 **New Zealand** 64 9 415 2280 **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 3170 **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 **Thailand** 66 2 6518311 **United Kingdom** 020 8328 2000