

AUTOMATED ELECTROPHORESIS

Experion™ Pro260 Analysis Kit

- For automated electrophoresis, detection, and basic analysis with the Experion system
- Analysis of 10 samples in as little as 30 minutes
- 33% higher resolution than other automated systems
- Protein quantitation, sizing, and analysis in a single step
- Innovative Caliper Life Sciences' LabChip technology



Sensitive, Reliable Protein Analysis in 30 Minutes

Introduction

Experion Pro260 analysis kits combine state-of-the-art chip design with high-quality reagents to perform reproducible, quantitative, and accurate protein analysis in a matter of minutes with the Experion automated electrophoresis system. Streamlined chip preparation methods and minimal sample requirements result in rapid experiments with minimal hands-on time.

The Experion Pro260 Advantage

Innovative chip design, chemistry, software, and hardware make protein separation and quantitation fast and easy.

- Analysis of 10 samples in as little as 30 minutes — electrophoresis, staining, destaining, detection, and basic analysis in a single step
- Resolution and quantitation of proteins between 10–260 kD
- 33% higher resolution than other automated systems
- Sensitivity comparable to colloidal Coomassie Blue staining of SDS-PAGE gels and over 3 times greater than other automated electrophoresis systems
- Linear dynamic range of 5–2,000 µg/ml BSA
- Accurate sizing with the Experion Pro260 protein ladder, part of the Precision Plus Protein™ family of protein standards

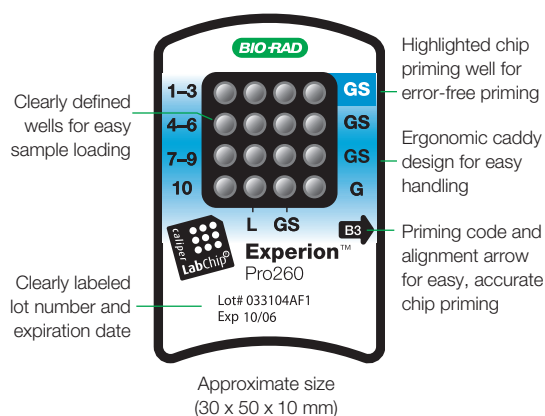
Experion Pro260 Chips

Experion Pro260 protein chips offer an optimized microchannel architecture within an ergonomic caddy for exceptional protein separation, chip handling, and results. Each chip holds up to ten samples and is clearly labeled with well number, priming code, lot number, and expiration date for easy, error-free runs.

Experion Pro260 Reagents

Bio-Rad has applied its expertise in protein separation and detection to the development of gel, buffer, and dye reagents for the Experion Pro260 analysis kits. Experion Pro260 analysis kits include the following reagents and supplies:

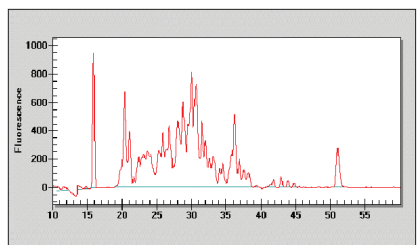
- High-quality gel matrix for separation and resolution similar to mini gels
- Fluorescent dye for high-sensitivity and accurate detection
- Experion Pro260 protein ladder for accurate sizing
- Optimized sample buffer for accurate quantitation and reproducible results
- Spin filters for convenient gel preparation



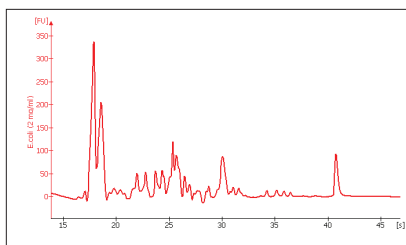
BIO-RAD



Experience Meets Innovation



Run time, sec



Run time, sec

Comparison of separation of proteins from equal amounts of *E. coli* lysate on automated electrophoresis systems.

Left panel, separation on the Experion system using the Pro260 kit; right panel, separation on a competitor's automated electrophoresis system using a competitor's kit. Note the greater number of peaks and increased resolution of the Experion Pro260 result.

Chemical Compatibility

The Experion Pro260 analysis kit delivers reproducible results with samples in all of the following buffers. For a complete chemical compatibility list, refer to the Experion Pro260 analysis kit instruction manual.

Buffer	Formulation	Buffer	Formulation
IEF cathode buffer	30 mM lysine, 30 mM arginine	IMAC elution buffer	50 mM potassium phosphate, 150 mM sodium chloride, 250 mM imidazole, pH 8
Tris/Tricine/SDS	150 mM Tris, 150 mM Tricine, 0.15% SDS, pH 8.3	Protein A resin wash buffer I	50 mM sodium acetate, 250 mM sodium chloride, pH 4
Tris/glycine/SDS	37.5 mM Tris, 288 mM glycine, 0.15% SDS, pH 8.3	Protein A resin wash buffer II	50 mM sodium bicarbonate, 250 mM sodium chloride, pH 8
Tris/CAPS	90 mM Tris, 60 mM CAPS, 0.15% SDS, pH 9.6	Tris-borate-EDTA (TBE)	133 mM Tris, 133 mM boric acid, 3 mM EDTA, pH 8.3
Tricine sample buffer	100 mM Tris-HCl, pH 6.8, 1% SDS, 20% glycerol, 0.02% Coomassie Blue	PBS-Tween	10 mM sodium phosphate, 150 mM sodium chloride, 0.05% Tween, pH 7.2
Zymogram sample buffer	31.3 mM Tris-HCl, pH 6.8, 2% SDS, 12.5% glycerol, 0.005% Bromophenol Blue	Commercial Buffers	
Phosphate-buffered saline (PBS)	10 mM sodium phosphate, 150 mM sodium chloride, pH 7.5	XT MOPS buffer (1.5x)	
		XT Tricine buffer (1.5x)	
		XT sample buffer	
		BugBuster (Novagen)	
		B-PER (Pierce)	

Ordering Information

Catalog #	Description
700-7101	Experion Pro260 Analysis Kit for 10 Chips, includes 10 Pro260 chips, 3 x 520 µl Pro260 gel, 45 µl Pro260 stain, 60 µl Pro260 ladder (10–260 kD), 400 µl Pro260 sample buffer, 3 spin filters
700-7102	Experion Pro260 Analysis Kit for 25 Chips, includes 25 Pro260 chips, 5 x 520 µl Pro260 gel, 2 x 45 µl Pro260 stain, 2 x 60 µl Pro260 ladder (10–260 kD), 2 x 400 µl Pro260 sample buffer, 5 spin filters
700-7151	Experion Pro260 Chips, 10
700-7152	Experion Pro260 Reagents and Supplies, for 10 chips, includes 3 x 520 µl Pro260 gel, 45 µl Pro260 stain, 60 µl Pro260 ladder (10–260 kD), 400 µl Pro260 sample buffer, 3 spin filters
700-7251	Experion Cleaning Chips, 10
700-7256	Experion Pro260 Ladder, 60 µl (10–260 kD)
700-7110	Experion Pro260 Starter Kit
700-7000	Experion System, 100/240 V, for protein analysis, includes electrophoresis station, priming station, software, USB2 cable, test chip, instructions (analysis kits sold separately)



LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only. These products are licensed under US Patent Nos. 5,863,753, 5,658,751, 5,436,134, and 5,582,977, and pending patent applications, and related foreign patents, for internal research and development use only in detecting, quantitating, and sizing macromolecules, in combination with microfluidics, where internal research and development use expressly excludes the use of this product for providing medical, diagnostic, or any other testing, analysis, or screening services, or providing clinical information or clinical analysis, in any event in return for compensation by an unrelated party.

B-PER is a registered trademark of Thermo Fisher Scientific, Inc. BugBuster is a registered trademark of EMD Chemicals, Inc. Coomassie is a trademark of BASF Aktiengesellschaft. Tween is a trademark of ICI Americas Inc.

BIO-RAD

**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