



# Experion™

Automated Electrophoresis System

## Focus on the results, not the method.

- Automated separation, staining, destaining, imaging, band detection, quantitation, and basic data analysis of protein and RNA samples in just 30 minutes
- Significantly reduced time to results, hands-on time, reagent usage, and sample consumption
- Single-step protein size determination (10–260 kD) or RNA purity checks at nanogram or picogram levels



**BIO-RAD**

# Experience Speed in Automation

The Experion system delivers affordable, accessible automation for electrophoresis. Combining Bio-Rad's expertise in electrophoresis with the innovation of Caliper Life Sciences' LabChip microfluidic technology, the Experion system delivers outstanding performance. With the Experion system, your routine protein and RNA separations will yield reproducible results in just 30 minutes.

## Rapid, Automated Results

The Experion system automatically performs the multiple steps of gel-based electrophoresis. You can walk away and do more with your time while the system separates and quantitates your protein and RNA samples.

## Integrated System Design

The components of the Experion system incorporate efficient and creative designs to deliver high-quality results. Optimized microfluidic chip design, electrophoresis-grade reagents, exclusive protein and RNA standards, easy-to-use automated priming and electrophoresis stations, and powerful software analysis tools all combine to form an integrated system that streamlines separation and analysis.

## Optimal Performance and User-Friendly Tools

- The highest protein resolution and sensitivity and best quantitation results available in an automated system
- Resolution and sensitivity comparable to colloidal Coomassie mini-gels
- Accurate single-step protein sizing from 10 to 260 kD
- Significantly reduced time to results, hands-on time, reagent usage, and sample consumption
- Accurate RNA quantitation at nanogram levels
- 2-in-1 process for RNA: integrity checks and quantitation
- Quick comparisons of protein or RNA components across all samples in a chip
- Convenient software tools for easy and efficient data analysis — sizing and quantitation calculations are performed automatically



### Experion automated electrophoresis system components.

1, operating software and data analysis tools; 2, electrophoresis station; 3, priming station; 4, vortex station; 5, RNA HighSens analysis chip (picogram sensitivity range); 6, RNA StdSens analysis chip (nanogram sensitivity range); 7, Pro260 protein analysis chip; 8, analysis kit reagents for protein and RNA samples.

## Ordering Information

Catalog #	Description
700-7000	<b>Experion System</b> , 100–120/220–240 V, for protein analysis, includes electrophoresis station, priming station, software, USB2 cable, test chip, instructions (analysis kits sold separately)
700-7001	<b>Experion System</b> , 100–120 V, for RNA analysis, includes electrophoresis station, priming station, vortex station, software, USB2 cable, test chip, instructions (analysis kits sold separately)
700-7002	<b>Experion System</b> , 220–240 V, for RNA analysis, includes electrophoresis station, priming station, vortex station, software, USB2 cable, test chip, instructions (analysis kits sold separately)
700-7102	<b>Experion Pro260 Analysis Kit for 25 Chips</b> , 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
700-7104	<b>Experion RNA StdSens Analysis Kit for 25 Chips</b> , includes 25 RNA StdSens chips, 2 x 1,250 µl RNA gel, 2 x 20 µl RNA StdSens stain, 2 x 20 µl RNA ladder, 2 x 900 µl RNA StdSens loading buffer, 4 spin filters
700-7106	<b>Experion RNA HighSens Analysis Kit for 25 Chips</b> , includes 25 RNA HighSens chips, 2 x 1,250 µl RNA gel, 2 x 20 µl RNA HighSens stain, 20 µl RNA ladder, 2 x 900 µl RNA HighSens loading buffer, 2 x 100 µl RNA sensitivity enhancer, 4 spin filters



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 U.S. 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.

**BIO-RAD**

**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

Web site [www.bio-rad.com](http://www.bio-rad.com) USA (800) 4BIORAD Australia 02 9914 2800 Austria (01)-877 89 01 Belgium 09-385 55 11 Brazil 55 21 2527 3454 Canada (905) 712-2771 China (86 21) 6426 0808 Czech Republic + 420 2 41 43 05 32 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)-2398112/3/4, 5018111, 6450092/93 Israel 03 951 4127 Italy 39 02 216091 Japan 03-5811-6270 Korea 82-2-3473-4460 Latin America 305-894-5950 Mexico 55-52-00-05-20 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 7 095 721 1404 Singapore 65-64153188 South Africa 00 27 11 4428508 Spain 34 91 590 52 00 Sweden 08 555 12700 Switzerland 061 717 95 55 Taiwan (886 2) 2578 7189/2578 7241 United Kingdom 020 8328 2000