

# Why choose the **Experion™ Automated Electrophoresis System** from Bio-Rad?

**These key features make the Experion system the obvious choice:**

### Reproducibility

- Automated chip priming station
  - you'll never waste a chip
  - intra- and inter-chip data are highly reproducible

### Standardized RNA Integrity Analysis

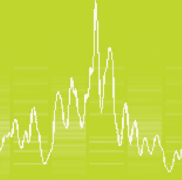
- Simplified RNA integrity assessment via RNA quality indicator (RQI)
- Customizable, color-coded RNA integrity categories
- 28S/18S ribosomal RNA ratio
- RNA sample electropherogram profile



### Unique and Flexible Software

- At-a-glance results display
- User-friendly graphical interface
- Exportable run summary tables
- Concise export options for Excel and database users
- Quick statistics without exporting
- Flexibility to assign and refine peak areas
- Quick and easy electropherogram annotation for presentations and publication
- 5 different methods for protein quantitation
  - relative to 260 kD internal standard
  - relative to user defined internal standard
  - calibration curve absolute quantitation
  - calibration curve and user defined internal standard
  - % total

**BIO-RAD**



### For Core Lab and Pharmaceutical Customers

- IQ/OQ validation software package available
- Security package (for 21 CFR Part 11 compliance)
- High reproducibility of intra- and inter-chip data

### Technical Support

Microfluidics may be a new area for your lab; let us make your transition smoother:

- Starter kits with control sample
- Field application specialists (FAS)
- Training videos
- Telephone calls answered by in-house technical support team: 1-800-4BIORAD (1-800-424-6723)

Excel is a trademark of Microsoft Corporation.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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