

AUTOMATED ELECTROPHORESIS

Experion™ RNA HighSens and RNA StdSens Analysis Kits

- For automated electrophoresis, detection, and basic analysis with the Experion system
- Analysis of 1–12 samples in as little as 10–30 minutes
- Accurate RNA quantitation
- Single-step RNA integrity and concentration determination
- Innovative Caliper Life Sciences' LabChip technology



Single-Step RNA Analysis in 30 Minutes

Introduction

Experion RNA HighSens and RNA StdSens analysis kits combine state-of-the-art chip design with high-quality reagents to perform reproducible, quantitative, and accurate RNA 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 RNA HighSens and RNA StdSens Advantage

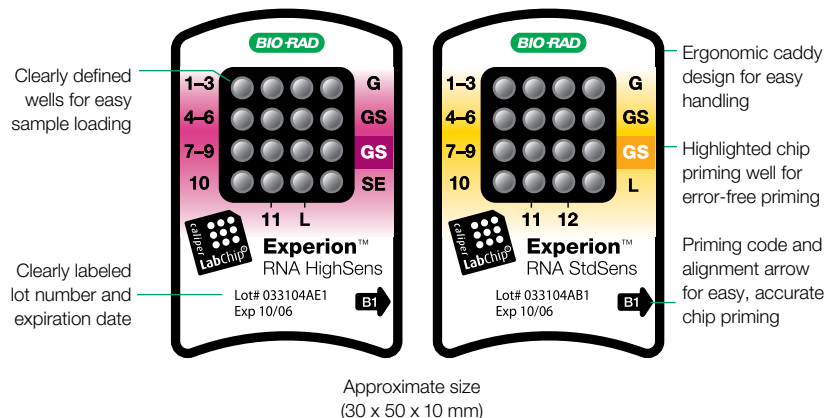
Innovative chip design, chemistry, software, and hardware make RNA separation and quantitation fast and easy.

- Single-step determination of RNA integrity (with the RQI-RNA quality indicator number)
- RNA ladder included in each kit
- Analysis of 1–12 samples in as little as 10–30 minutes

- Accurate quantitation of nanogram amounts of RNA
- Sensitivity down to 100 pg/μl with the RNA HighSens kit, dynamic range of 100–5,000 pg/μl
- Sensitivity down to 5 ng/μl with the RNA StdSens kit, dynamic range of 5–500 ng/μl

Experion RNA HighSens and RNA StdSens Chips

Experion RNA chips offer an optimized microchannel architecture within an ergonomic caddy for exceptional RNA separation, chip handling, and results. Each RNA HighSens chip holds up to 11 samples, and each RNA StdSens chip holds up to 12. The chips are clearly labeled with well number, priming code, lot number, and expiration date for easy, error-free runs.



BIO-RAD



Experience Meets Innovation

Experion RNA HighSens and RNA StdSens Reagents

Accurate assessment of RNA integrity is essential to the success of quantitative RT-PCR and microarray experiments. Experion RNA analysis kits offer single-step RNA integrity determination for confidence without delay. Experion RNA analysis kits include the following reagents and supplies:

- High-quality gel matrix for improved separation compared to agarose RNA gels, with significantly lower sample requirements
- Fluorescent dye for high-sensitivity and accurate detection
- Exclusive Experion RNA ladder for accurate sizing and quantitation
- Optimized loading buffer for reproducible results
- Sensitivity enhancer in RNA HighSens kit for accurate analysis at picogram levels
- Spin filters for convenient gel preparation

Ordering Information

Catalog #	Description
700-7103	Experion RNA StdSens Analysis Kit for 10 Chips, includes 10 RNA StdSens chips, 1,250 µl RNA gel, 20 µl RNA StdSens stain, 20 µl RNA ladder, 900 µl RNA StdSens loading buffer, 2 spin filters
700-7104	Experion RNA StdSens Analysis Kit for 25 Chips, 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
700-7105	Experion RNA HighSens Analysis Kit for 10 Chips, includes 10 RNA HighSens chips, 1,250 µl RNA gel, 20 µl RNA HighSens stain, 20 µl RNA ladder, 900 µl RNA HighSens loading buffer, 100 µl RNA sensitivity enhancer, 2 spin filters
700-7106	Experion RNA HighSens Analysis Kit for 25 Chips, 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
700-7111	Experion RNA StdSens Starter Kit
700-7001	Experion System, 100–240 V, for RNA analysis, includes electrophoresis station, priming station, vortex station, software, USB2 cable, test chip, instructions (analysis kits sold separately)
700-7252	Experion Electrode Cleaner, 250 ml
700-7253	Experion DEPC-Treated Water, 100 ml



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.



**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