

Storage Conditions


Product	Storage
Pro260 chips	Room temperature
Pro260 reagents	4°C (protected from light)

Essential Practices

- Always wear gloves when handling reagents and chips
- Handle chips by the edges; do not touch the glass
- Remove chip from packaging only immediately before use
- Avoid sources of dust and contaminants when preparing samples and loading the chip.  
Foreign particles in reagents, samples, and the wells of the chip can interfere with results
- Deep-clean the electrodes if contamination is suspected or if a chip was left in the instrument overnight
- Use of colored or coated (for example, siliconized polypropylene) tubes when preparing kit reagents or samples is not recommended; such tubes may cause artifacts during the separation
- Refer to the instruction manual for more details

Ordering Information

Catalog #	Description	Catalog #	Description
700-7000	<b>Experion System</b> , 100–240 V, for protein analysis, includes electrophoresis station, priming station, software, USB2 cable, instructions (analysis kits sold separately)	700-7110 700-7151 700-7152	<b>Experion Pro260 Starter Kit</b> <b>Experion Pro260 Chips</b> , 10 <b>Experion Pro260 Reagents and Supplies</b> , 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
700-7101	<b>Experion Pro260 Analysis Kit for 10 Chips</b> , includes 10 Pro260 chips, Experion Pro260 reagents and supplies for 10 chips	700-7251	<b>Experion Cleaning Chips</b> , 10
700-7102	<b>Experion Pro260 Analysis Kit for 25 Chips</b> , includes 25 Pro260 chips, Experion Pro260 reagents and supplies for 25 chips	700-7254 163-2091 161-0610	<b>Experion Spin Filters</b> , 10 <b>ReadyPrep™ Proteomics Grade Water</b> , 500 ml <b>Dithiothreitol (DTT) reducing agent</b> , 1g
701-7000	<b>Experion System</b> , 100–240 V, for protein analysis (700-7000), Experion Pro260 starter kit (700-7110)	161-0710 700-7256	<b>2-Mercaptoethanol</b> , 25 ml <b>Experion Pro260 Ladder</b>

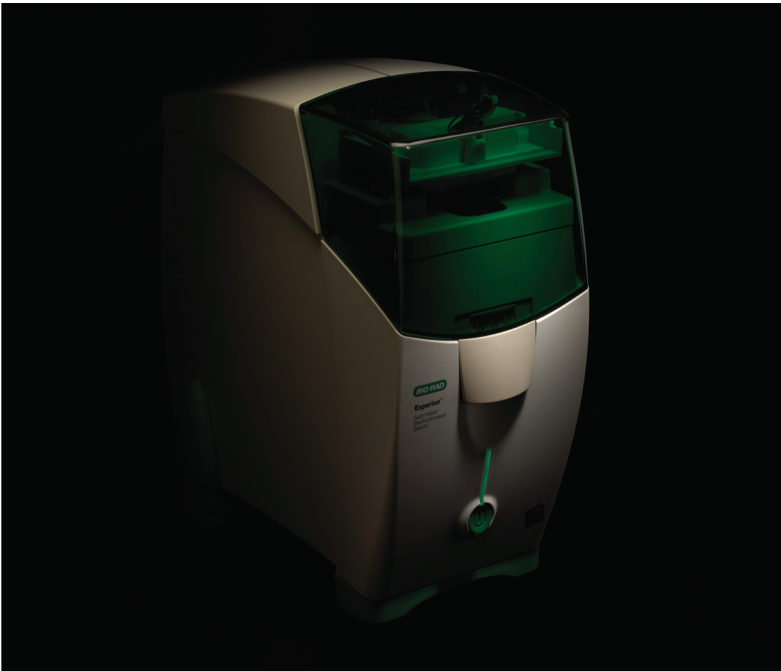
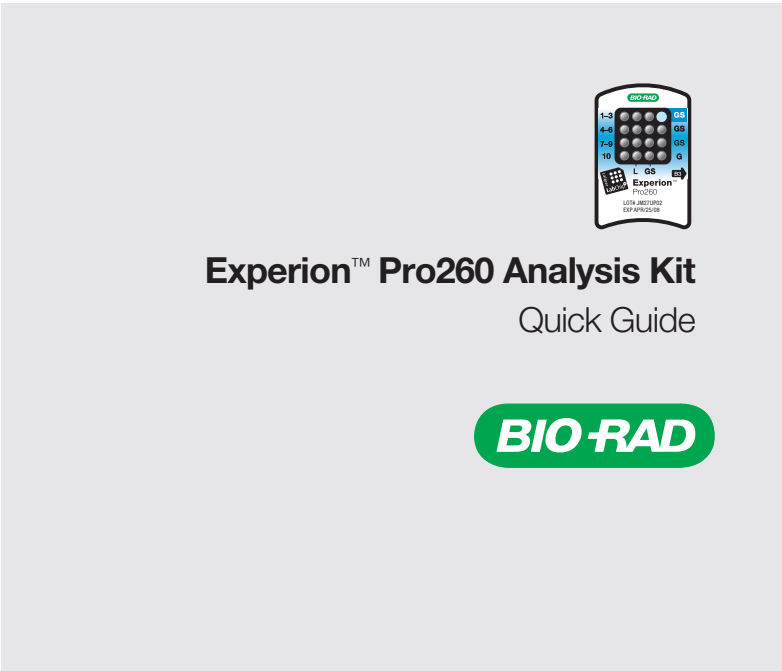
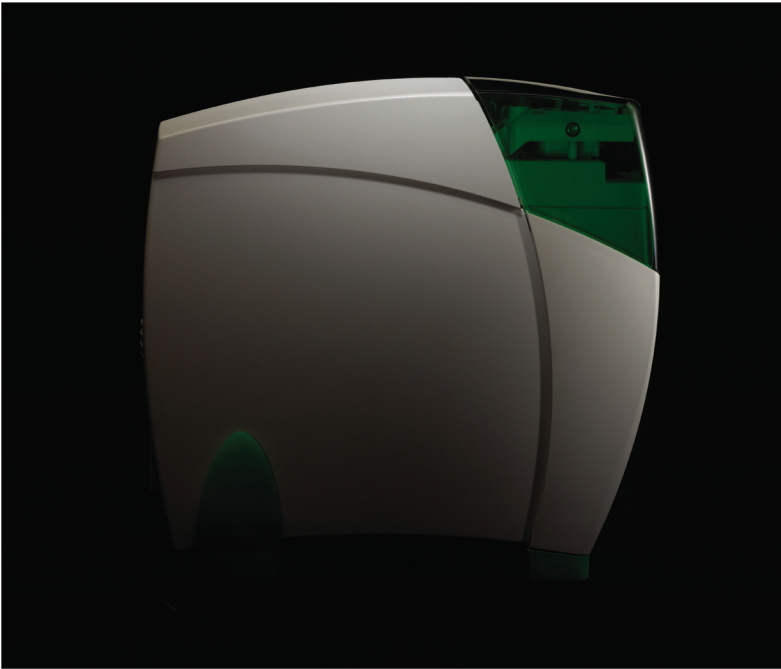
 LabChip and the LabChip logo are trademarks of Caliper Life Sciences, Inc. Bio-Rad Laboratories, Inc. is licensed by Caliper Life Sciences, Inc. to sell products using the LabChip technology for research use only. These products are licensed under U.S. Patent Nos. 5,863,753, 5,658,751, 5,436,134, and 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** 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



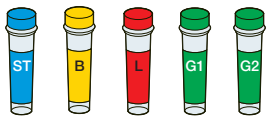
# Experion Pro260 Analysis Kit Quick Guide

For complete instructions, refer to the Experion Pro260 analysis kit instruction manual. Full manuals are available online at [www.bio-rad.com](http://www.bio-rad.com) or contact us by phone at 1 800 424 6723 for an electronic copy. Read the full protocol and essential practices sections if using for the first time.

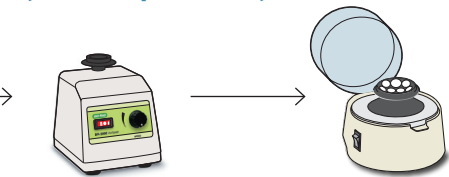


1

## Equilibrate Kit Reagents (ST = stain, B = sample buffer, L = ladder, G = gel)



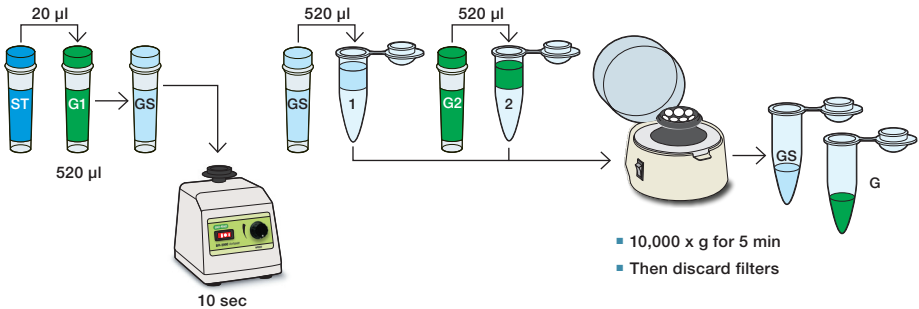
- Equilibrate reagents to room temperature (~15 min). Protect them from light



- Briefly vortex and spin down reagents

2

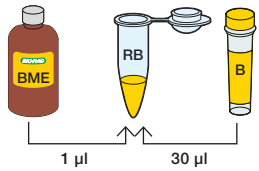
## Prepare the Gel (G) and Gel-Stain (GS) Solution



**Note:** Unused filtered GS solution and filtered G may be stored for up to 30 days at 4°C protected from light; each can be used as is within 30 days. After 30 days, GS solution and G must be refiltered before use. GS solution and G may be stored (4°C protected from light) for another 30 days and refiltered before use. Discard GS solution and G after the second refiltering.

3

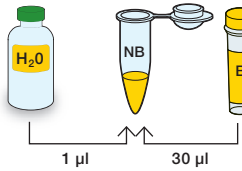
## Prepare the Sample Buffer Reducing Conditions



- Prepare 1 tube of reducing buffer (RB) for each chip to be run
- The ladder must be prepared with RB
- Prepare a fresh solution daily

**Note:** β-mercaptoethanol (BME) or dithiothreitol (DTT) can be used.

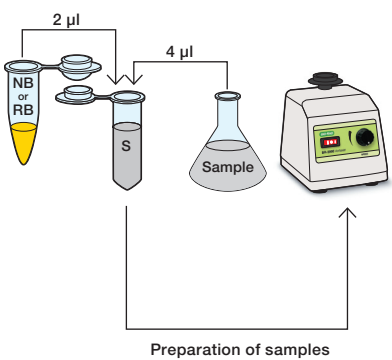
## Nonreducing Conditions



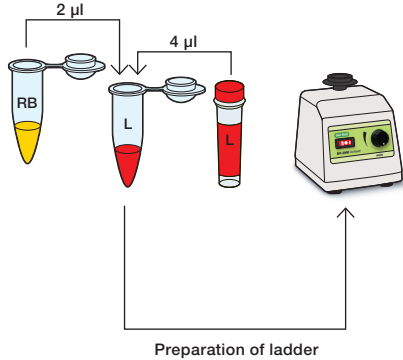
- Prepare 1 tube of nonreducing buffer (NB) for each chip to be run
- Samples prepared with RB and NB can be run on the same chip
- Only use ultrapure water (0.2 µm, filtered)

4

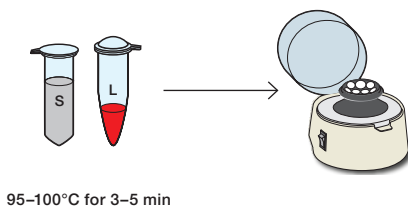
## Prepare the Samples and Pro260 Ladder



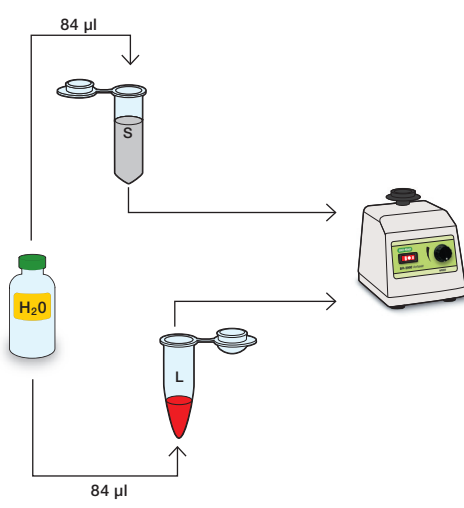
Preparation of samples



Preparation of ladder



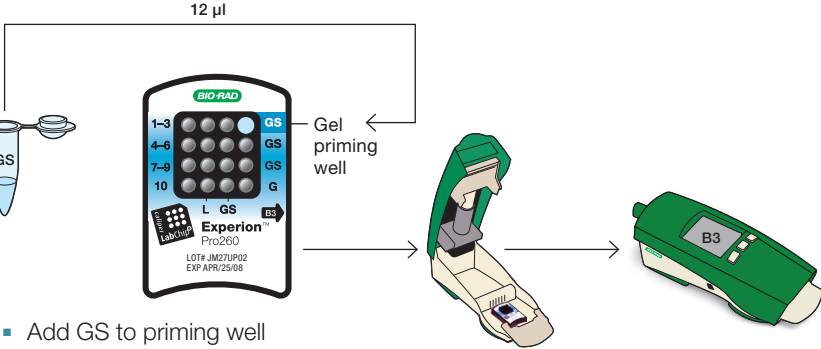
95–100°C for 3–5 min



**Note:** Use only ultrapure water for the 84 µl dilution step.

5

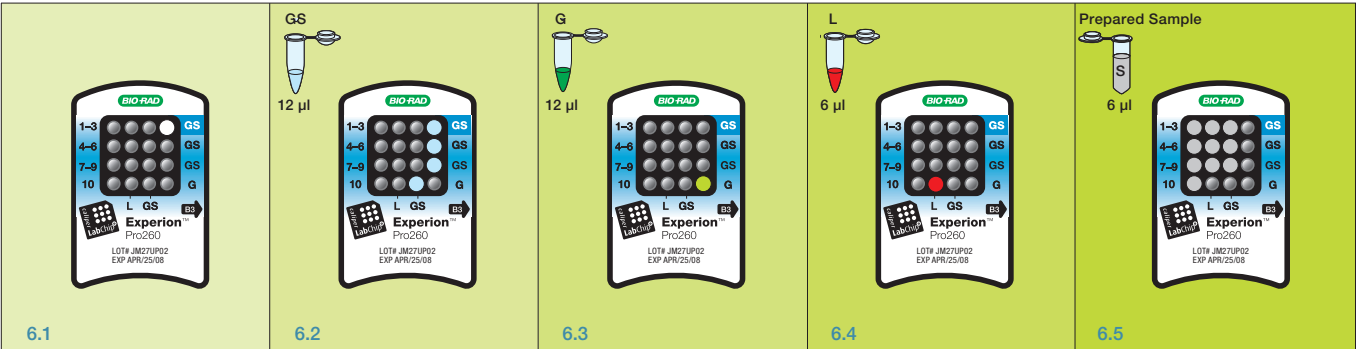
## Prime the Chip



- Add GS to priming well
- Select B3 on priming station
- Place chip in station and press **Start**
- Remove chip after priming is complete
- Flip chip over and visually inspect the microchannels for trapped air bubbles or incomplete priming

6

## Load the Prepared Samples and Pro260 Ladder Onto the Chip



- Remove remaining GS solution from the **GS** well (gel priming well)
- Pipet 12 µl filtered GS solution into all 4 **GS** wells
- Pipet 12 µl filtered G into well **G**
- Pipet 6 µl diluted protein ladder into well **L**
- Pipet 6 µl diluted sample into wells 1–10

**Note:** Do not leave any sample well empty. If necessary, pipet a replicate into any empty sample wells; run the chip in the Experion electrophoresis station within 5 min of loading.

7

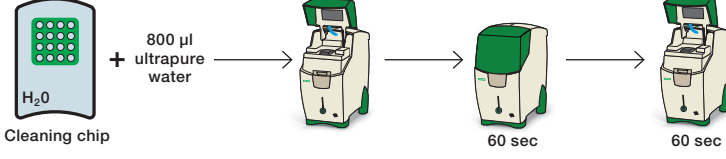
## Run the Pro260 Analysis



- Select **New Run**, then **Protein 260** assay
- Click the **Start** (▶) button onscreen
- Select number of samples to run
- When the run is complete, remove and discard the used chip

8

## Clean the Electrodes



**Note:** If using the instrument for RNA analysis after a protein run, use the deep-cleaning procedure instructions in the RNA analysis kit manual or the software version 3.0 Help section (use search term: electrodes).