

Storage Conditions

Product	Storage
DNA 1K chips	Room temperature
DNA 1K 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 (Bulletin 10006453) for more details

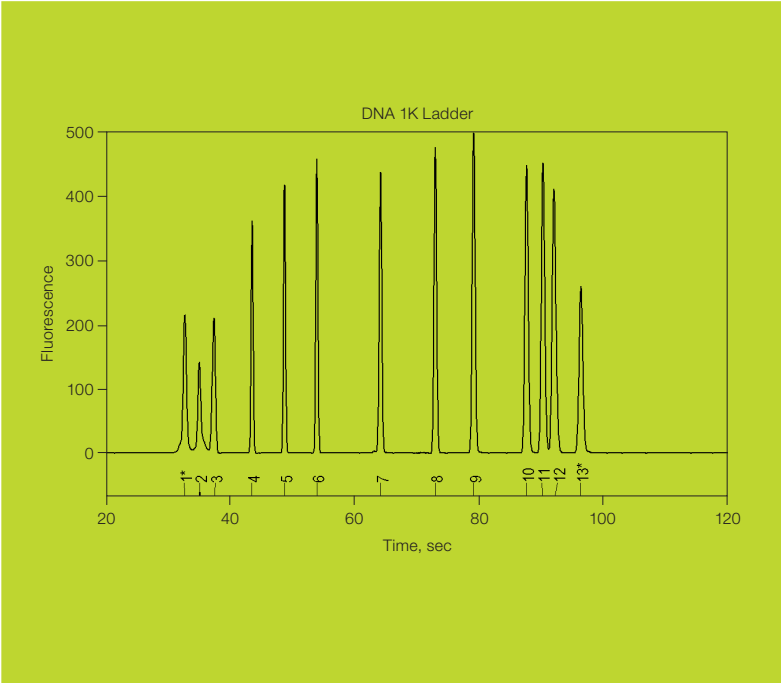
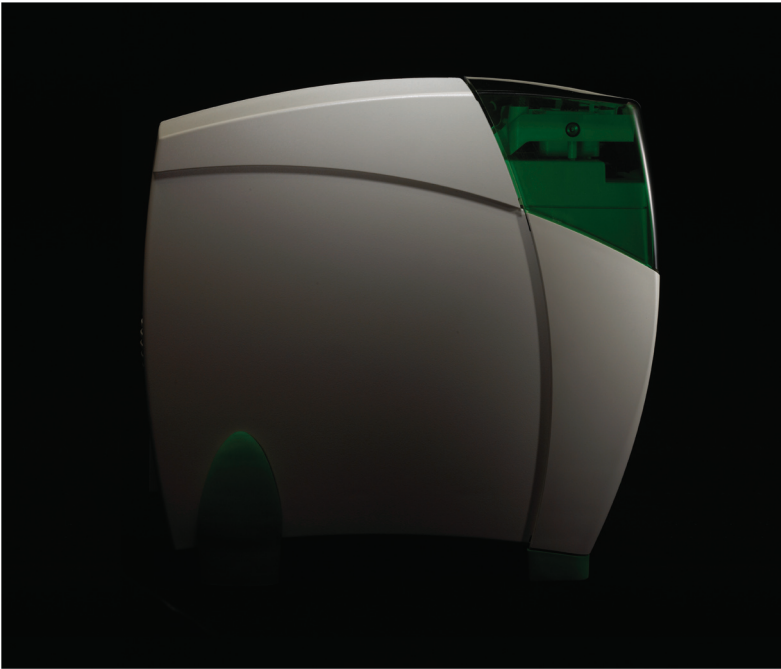
Ordering Information

Catalog #	Description
700-7001	Experion System, 100–240 V , for RNA and DNA analyses, includes electrophoresis station, priming station, vortex station, software, USB2 cable
700-7107	Experion DNA 1K Analysis Kit for 10 Chips , includes 10 DNA chips, 1 cleaning chip, 3 x 250 µl DNA 1K gel, 40 µl DNA 1K stain, 20 µl DNA 1K ladder, 750 µl DNA 1K loading buffer, 3 spin filters
700-7163	Experion DNA Chips , 10, for DNA 1K and 12K analyses, plus 1 cleaning chip
700-7164	Experion DNA 1K Reagents and Supplies for 10 Chips , includes 3 x 250 µl DNA 1K gel, 40 µl DNA 1K stain, 20 µl DNA 1K ladder, 750 µl DNA 1K loading buffer, 3 spin filters
700-7253	Experion DEPC-Treated Water , 100 ml
700-7261	Experion DNA 1K Ladder , 20 µl
700-7251	Experion Cleaning Chips , 10
700-7254	Experion Spin Filters , 10
700-7264	Cleaning Swabs , lint-free, for Experion system electrode deep cleaning, 25
700-7252	Experion Electrode Cleaner , 250 ml
732-6300	PCR Kleen Spin Columns , 25
732-6223	Micro Bio-Spin 30 Columns , includes 25 columns in Tris buffer, 50 collection tubes



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

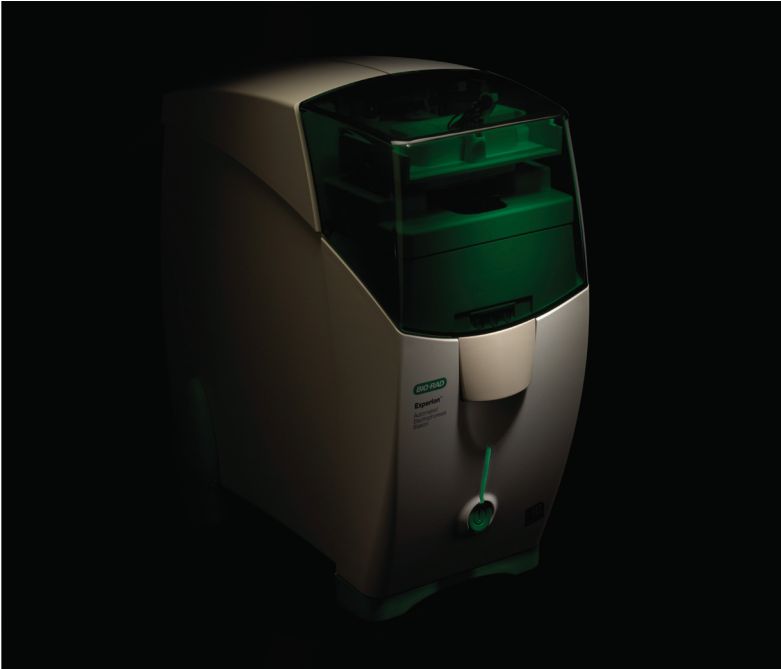
Bio-Rad Laboratories, Inc.

Life Science Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 31 3689 6600 **Canada** 905 364 3435 **China** 86 20 8732 2339 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00 **France** 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 777 4396 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **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™ DNA 1K Analysis Kit
Quick Guide

BIO-RAD

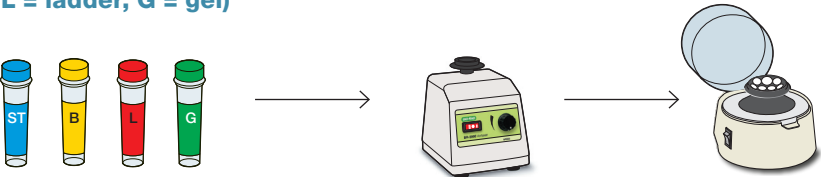


Experion DNA 1K Analysis Kit Quick Guide

For complete instructions, refer to the Experion DNA analysis kit instruction manual. Full manuals are available online at www.bio-rad.com. You can contact us by phone at 1-800-424-6723 or call your local Bio-Rad office for an electronic copy. Read the full protocol and essential practices sections if using for the first time.

1

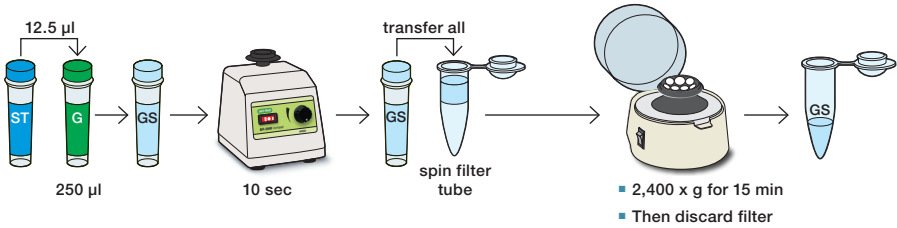
Equilibrate Kit Reagents (ST = stain, B = loading buffer, L = ladder, G = gel)



- Equilibrate reagents to room temperature (~15 min). Protect them from light
- Briefly vortex and spin down reagents

2

Prepare Gel-Stain (GS) Solution



Note: After use, store filtered GS (protected from light) at 4°C. Discard any remaining filtered GS after 1 month. Do not refilter.

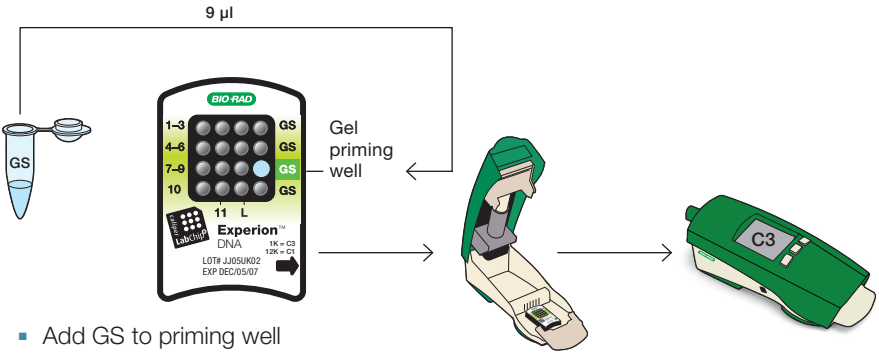
3

Prepare DNA Samples (If Needed)

- If samples are very concentrated, dilute (with TE buffer or DNase-free water) to approximately 50 ng/µl. Buffer exchange or clean up samples (for example, using PCR clean up kits) if samples are in salt concentrations ≥ 200 mM KCl (or NaCl) and 15 mM $MgCl_2$

4

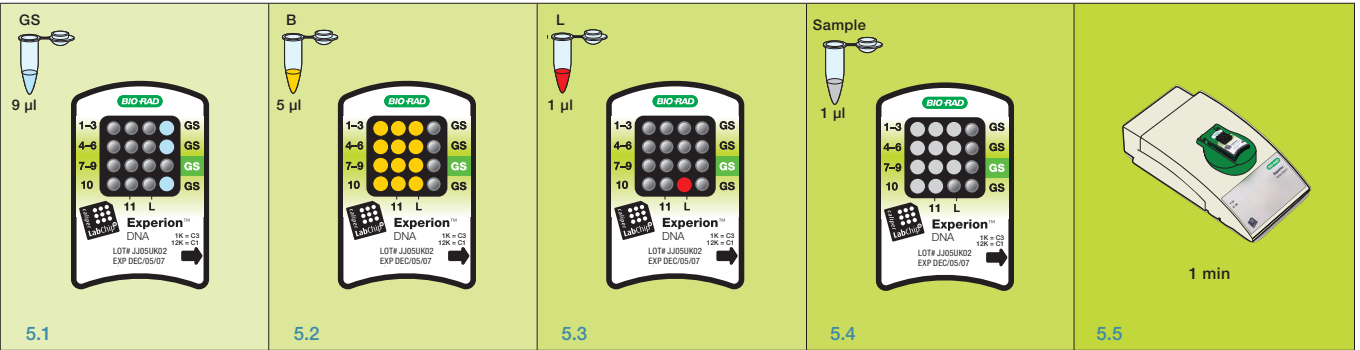
Prime the Chip



- Add GS to priming well
- Select C3 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

5

Load the Prepared Samples and DNA 1K Ladder Onto the Chip



- Pipet 9 µl filtered GS solution into other 3 **GS** wells
- Pipet 5 µl loading buffer into the **L** well and each sample well (1–11); do not leave any well empty
- Pipet 1 µl DNA 1K ladder into well **L**
- Pipet 1 µl of DNA sample into each of the 11 sample wells. Pipet 1 µl of TE or DNase-free water into any unused wells
- Place the chip in the Experion vortex station and vortex (press Mix; vortex station automatically stops after 1 min)
- Run the chip in the Experion electrophoresis station within 5 min of loading

6

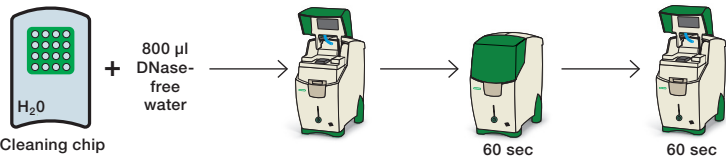
Run the DNA 1K Analysis



- Select **New Run**, then **DNA 1K** assay
- Click the **Start** (▶) button onscreen
- Select number of samples to run and start
- When the run is complete, remove and discard the used chip

7

Clean the Electrodes



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