

AUTOMATED ELECTROPHORESIS

Experion™ DNA 1K and 12K Analysis Kits

- For automated electrophoresis detection and basic analysis with the Experion system
- Accurate sizing and quantitation of DNA fragments
- Analysis of 1–11 samples in 30 to 40 minutes
- Innovative Caliper Life Sciences' LabChip technology



Single-Step DNA Analysis in 40 Minutes

Introduction

Experion DNA 1K and DNA 12K analysis kits combine state-of-the-art chip design with high-quality reagents to perform reproducible, quantitative, and accurate analysis of DNA fragments in minutes with the Experion system. Streamlined chip preparation methods and minimal sample requirements result in rapid experiments with minimal hands-on time.

The Experion DNA 1K and DNA 12K Advantage

Innovative chip design, chemistry, software, and hardware make DNA sizing and quantitation fast and easy.

- Analysis of 1–11 samples in 30 to 40 minutes
- Single-step determination of DNA purity
- Accurate sizing and quantitation of DNA fragments

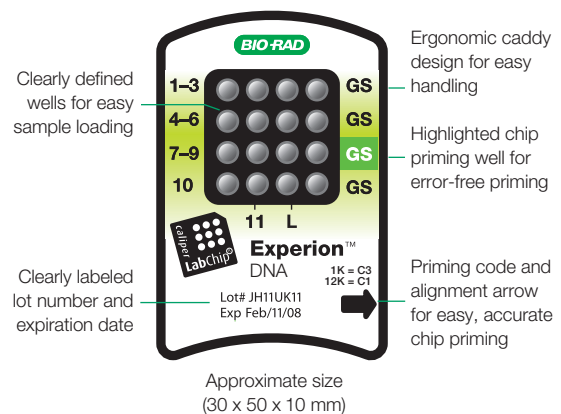
Experion DNA 1K and 12K Chips

Experion DNA 1K and DNA 12K chips offer an optimized microchannel architecture within an ergonomic caddy for exceptional DNA separation, chip handling, and results. Each chip holds up to 11 samples and is clearly labeled with well number, priming code, lot number, and expiration date for easy, error-free runs.

Experion DNA 1K and 12K Reagents

Both the DNA 1K and DNA 12K assays utilize the same chips and similar protocols, which makes it easy to switch between DNA assays. The DNA assays yield good resolution over the separation range, either between 15–1,500 bp or between 50–17,000 bp, thereby allowing accurate quantitation and sizing. Experion DNA analysis kits include the following reagents and supplies:

- High-quality gel matrix for separation comparable to agarose DNA gels
- Fluorescent dye for high sensitivity and accurate detection
- Experion DNA ladder for accurate sizing and quantitation
- Optimized loading buffer for reproducible results
- Spin filters for convenient gel preparation



BIO-RAD



Experience Meets Innovation

Specifications

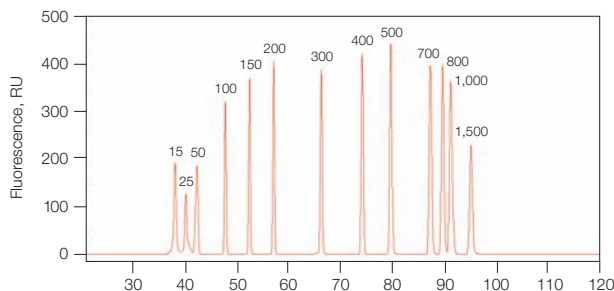
	DNA 1K Assay	DNA 12K Assay
Analytical		
Separation sizing range	15–1,500 bp	50–17,000 bp
Resolution	25–100 bp: 5 bp 100–700 bp: 5% 700–1,000 bp: 10%	100–1,500 bp: 10% 1,500–12,000 bp: 20%
Quantitation range*	0.5–50 ng/μl	0.5–50 ng/μl
Limits of detection**	0.1 ng/μl	0.1 ng/μl
Sizing accuracy***	±10%	±15%
Sizing reproducibility***	CV ≤ 5%	CV ≤ 5%
Quantitation accuracy*** (in TE buffer)	±25%	±20%
Quantitation reproducibility*** (in TE buffer)	CV ≤ 20%	CV ≤ 20%
Maximum salt concentration in samples	200 mM KCl/NaCl 15 mM MgCl ₂	250 mM KCl/NaCl 15 mM MgCl ₂
Physical		
Sample volume	1 μl	1 μl
Number of samples/chip	11	11

* 100 bp and larger.

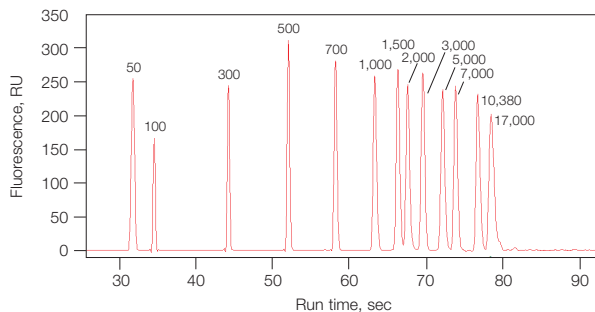
** Outside recommended quantitation range.

*** Applies to interchip and intrachip variation.

DNA 1K ladder



DNA 12K ladder



The DNA 1K and 12K ladders. The DNA fragment sizes are indicated in base pairs above the peaks.

Experion Software Requirements

The DNA 1K and DNA 12K analysis kits require Experion software version 2.1 or higher. Current version is 3.0.

Ordering Information

Catalog #	Description
700-7107	Experion DNA 1K Analysis Kit for 10 Chips, includes 10 DNA chips, Experion DNA 1K reagents and supplies for 10 chips
700-7108	Experion DNA 12K Analysis Kit for 10 Chips, includes 10 DNA chips, Experion DNA 12K reagents and supplies for 10 chips
700-7050	Experion Software, version 3.0, system operation and standard data analysis tools, PC
700-7163	Experion DNA Chips, 10
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
700-7165	Experion DNA 12K Reagents and Supplies, for 10 chips, includes 650 μl DNA 12K gel, 40 μl DNA 12K stain, 20 μl DNA 12K ladder, 750 μl DNA 12K loading buffer, 3 spin filters
700-7251	Experion Cleaning Chips, 10
700-7252	Experion Electrode Cleaner, 250 ml
700-7253	Experion DEPC-Treated Water, 100 ml
700-7261	Experion DNA 1K Ladder
700-7262	Experion DNA 12K Ladder



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

**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