



dsDNA 1000 Fluorescent Detection and Analysis Kit

Specifications

Fragment size range	100–1,000 bp
Resolution	10–20 bp
Number assays	400 runs
Capillary	BioCAP 75 μ m ID, coated, 50 cm length, 2
Dye	Fluorescent nucleic acid stain
Run buffer	Hydrophilic polymer in TBE
DNA standard	100 bp ladder
Dye excitation wavelength	488 nm
Dye emission wavelength	520 nm

Fluorescent DNA analysis.

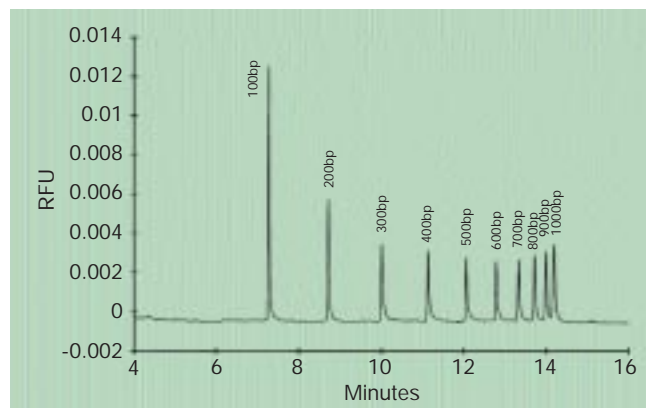
Bio-Rad now offers a completely automated DNA analysis system that's extremely fast and convenient. Our new dsDNA 1000 Fluorescent Detection and Analysis Kit is designed to complement the new BioFocus[®] capillary electrophoresis laser Induced Fluorescence LIF² Detection System. These two comprise an excellent choice for automated fluorescent analysis of DNA fragment size polymorphisms, DNA point mutations, microsatellite DNA analysis, DNA deletion analysis, and PCR[®] product analysis.

Easy labeling, high resolution.

The new dsDNA 1000 Fluorescent Detection and Analysis Kit contains all the reagents you need to quickly label and resolve double stranded DNA fragments ranging from 100 to 1,000 base pairs. The kit utilizes a special hydrophilic polymer buffer which creates a sieving effect analogous to acrylamide. The ready-to-use buffer is replenished prior to each analysis, insuring the highest reproducibility of DNA fragment migration time and peak area. A 100 base-pair DNA ladder is included so you can accurately determine fragment sizes.

Superior fluorescence sensitivity.

The BioFocus LIF² detector is at least 1,000 times more sensitive than existing UV absorbance detection. The key to laser induced fluorescence sensitivity is focusing a beam of high intensity, single wavelength laser light on a capillary. The BioFocus LIF² detector is capable of detecting picomolar concentrations of DNA, which is as sensitive as autoradiography, without the hazardous radioisotopes.



Bio-Rad 100 bp molecular ruler (170-8202) resolved with the dsDNA Fluorescent Detection and Analysis Kit.

Focus on Performance
The BioFocus[®] System

BIO-RAD



dsDNA 1000 Fluorescent Detection and Analysis Kit

Electrophoretic analysis that's completely automated.

Simply load the carousels with buffers and samples, program the run conditions, and the BioFocus system does it all – real time detection, automatic data integration, and report printing for each DNA analysis. Not only are capillary separations fast, but the resolution is outstanding.

The BioFocus system virtually eliminates the need for conventional separation techniques including gel casting, sample loading, staining, exposing film, and film imaging.

BioFocus LIF² is the most advanced capillary electrophoresis system. Let the BioFocus LIF² system work for you! Contact your local Bio-Rad representative today.

* The Polymerase Chain Reaction (PCR) Process is covered by patents owned by Hoffman-LaRoche. Use of the PCR process requires a license.

Ordering Information

Catalog No.

148-4132

Product Description

dsDNA 1000 Fluorescent Detection and Analysis Kit, includes 2 capillaries, run buffer, DNA ladder, fluorescent nucleic acid stain

148-5100

dsDNA 1000 Fluorescent Nucleic Acid Stain

Focus on Performance
The BioFocus[®] System

BIO-RAD

**Bio-Rad
Laboratories**

Life Science
Group

U.S. (200) 620-8740 • Canada (604) 761-1000 • Australia 08 2214 2200 • Austria (1) 277 2501 • Belgium 09 226 66 11 • Canada (905) 713 2771 • China (20) 01 2049 922 • Denmark 22 17 2247 • Finland 20 246 2207 • France (1) 62 20 49 20 • Germany 029 2 62 35 0 • India 21 11 49 0 02 • Italy 02 2 602 1 • Japan 03 62 11 62 0 • Hong Kong 22 222 00 • The Netherlands 020 2 62 40 99 • New Zealand 09 442 2222 • Singapore (65) 273 2227 • Spain (91) 661 70 25 • Sweden 08 01 2 62 7 0 00 • Switzerland 01 2 62 65 66 • United Kingdom 020 12 11 35