



## Capillary Electrophoresis Oligonucleotide Analysis Kit

### Determining the purity of synthetic oligonucleotides has never been so fast, easy, and reliable.

Before using oligonucleotides in most molecular biology experiments, it's important to check their quality. But to perform a routine analysis with HPLC or gel electrophoresis, you need large volumes of reagents, as well as plenty of spare time.

However, capillary electrophoresis makes purity screening of synthetic DNA much faster. And Bio-Rad's Capillary Electrophoresis Oligonucleotide Analysis Kit makes analysis of synthetic oligonucleotides simple and reliable.

### One complete kit.

The Oligonucleotide Analysis Kit from Bio-Rad contains standards, analysis buffer, two lengths of BioCAP™ oligonucleotide analysis capillaries, and detailed instructions – everything you need for purity analysis.

### Unique polymer-coated capillary.

BioCAP coated capillaries feature a patented internal coating that's covalently bound to the capillary wall. Because this coating virtually eliminates electro-osmotic flow in the capillary, you get rapid and reproducible high resolution separations. In addition, the high stability of the coating provides excellent capillary lifetime.

### Fast, automatic results.

You want fast, efficient automation to maximize productivity. And that's just what the BioFocus® capillary electrophoresis system has to offer. Once you've loaded the samples into the carousel, our system automatically analyzes each sample, and performs real time data collection, data integration, and report generation.

### Accurate, reproducible oligonucleotide analysis.

The BioFocus CE oligonucleotide analysis kit features a hydrophilic polymer sieving buffer that provides single-base resolution of oligonucleotides in the 4 to 40 mer range. The polymer containing buffer is replenished prior to each new run, providing high reproducibility of migration time and peak area.

Focus on Performance  
The BioFocus® System

**BIO-RAD**

# Information Sheet

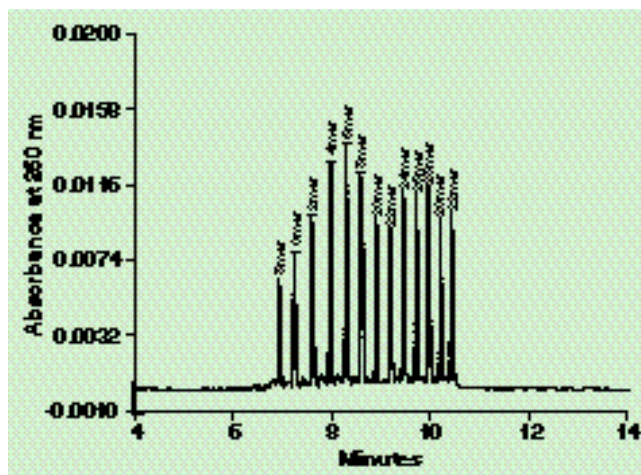


## Capillary Electrophoresis Oligonucleotide Analysis Kit

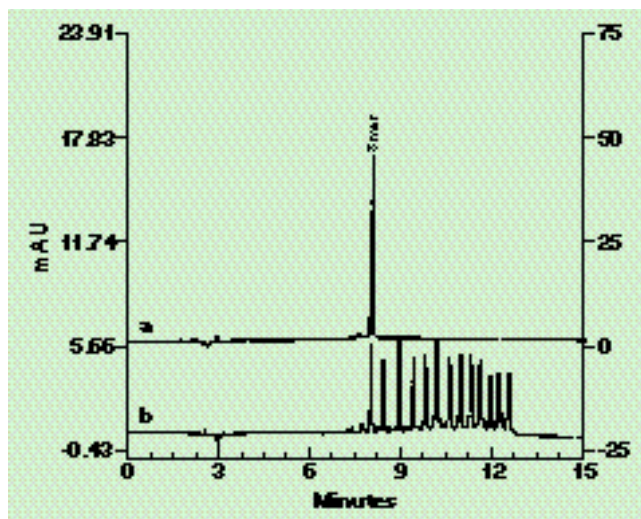
### Ordering Information

Catalog No. Product Description  
148-4140 CE Oligonucleotide Analysis Kit, includes 2 capillaries, run buffer, oligonucleotide standard, and protocol for use on BioFocus systems

148-5026 CE Oligonucleotide Buffer



Electropherogram of the CE oligonucleotide size standards.



Overlay of 8 mer sample (a) and the Oligonucleotide size standard (b).

Focus on Performance  
The BioFocus® System



Bio-Rad  
Laboratories

Life Science  
Group

U.S. (800) 4BIORAD • California (510) 741-1000 • Australia 02-9914-2800 • Austria (1)-877 89 01 • Belgium 09-385 55 11 • Canada (905) 712-2771 • China (86-10) 2046622 • Denmark 39 17 9947 • Finland 90 804 2200 • France (1) 43 90 46 90 • Germany 089 318 84-0 • India 91-11-461-0103 • Italy 02-21609 1 • Japan 03-5811-6270 • Hong Kong 7893300 • The Netherlands 031318-540666 • New Zealand 09-443 3099 • Singapore (65) 272-9877 • Spain (91) 661 70 85 • Sweden 46 (0) 8 627 50 00 • Switzerland 01-809 55 55 • United Kingdom 0800 181134