



BioFocus® Verification Kit

Keep it in check.

Here's a simple way to check the performance of your BioFocus capillary electrophoresis system – the BioFocus verification kit. It contains everything you need to quickly and accurately run a performance diagnosis of important BioFocus parameters:

- High voltage power supply
- Temperature control
- Detector
- Injection pneumatics

The BioFocus system will automatically perform the verification with software versions 5.2 or higher. For earlier versions, you can simply program the system to perform these functions.

Simple chemistry.

Benzene sulfonic acid was selected as the test sample because it is chemically stable, offers minimal interaction with the capillary, and under test conditions, has proven to be exceptionally reproducible. The acidic run buffer eliminates any effects of electro-osmotic flow in the capillary. And this kit takes advantage of our BioCAP™ fused silica capillary which provides a consistent test environment for evaluating important system parameters.

Simple procedure.

After you've loaded your sample and buffer, simply select the automatic sequence 'verify', and the BioFocus system automatically performs ten test runs, storing all data. To analyze the data, select 'verification', and the system automatically analyzes migration time and peak area for each run. The relative standard deviation of the migration times and peak areas are calculated, and the system will either 'pass' or 'fail'. A detailed verification report is automatically generated.

Focus on Performance
The BioFocus® System

BIO-RAD

Information Sheet



BioFocus® Verification Kit

Bio-Rad Laboratories BioFocus Integrator v 5.20

BioFocus System VERIFICATION REPORT

Report Date Mon Jan 20 14:34:17 1997
Instrument Model BioFocus 3000TC
Instrument Serial # 951BR0038
Instrument Hardware ID 0101
Z180 Firmware Rev. J
HC11 Firmware Rev. F
Operating Software Version 5.20

Operator Name Rowena Ng
Institution CE Chemistry R&D
Life Science Group-
Bio-Rad Labs, Hercules, CA

Comments CE Applications Laboratory

OPERATING PARAMETERS

Capillary 24 cm x 50 µm uncoated
Capillary Temp 22 deg C
Sample Verification Test Sample
Inlet 148-5010 phosphate, 0.1M
Outlet 148-5010 phosphate, 0.1M
Prep Cycle 1 Pre-Inject: Inlet, PM_PREP; Outlet,
Waste; 30 sec HI Pressure
Injection 8 psi*sec
Run Voltage 8.00 kV - to +
Detector Single wavelength at 200 nm
Carousel Temp not set

TEST DATA

File Name	Migration Time	Peak Area	Date/Time
VER_001	4.06	1332918	Mon Jan 20 12:56:46 1997
VER_002	3.96	1306927	Mon Jan 20 13:04:50 1997
VER_003	3.87	1310403	Mon Jan 20 13:12:56 1997
VER_004	4.01	1353899	Mon Jan 20 13:21:00 1997
VER_005	4.01	1359515	Mon Jan 20 13:29:04 1997
VER_006	4.02	1361984	Mon Jan 20 13:37:10 1997
VER_007	3.99	1357714	Mon Jan 20 13:45:16 1997
VER_008	4.02	1363959	Mon Jan 20 13:53:20 1997
VER_009	4.00	1367551	Mon Jan 20 14:01:26 1997
VER_010	4.00	1366857	Mon Jan 20 14:09:30 1997

CALCULATED MEAN 4.003 1348162
CALCULATED STD DEV 0.028 23059
CALCULATED RSD 0.70% 1.71%
ACCEPTABLE RSD 1.50% 2.50%

RESULTS PASS PASS

Signature: _____ Date: _____

A detailed verification report is automatically printed.

Ordering Information

Catalog No. Product Description

148-4170 **BioFocus System Verification Kit**, includes 2 uncoated capillaries (40 cm x 50 µm ID x 375 µm OD, window at 8 cm), run buffer, test sample, and instructions for use on BioFocus systems

Buffers and Accessories

148-5010 **Phosphate Buffer**, 01 M, pH 2.5, 4 x 60 ml
148-5011 **Phosphate Buffer**, 0.1 M, pH 2.5, 250 ml
148-3060 **BioCAP Bare Silica Capillary**, 50 µm ID x 375 µm OD x 40 cm, window at 8 cm
148-3053 **Capillary Cartridge**, 24 cm x 50 µm, uncoated, capillary installed

Focus on Performance
The BioFocus® System

BIO-RAD

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