



**BIO-RAD**

## Declaration of Conformity

Bio-Rad Laboratories, Inc., 1000 Alfred Nobel Drive, Hercules, California, 94547, U. S. A., declares that the product

Econo Gradient Pump (731-9002)  
SPL3-2 Splitter Valve (731-9011)  
SV-3 Diverter/Bypass Valve (731-8322)  
Gradient Mixer (731-8323)

to which this declaration relates, is in conformity to the following standards or normative documents.

EN50082-1: 1997

EN61000-3-2: 1995 Class A

EN61000-3-3: 1995

EN61000-4-2: 1995

EN61000-4-3: 1996

EN61000-4-4: 1995

EN61000-4-5: 1995

EN61000-4-6: 1996

EN61000-4-8: 1994

EN61000-4-11: 1994

ENV50204: 1995

EN50081-2: 1994

EN55011: 1991 Class A

EN61010-1: 1993 + A2

IEC61010-1: 1990 + A1 + A2

following the provisions of the 89/336/EEC & 73/23/EEC Directive.

The product is imported into the EU by Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead (London area), Hertfordshire HP2 7TD England.

July 12, 1999  
date of issue

Norman Schwartz  
Vice President