



**BIO-RAD**

## Declaration of Conformity

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

iCycler Thermal Cycler  
iCycler iQ Module

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

EN55011: 1994 with A1:1995 Class A Group 1

EN50082-1: 1997

EN61000-4-2: 1995

EN61000-4-3: 1997

ENV50204: 1995

EN61000-4-4: 1995

EN61000-4-5: 1995

EN61000-4-6: 1996

EN61000-4-8: 1994

EN61000-4-11: 1994

EN61010-1: 1993 + 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.

October 18, 1999

date of issue

Norman Schwartz  
Vice President