

<u>┇╱╱╱╱┧╒╱╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙</u>┋╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙┋╱┈╙

Declaration of Conformity

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

iCycler Thermal Cycler

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.

June 30,1999 date of issue

Norman Schwartz Vice President

C1708701 Rev A