

BIO-RAD

Declaration of Conformity

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

Buffer Recirculation Pump 220-240 V

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

EN55011:1991
EN50082-2:1994
EN61000-4-2:1995
ENV50141:1994
ENV50140:1994
EN61000-4-4:1995
EN61000-4-8:1995
EN61010-1:1993

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.

29 November, 1995
date of issue

Burton A. Zabin

Burton Zabin
Vice President

DC1702930 Rev A