



BIO-RAD

Declaration of Conformity

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

BioLogic LP System (731-8301)
BioLogic LP Controller (731-8349)
BioLogic LP Buffer Select Valve (731-8321)
BioLogic LP Div/Bypass Valve (731-8322)
BioLogic LP Mixer (731-8323)
BioLogic LP Optics Module (731-8324)
Model EG-1 Flow Cell (731-8155)

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

EN55011:1991 Class A
EN50082-2:1994
EN61000-4-2:1995
EN61000-4-4:1995
EN61000-4-8:1995
ENV54140:1994
ENV54141:1994
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.

22 March, 1996
date of issue



Burton Zabin
Vice President

DC7318301 Rev A