

BIO-RAD

Declaration of Conformity

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

BioLogic QuadTec (7601305)

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

Low Voltage Ordinance (73/23/EWG); EN 61010-1; 1993

Engineering Guidelines (89/392/EWG)

EMV Ordinance (89/336/EWG)

EN 50081-1; 1992

EN 55011; 1991 Class B

EN 55022; 1987 Class B

EN 50082-1; 1992

IEC 801-1; 1984

includes IEC 41 B (sec) 81; 1992

IEC 801-3; 1984

IEC 801-4; 1988

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 6, 2000
date of issue

Norman Schwartz
Vice President