



BIO-RAD

Declaration of Conformity

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

Western Processor

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

EN61010-1: 1993

EN55014: 1993

EN55014: 1995

EN61000-4-2: 1995

EN61000-4-4: 1995

EN61000-4-5: 1995

EN61000-4-6: 1996

EN61000-4-11: 1994

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.

April 15, 1998
date of issue

CONTROLLED NETWORK COPY
This Document is Approved
for Tooling, Production and
General Distribution

Norman Schwartz
Vice President