



BIO-RAD

Declaration of Conformity

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

PROTEAN® Plus Cell

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

EN61010-1

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.

8 September, 2000
date of issue

Norman Schwartz
Vice President

DC1654150 Rev A