

PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

MHS-15 MERCURY HYDRIDE SYSTEM SERIAL NO.

This is to certify that this PerkinElmer product was tested and verified to be in conformance with all applicable quality requirements, including specifications, drawings, calibration, preservation, packing, marking requirements and part identification.

Declaration of Safety Code Compliance

This PerkinElmer product conforms to the applicable clauses stipulated in the CE Mark requirements for the LVD Directive (73/23/EEC and 93/68/EEC):

EN 61010-1:1993, Safety Requirements for Electrical Equipment for Laboratory Use
IEC 1010-1:1990 + A1:1992 + A2:1995, Safety Requirements for Electrical Equipment
for Laboratory Use
IEC 61010-1, 2nd Edition (draft), Safety Requirements for Electrical Equipment for Laboratory Use
EN 61010-2-061:1996, Safety Requirements for Laboratory Spectrometers with Thermal Atomization and Ionization



NOTE: The operation of certain types of equipment (e.g., signal generators) may be subject to given restrictions. Please refer to the appropriate information in the respective user documentation.

Declaration of System Validation

The product was found to meet its functional and performance specification prior to shipment. To support this declaration, the following Engineering, Assembly and Test documents are held by PerkinElmer and are available for reference upon request in justified cases and to an appropriate extent:

The Product Description	The System Design Documentation
The Functional Specification	The Source Code Documentation
The User Interface Definition	The Evaluation Documentation

NOTE: PerkinElmer will maintain possession of all documents and controls their reproduction, including parts of them.

The existence of these documents and the procedures used in their production are formal requirements of the PerkinElmer Quality System. The integrity of the PerkinElmer Quality System is routinely audited and is certified by the British Standards Institution as meeting the requirements of ISO 9001, the internationally recognized standard for Quality Assurance.

A handwritten signature in black ink, appearing to read 'Zhang Fu', located below the text of the signature.

Zhang Fu

as

Manager, Quality and Compliance (Singapore)