

PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

**MODEL 2400 SERIES II
CHNS/O ANALYZER**

SERIAL NO. 241N

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 EMC & Safety Code Compliance

This PerkinElmer product conforms to the regulations stipulated in the CE Mark requirements for the EMC Directive (2004/108/EC) and the LVD Directive (2006/95/EC):

EN 55011:1998 Group 1, Class B, EMC -- RF Characteristics of ISM Equipment
EN 61326:1997 + A1:1998 + A2:2001, EMC -- Requirements for Electrical Equipment for Laboratory Use
IEC 61000-4-2:1995, EMC -- Electrostatic Discharge Requirements
IEC 61000-4-3:1995, EMC -- Radiated Electromagnetic Field Requirements
IEC 61000-4-4:1995, EMC -- Electrical Fast Transient/Burst Requirements
IEC 61000-4-5:1995, EMC -- Surge Immunity Requirements
IEC 61000-4-6:1996, EMC -- Conducted Disturbances (induced by RF fields) Requirements
IEC 61000-4-11:1994, EMC -- Voltage Dips, Short Interruptions, Voltage Variations Requirements
EN 61000-3-2:2000, EMC -- Harmonic Current Emissions
EN 61000-3-3:1995 + A1:2001, EMC -- Voltage Fluctuations and Flicker
EN 61010-1:2001, Safety Requirements for Electrical Equipment for Laboratory Use



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 Management System. The integrity of the PerkinElmer Quality Management System is routinely audited and has been certified to ISO 9001 since 1992.

John C. Bouse
Staff Engineer – Compliance
R&D Shared Services