

PRODUCT CERTIFICATION AND DECLARATION OF CONFORMITY

SIMULTANEOUS THERMAL ANALYZER SERIAL NO. (STA) 6000

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 55022:1998 + A1:2000 + A2:2003 Group 1, Class B (radiated), EMC -- RF Characteristics of ITE Equipment
EN 55022:2006 Group 1, Class B (conducted), EMC -- RF Characteristics of ITE Equipment
EN 61326:2006, EMC -- Requirements for Electrical Equipment for Laboratory Use
EN 61000-4-2:1995 + A1:1998 + A2:2001, EMC -- Electrostatic Discharge Requirements
EN 61000-4-3:2006, EMC -- Radiated Electromagnetic Field Requirements
ENV 50204:1995, EMC -- Radiated Electromagnetic Field from Digital Radio Telephones Requirements
EN 61000-4-4:2005, EMC -- Electrical Fast Transient/Burst Requirements
EN 61000-4-5:2007, EMC -- Surge Immunity Requirements
EN 61000-4-6:2007, EMC -- Conducted Disturbances (induced by RF fields) Requirements
EN 61000-4-8:1994 + A1:2001, EMC -- Power Frequency Magnetic Field Immunity Requirements
EN 61000-4-11:2004, EMC -- Voltage Dips, Short Interruptions, Voltage Variations Requirements
EN 61000-3-2:2000 + A2:2005, 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
EN 61010-2-010:2003, Safety Requirements for Laboratory Equipment for the Heating of Materials
EN 61000-2-081:2002, Particular Requirements for Automatic and Semi-automatic Laboratory Equipment for Analysis and Other Purposes



FCC Part 15, Class A, Radiated and Conducted Emissions

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.

John C. Bouse
Staff Engineer – Compliance
R&D Electrical Engineering