

Declaration of Conformity

STA 8000 Simultaneous Thermal Analyzer **Serial No:**

We, PerkinElmer, declare under our sole responsibility that the product to which this declaration relates is in conformity with the following European Directives:

EMC Directive 2004/108/EC
Low Voltage Directive 2006/95/EC
RoHS Directive 2011/65/EU

Standards to which conformity is declared:

EMC:	EN 61326-1:2006	Electrical equipment for measurement, control and laboratory use - EMC requirements. Including the underlying standards;
	EN 61000-4-2	
	EN 61000-4-3	
	EN 61000-4-4	
	EN 61000-4-5	
	EN 61000-4-6	
	EN 61000-4-11	
	EN 55011 Class A	
	EN 61000-3-2	
	EN 61000-3-3	
LVD:	EN 61010-1:2010	Safety Requirements for Electrical Equipment for Measurement, Control and Laboratory Use.
	EN 61010-2-010:2003	

CE marking affixed in 2012

The integrity of the PerkinElmer Quality System is routinely audited and is certified as meeting the requirements of ISO 9001, the internationally recognized standard for Quality Assurance.



J. Dadswell
Quality & Compliance Manager R&D
PerkinElmer Ltd. UK

Issue date:
June 2013