

## Declaration of Conformity

**TMA 4000 Thermomechanical Analyzer**

**Serial No:**

We, PerkinElmer, declare 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**

Standards to which conformity is declared:

**EMC:**        **EN 55011:2009**                      Industrial, scientific and medical (ISM) radio-frequency equipment -  
Electromagnetic disturbance characteristics - Limits and methods of  
measurement.

**EN 55011:2009**                      Conducted and Radiated Emissions ClassA  
**EN 61000-4-2**                      Electrostatic Discharge  
**EN 61000-4-3**                      Radiated Susceptibility  
**EN 61000-4-4**                      Transient Susceptibility  
**EN 61000-4-5**                      Surge Susceptibility  
**EN 61000-4-6**                      Conducted Immunity  
**EN 61000-4-8**                      Magnetic Immunity  
**EN 61000-4-11**                      Voltage Fluctuations  
**EN 61000-3-2**                      Harmonics  
**EN 61000-3-3**                      Flicker

### **CFR47 FCC Part 15, SubPart B, Class B**

**LVD:**        **IEC 61010-1:2010**                      Safety Requirements for Electrical Equipment for Measurement,  
Control and Laboratory Use.  
**IEC 61010-2-010:2005**                      Particular requirements for laboratory equipment for the heating of  
materials  
**IEC 61010-2-081:2009**                      Particular requirements for automatic and semi-automatic laboratory  
equipment for analysis and other purposes

### **CE marking affixed in 2013**

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:**  
**December 2013**