

DECLARATION OF CONFORMITY



The manufacturer **PerkinElmer Singapore Pte Ltd**

hereby declares that the **Laboratory Equipment (Class 8721-06 & 8721-86)**

of equipment models **MicroBeta2
2450-0010, 2450-0020, 2450-0060, 2450-0120,
2460-0010, 2460-0020, 2460-0060, 2460-0120**

specifically **Model No.: _____ Serial No.: _____ Mfg Date.: _____**

*conforms to the
and its standards* **EMC Directive (2014/30/EU)
EN 61326-1:2013 Class A
EN 61000-3-2:2006 +A1: 2009 +A2 :2009, EN 61000-3-3:2008,
EN 55011:2009 +A1:2010
IEC 61000-4-2:2008, IEC 61000-4-3:2010, IEC 61000-4-4:2004 +A1:2010
IEC 61000-4-5:2005, IEC 61000-4-6:2008, IEC 61000-4-8:2009,
IEC 61000-4-11:2004**

as well as the **Low Voltage Safety Directive (2014/35/EU)**

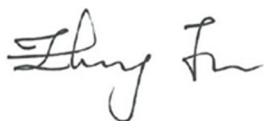
and its standards **IEC 61010-1:2010 (3rd Ed); CAN/CSA C22.2 61010-1-12, UL 61010-1(3rd ED)
Canada & US National Differences**

and most recently, the **Machinery Directive (2006/42/EC), described in Risk Assessment: MB2-RA,
05/12**

and also the **Restriction of Hazardous Substances in Electrical and Electronic
Equipment Directive 2011/65/EU**

CE Marking Affixed in 2013

This declaration is issued under the responsibility of



Zhang Fu
as
Senior Manager Quality Assurance

PerkinElmer Document: 8845401, Rev F