
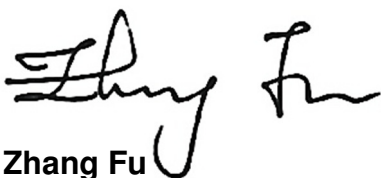


## DECLARATION OF CONFORMITY

<i>The manufacturer</i>	<b>PerkinElmer Singapore Pte Ltd</b>	 Since 2013
<i>hereby declares that the</i>	<b>Laboratory Equipment (Class 8721-06 &amp; 8721-86)</b>	
<i>of equipment models</i>	<b>MicroBeta2 2450-0010, 2450-0020, 2450-0060, 2450-0120, 2460-0010, 2460-0020, 2460-0060, 2460-0120</b>	
<i>specifically</i>	<b>Model No.:</b> _____ <b>Serial No.:</b> _____ <b>Mfg Date.:</b> _____	
<i>conforms to the and its standards</i>	<b>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</b>	
<i>as well as the</i>	<b>Low Voltage Safety Directive (2014/35/EU)</b>	
<i>and its standards</i>	<b>IEC 61010-1:2010 (3<sup>rd</sup> Ed); CAN/CSA C22.2 61010-1-12, UL 61010-1(3<sup>rd</sup> ED) Canada &amp; US National Differences</b>	
<i>and most recently, the</i>	<b>Machinery Directive (2006/42/EC), described in Risk Assessment: MB2-RA, 05/12</b>	
<i>and also the</i>	<b>Restriction of Hazardous Substances in Electrical and Electronic Equipment Directive 2011/65/EU</b>	

*This declaration is issued under the responsibility of*



Zhang Fu  
as  
Senior Manager Quality Assurance