

## DECLARATION OF CONFORMITY



Since  
2013

*The manufacturer* **PerkinElmer Singapore Pte Ltd**

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

*of equipment models* **WIZARD2**  
**2470-0010, 2470-0020, 2470-0050, 2470-0100, 2470-0150, 2470-0200,**  
**2480-0010, 3470-0050, 3470-0100, 3470-0150, 3470-0200**

*specifically* **Model No.:** \_\_\_\_\_ **Serial No.:** \_\_\_\_\_ **Mfg Date:** \_\_\_\_\_

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

*as well as the and its standards* **Low Voltage Safety Directive (2014/35/EU),**  
EN 61010-1:2010  
EN 61010-2-101:2017

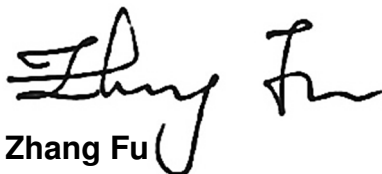
IEC 61010-1:2010; IEC 61010-2-101:2015; CAN/CSA- C22.2 61010-1-12; CAN/CSA-  
C22.2 61010-2-101-15; UL 61010-1 (3<sup>rd</sup> Ed)

*and the* **In Vitro Diagnostic Directive IVDD 98/79/EC, Annex III**  
Other device (all devices except Annex II and self-testing devices)

*and also the* **Restriction of Hazardous Substances in Electrical and Electronic Equipment Directive 2011/65/EU**  
Beside an application exempted from the restriction in Article 4(1)  
Specific to medical devices by Annex IV of RoHS Directive as below:  
5. Lead in shielding for ionizing radiation.

**Authorized Representative**  
EMERGO EUROPE  
Prinsessegracht 20  
2514 AP The Hague  
The Netherlands

***This declaration is issued under the responsibility of***



**Zhang Fu**  
as  
**Senior Manager Quality Assurance**