

# DECLARATION OF CONFORMITY

---

**The manufacturer located in** **PerkinElmer Singapore Pte Ltd**  
2 Tukang Innovation Grove  
#04-01 JTC MedTech Hub @ MedTech Park  
Singapore 618305

**hereby declares under sole responsibility that the product**  
Instrument: FIMS 100  
Serial Number:

**conforms to the and its standards** EMC Directive (2014/30/EU)  
EN 61326-1:2013  
EN 61326-2-2:2013  
IEC 61326-1:2012  
IEC 61326-2-2:2012

**as well as the and its standards** Low Voltage Directive (2014/35/EU)  
EN 61010-1:2010  
EN 61010-2-081:2015  
IEC 61010-1:2010  
IEC 61010-2-081:2015

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



*This declaration is issued under the responsibility of*



Zhang Fu  
**as**  
Manager, Quality and Compliance (Singapore)

2 Tukang Innovation Grove #04-01 JTC MedTech Hub @ MedTech Park Singapore 618305  
[in the EU: PerkinElmer, Chalfont Road, Buckinghamshire, Seer Green, HP9 2FX, United Kingdom]