

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: S10 AUTOSAMPLER
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)

