

Manufacturer Sartorius Liquid Handling oy
Tulppatie 1, 00880 Helsinki, Finland

We hereby declare under our sole responsibility that the following products

Model & Name

710969	Rline
710993	Rline
710995	Rline
LH-710969KIT	Rline Development Kit
LH-710993KIT	Rline Development Kit
LH-710995KIT	Rline Development Kit

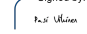
Device type Electronic laboratory equipment

in the form as delivered fulfil all the relevant provisions of the following British Regulations and meet the applicable requirements of the British Designated Standards including any amendments valid at the time this declaration was signed listed below:

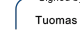
The Electromagnetic Compatibility Regulations 2016
UK Statutory Instruments 2016 No. 1091
BS EN 61326-1:2013
BS EN 61326-1:2021

The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012 (RoHS)
UK Statutory Instruments 2012 No. 3032
BS EN 50581:2012, BS EN 61010-1:2010

Sartorius Liquid Handling oy
Helsinki, 14.10.2024

Signed by:

Signer Name: Pasi Vihinen
Signing Reason: I approve this document
Signing Time: 2024-10-16 | 08:43:52 CEST
B63019D13BB549149E052159F2E6C344

Pasi Vihinen
Manager of Sustaining
Engineering, EMEA

Signed by:

Tuomas Huhmarniemi
Signer Name: Tuomas Huhmarniemi
Signing Reason: Hyväksyn tämän asiakirjan
Signing Time: 2024-10-15 | 16:16:17 CEST
CC6DF5A424E6442D9507ABA81B686FAC

Tuomas Huhmarniemi
Head of Quality