

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

We hereby declare under our sole responsibility that the following product

Name Midi Plus

Model(s) 710930
710931

Device type Electronic laboratory equipment

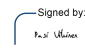
in the form as delivered fulfils all the relevant provisions of the following British Regulations and meets the applicable requirements of the British Designated Standards including any amendments valid at the time this declaration was signed listed below:


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

Accessories AC Adapter 9V 5, 5mm 800mA Universal LH-731608

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 | 08:42:41 CEST
B63019D13BBS49149E052159F2E6C344

Signed by:

Signer Name: Tuomas Huhmarniemi
Signing Reason: Hyväksyn tämän asiakirjan
Signing Time: 2024-10-15 | 16:14:25 CEST
CC6DF5A42E6442D9507ABA61B686FAC

Pasi Vihinen
Manager of Sustaining
Engineering, EMEA

Tuomas Huhmarniemi
Head of Quality