



Original

SARTORIUS

EU/EC Declaration of Conformity

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

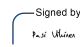
in the form as delivered fulfils all the relevant provisions of the following European Directives - including any amendments valid at the time this declaration was signed - and meets the applicable requirements of the harmonized European Standards including any amendments valid at the time this declaration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


2014/30/EU Directive on Electromagnetic Compatibility
EN 61326-1:2013
EN 61326-1:2021

2011/65/EU Restriction of the use of certain hazardous substances in electrical and electronic equipment (RoHS) (incl. (EU) 2015/863)
EN IEC 63000:2018

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 | 08:42:54 CEST
B63019D13BB549149E052159F2E6C344

Signed by:

Signer Name: Tuomas Huhmarniemi
Signing Reason: Hyväksyn tämän asiakirjan
Signing Time: 2024-10-15 | 16:14:43 CEST
CC6DF5A424E6442D9507ABA61B686FAC

Pasi Vihinen
Manager of Sustaining
Engineering, EMEA

Tuomas Huhmarniemi
Head of Quality