



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 products

Model & Name 730981 Charging Stand
730991 Charging Carousel

Device type Electronic laboratory equipment

in the form as delivered fulfil all the relevant provisions of the following European Directives – including any amendments valid at the time this declaration was signed – and meet 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

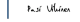
EN 61326-1:2013

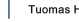
EN 61326-1:2021

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

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 | 08:44:25 CEST
B63019D13BB549149E052159F2E6C344

Signed by:

Signer Name: Tuomas Huhmarniemi
Signing Reason: Hyväksyn tämän asiakirjan
Signing Time: 2024-10-15 | 16:17:13 CEST
CC6DF5A424E6442D9507ABA81B686FAC

Pasi Vihinen
Manager of Sustaining
Engineering, EMEA

Tuomas Huhmarniemi
Head of Quality