



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 Picus® 2

Model(s) LH-747021, LH-747041, LH-747061, LH-747081, LH-747101, LH-747111,
LH-747321, LH-747341, LH-747361, LH-747391, LH-747421, LH-747441,
LH-747461, LH-747491

Device type Electronic laboratory equipment, radio equipment for laboratory use

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:

2014/53/EU Directive on Radio Equipment

EN 300 328 V2.2.2
EN 301 489-17 V3.2.4

2014/30/EU Directive on Electromagnetic Compatibility

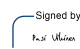
EN 61326-1:2013
EN 61326-1:2021
EN 301 489-1 V2.2.3

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 Charging Stand 730981
Charging Carousel 730991
Picus® AC Adapter LH-735001
Sartorius Pipetting mobile application

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 | 08:43:18 CEST
B63019D13BB549149E052159F2E6C344

Pasi Vihinen
Manager of Sustaining
Engineering, EMEA

Signed by:

Signer Name: Tuomas Huhmarniemi
Signing Reason: Hyväksyn tämän asiakirjan
Signing Time: 2024-10-15 | 16:15:21 CEST
CC6DF5A424E6442D9507ABA61B686FAC

Tuomas Huhmarniemi
Head of Quality