

**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 UK Regulations and meets the applicable requirements of the UK Designated Standards including any amendments valid at the time this declaration was signed listed below:


**The Radio Equipment Regulations 2017**  
**UK Statutory Instruments 2017 No. 1206**  
EN 300 328 v. 2.2.2  
EN 301 489-17 V3.2.4

**The Electromagnetic Compatibility Regulations 2016**  
**UK Statutory Instruments 2016 No. 1091**  
BS EN 61326-1:2013  
BS EN 61326-1:2021  
EN 301 489-1 V2.2.3


**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** 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 | 08:43:05 CEST  
B63019D13BB549148E052159F2E6C344

Pasi Vihinen  
Manager of Sustaining  
Engineering, EMEA

Signed by:  
  
Signer Name: Tuomas Huhmarniemi  
Signing Reason: Hyväksyn tämän asiakirjan  
Signing Time: 2024-10-15 | 16:15:03 CEST  
CC6DF5A424E6442D9507ABA61B686FAC

Tuomas Huhmarniemi  
Head of Quality