

UK Declaration of Conformity

Manufacturer Sartorius Biohit Liquid Handling Oy
Laippatie 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 pipette

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
EN 301 489-17 V3.2.4

The Electromagnetic Compatibility Regulations 2016
UK Statutory Instruments 2016 No. 1091
BS EN 61326-1:2013
EN 301 489-1 V1.9.2

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 Biohit Liquid Handling Oy
Helsinki, 29 June 2023

Hannes Jantunen
Head of Development

Tuomas Huhmarniemi
Head of Quality