



UK Declaration of Conformity

Product: **THERMOstar (0300)**

BMG LABTECH GmbH declares the conformity of the product THERMOstar (type 0300) to the following UK Regulation(s):

The Electrical Equipment (Safety) Regulations 2016
The Electromagnetic Compatibility Regulations 2016
The Restriction of the Use of Certain Hazardous Substances in Electrical and Electronic Equipment Regulations 2012

Applied harmonised standards:

Safety:

BS EN 61010-1: 2010 Safety requirements for electrical equipment for measurement, control and laboratory use

EMC:

BS EN 61326-1: 2013 Electrical equipment for measurement, control and laboratory use – EMC requirements

BS EN 61000-3-2: 2014 Limits for harmonic current emissions

BS EN 61000-3-3: 2013 Limitation of voltage fluctuations and flicker in low-voltage systems...

BS EN 55011: 2016 Industrial, scientific and medical equipment – Limits CISPR 11

Restriction of the use of certain hazardous substances in electronic equipment (RoHS):

BS EN IEC 63000: 2018 Technical documentation for the assessment of electrical and electronic products with respect to the restriction of hazardous substances

Ortenberg, 2022-09-19

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)