



UK Declaration of Conformity

Product: **CLARIOstar^{PLUS} (0430)**

BMG LABTECH GmbH declares the conformity of the product CLARIOstar^{PLUS} (type 0430) 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
BS EN 62471: 2008	Photobiological safety of lamps and lamp systems
BS EN 60825-1: 2014	Safety of Laser Products

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. Rübiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)