



# UK Declaration of Conformity

**Product: PHERAstar FSX (0472)**

BMG LABTECH GmbH declares the conformity of the product PHERAstar FSX (type 0472) 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:

<b>BS EN 61010-1: 2010</b>	Safety requirements for electrical equipment for measurement, control and laboratory use
<b>BS EN 62471: 2008</b>	Photobiological safety of lamps and lamp systems
<b>BS EN 60825-1: 2014</b>	Safety of laser products

### EMC:

<b>BS EN 61326-1: 2013</b>	Electrical equipment for measurement, control and laboratory use – EMC requirements
<b>BS EN 61000-3-2: 2014</b>	Limits for harmonic current emissions
<b>BS EN 61000-3-3: 2013</b>	Limitation of voltage fluctuations and flicker in low-voltage systems...
<b>BS EN 55011: 2016</b>	Industrial, scientific and medical equipment – Limits CISPR 11

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

<b>BS EN IEC 63000: 2018</b>	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)