



# UK Declaration of Conformity

Product: **VANTAstar (0421)**

BMG LABTECH GmbH declares the conformity of the product VANTAstar (type 0421) including option pump module (type 0270) 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

### 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, 2022-09-15

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)