



# EU Declaration of Conformity

**Product: NEPHELOstar<sup>Plus</sup> (0506)**

BMG LABTECH GmbH declares the conformity of the product NEPHELOstar<sup>Plus</sup> (type 0506) to the following standards or directive documents:

## Safety:

**EN 61010-1: 07-2011** Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements  
**2014/35/EU: 02-2014** EU Low Voltage Directive  
**EN 60825-1: 07-2015** Safety of Laser Products

## EMC:

**2014/30/EU: 02-2014** EMC Guideline  
**EN 61326-1: 07-2013** Electrical equipment for measurement, control and laboratory use...  
**EN 55011: 04-2011** Industrial, scientific and medical (ISM) radio-frequency equipment...  
**EN 61000-3-2: 03-2015** Limits for harmonic current emissions  
**EN 61000-3-3: 03-2014** Limitation of voltage fluctuations and flicker in low-voltage systems...  
**EN 61000-4-2: 12-2009** Electrostatic discharge immunity test  
**EN 61000-4-3: 04-2011** Radiated, radio-frequency, electromagnetic field immunity test  
**EN 61000-4-4: 04-2013** Electrical fast transient/burst immunity test  
**EN 61000-4-5: 03-2015** Surge immunity test  
**EN 61000-4-6: 08-2014** Immunity to conducted disturbances, induced by radio-frequency fields  
**EN 61000-4-11: 02-2005** Voltage dips, short interruptions and voltage variations immunity tests  
**EN 62471: 03-2009** Photobiological safety of lamps and lamp systems

## Restriction of the use of certain hazardous substances in electronic equipment

**2011/65/EU 2011-06-08** RoHS-Directive

Ortenberg, 2015-12-14

Quality Representative N. Herzog

Managing Director T. Rübiger

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)