



## EU Declaration of Conformity

**Product: NEPHELOstar (0504)**

BMG LABTECH GmbH declares the conformity of the product NEPHELOstar (type 0504) to the following standards or normative documents:

### Safety:

**EN 61010-1:2001** (Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements), **Classification VDE 0411 Part 1 + A1:2002**  
**EU Low Voltage (73/23/EWG) 1993**  
**EN 60825-1:10-2003 Safety of Laser Products**

### EMC:

**DIN EN 61326-1/1997 + DIN EN 61326-A1:1998 + DIN EN 61326-A2:2001 + DIN EN 61326-A3:2003**  
(Electrical equipment for measurement, control and laboratory use...)  
**DIN EN 55011/2003** (Industrial, scientific and medical (ISM) radio-frequency equipment...)  
**DIN EN 61000-3-2/2001** (Limits for harmonic current emissions)  
**DIN EN 61000-3-3/2002** (Limitation of voltage fluctuations and flicker in low-voltage systems... $\leq 16A$ )  
**DIN EN 61000-4-2/2001** (Electrostatic discharge immunity test)  
**DIN EN 61000-4-3/2003** (Radiated, radio-frequency, electromagnetic field immunity test)  
**DIN EN 61000-4-4/2002** (Electrical fast transient/burst immunity test)  
**DIN EN 61000-4-5/2001** (Surge immunity test)  
**DIN EN 61000-4-6/2001** (Immunity to conducted disturbances, induced by radio-frequency fields)  
**DIN EN 61000-4-11/2001** (Voltage dips, short interruptions and voltage variations immunity tests)

Offenburg, 2007-08-06

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH  
Hanns-Martin-Schleyer-Str. 10  
77656 Offenburg (GERMANY)