



# EU Declaration of Conformity

**Product:** **RUBYstar (0460)**

BMG LABTECH GmbH declares the conformity of the product RUBYstar (type 0460) to the following standards or normative documents:

## Safety:

**EN 61010-1:08-2002** (Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements), **Classification VDE 0411 Part 1 + A1:11-2002 + A2:01-2004 2006/95/EG:12-2006** (EU Low Voltage Guideline)  
**EN 60825-1:10-2003 Safety of Laser Products**

## EMC:

**2004/108/EG:12-2004** (EMC Guideline)  
**EN 61326-1:10-2006** (Electrical equipment for measurement, control and laboratory use...)  
**EN 55011:08-2003** (Industrial, scientific and medical (ISM) radio-frequency equipment...)  
**EN 61000-3-2:10-2006** (Limits for harmonic current emissions)  
**EN 61000-3-3:07-2006** (Limitation of voltage fluctuations and flicker in low-voltage systems...<=16A)  
**EN 61000-4-2:12-2001** (Electrostatic discharge immunity test)  
**EN 61000-4-3:12-2006** (Radiated, radio-frequency, electromagnetic field immunity test)  
**EN 61000-4-4:07-2005** (Electrical fast transient/burst immunity test)  
**EN 61000-4-5:07-2006** (Surge immunity test)  
**EN 61000-4-6:12-2001** (Immunity to conducted disturbances, induced by radio-frequency fields)  
**EN 61000-4-11:02-2005** (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)