



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)