



EU Declaration of Conformity

Product: **Omega (0415)**

BMG LABTECH GmbH declares the conformity of the product Omega (type 0415) 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

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. Räßler

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)