



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:

2014/35/EU: 02-2014 EU Low Voltage Directive
EN 61010-1: 03-2020 Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements

EMC:

2014/30/EU: 02-2014 EU EMC Guideline
EN 61326-1: 07-2013 Electrical equipment for measurement, control and laboratory use – EMC requirements
EN 61000-3-2: 12-2019 Limits for harmonic current emissions
EN 61000-3-3: 07-2020 Limitation of voltage fluctuations and flicker in low-voltage systems...
EN 61000-4-2: 12-2009 Electrostatic discharge immunity test
EN 61000-4-3: 11-2021 Radiated, radio-frequency, electromagnetic field immunity test
EN 61000-4-4: 04-2013 Electrical fast transient/burst immunity test
EN 61000-4-5: 03-2019 Surge immunity test
EN 61000-4-6: 08-2014 Immunity to conducted disturbances, induced by radio-frequency fields
EN 61000-4-11: 10-2021 Voltage dips, short interruptions and voltage variations immunity tests
EN 55011: 05-2022 Industrial, scientific and medical equipment – Limits CISPR 11
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 2 -Directive, including 2015/863/EU (RoHS 3)
1907/2006/EU: 2007-07-01 REACH Regulation

Ortenberg, 2022-09-19

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)