



EU Declaration of Conformity

Product: **PHERAS[®] FSX (0472)**

BMG LABTECH GmbH declares the conformity of the product PHERAS[®] FSX (type 0472) 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
EN 60825-1: 07-2015 Safety of Laser Products

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. Räßiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)