



EU Declaration of Conformity

Product: SPECTROstar^{Nano} (0601)

BMG LABTECH GmbH declares the conformity of the product SPECTROstar^{Nano} (type 0601) 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, 2022-09-20

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)