



EU Declaration of Conformity

Product: **VANTAstar (0421)**

BMG LABTECH GmbH declares the conformity of the product VANTAstar (type 0421) including option pump module (type 0270) to the following standards or directive documents:

Safety:

2014/35/EU: 02-2014
EN 61010-1: 03-2020

EU Low Voltage Directive
Safety requirements for electrical equipment for measurement, control and laboratory use

EMC:

2014/30/EU: 02-2014
EN 61326-1: 07-2013

EU EMC Guideline
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, 2022-09-13

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH
Allmendgrün 8
77799 Ortenberg (GERMANY)