



# EU Declaration of Conformity

Product: **Atmospheric Control Unit  
ACU (0221)**

BMG LABTECH GmbH declares the conformity of the product Atmospheric Control Unit (ACU type 0221) 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

**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. Rübiger

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)