



# 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:**

**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

**EMC:**

<b>2014/30/EU: 02-2014</b>	EMC Guideline
<b>EN 61326-1: 10-2013</b>	Electrical equipment for measurement, control and laboratory use...
<b>EN 55011: 04-2011</b>	Industrial, scientific and medical (ISM) radio-frequency equipment...
<b>EN 61000-3-2: 03-2010</b>	Limits for harmonic current emissions
<b>EN 61000-3-3: 06-2009</b>	Limitation of voltage fluctuations and flicker in low-voltage systems...
<b>EN 61000-4-2: 12-2009</b>	Electrostatic discharge immunity test
<b>EN 61000-4-3: 04-2011</b>	Radiated, radio-frequency, electromagnetic field immunity test
<b>EN 61000-4-4: 11-2010</b>	Electrical fast transient/burst immunity test
<b>EN 61000-4-5: 06-2007</b>	Surge immunity test
<b>EN 61000-4-6: 12-2009</b>	Immunity to conducted disturbances, induced by radio-frequency fields
<b>EN 61000-4-11: 02-2005</b>	Voltage dips, short interruptions and voltage variations immunity tests

**Restriction of the use of certain hazardous substances in electronic equipment**

**2011/65/EU 2011-06-08** RoHS-Directive

Ortenberg, 2016-04-20

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)