



# UK Declaration of Conformity

**Product: LVis Plate (0680)**

BMG LABTECH GmbH declares the conformity of the product LVis Plate (type 0680) to the following UK Regulation(s):

**The Electrical Equipment (Safety) Regulations 2016  
The Restriction of the Use of Certain Hazardous Substances  
in Electrical and Electronic Equipment Regulations 2012**

**Applied harmonised standards:**

**Safety:**

**BS EN 61010-1: 2010**

Safety requirements for electrical equipment for measurement, control and laboratory use

**Restriction of the use of certain hazardous substances in electronic equipment (RoHS):**

**BS EN IEC 63000: 2018**

Technical documentation for the assessment of electrical and electronic products with respect to the restriction of hazardous substances

Ortenberg, 2022-09-19

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)