



EU Declaration of Conformity

Product: Evaluation Plate (0681)

BMG LABTECH GmbH declares the conformity of the product Evaluation Plate (type 0681) 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

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)