



# EU Declaration of Conformity

**Product: LVis Plate (0680)**

BMG LABTECH GmbH declares the conformity of the product LVis Plate (type 0680) to the following standards or directive documents:

## Safety:

**EN 61010-1:2001** (Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements), **Classification VDE 0411 Part 1 + A1:11-2002 + A2:01-2004 2006/95/EG:12-2006** (EU Low Voltage Guideline)

Ortenberg, 2013-02-01

Quality Representative N. Herzog

Managing Director T. Räßler

BMG LABTECH GmbH  
Allmendgrün 8  
77799 Ortenberg (GERMANY)