



# EU Declaration of Conformity

**Product:** **NOVOstar (0700)**

BMG LABTECH GmbH declares the conformity of the product NOVOstar (type 0700) to the following standards or directive documents:

## Safety:

**EN 61010-1:08-2002** (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)

## EMC:

**2004/108/EG:12-2004** (EMC Guideline)

**EN 61326-1:10-2006** (Electrical equipment for measurement, control and laboratory use...)

**EN 55011:08-2003** (Industrial, scientific and medical (ISM) radio-frequency equipment...)

**EN 61000-3-2:10-2006** (Limits for harmonic current emissions)

**EN 61000-3-3:07-2006** (Limitation of voltage fluctuations and flicker in low-voltage systems...<=16A)

**EN 61000-4-2:12-2001** (Electrostatic discharge immunity test)

**EN 61000-4-3:12-2006** (Radiated, radio-frequency, electromagnetic field immunity test)

**EN 61000-4-4:07-2005** (Electrical fast transient/burst immunity test)

**EN 61000-4-5:07-2006** (Surge immunity test)

**EN 61000-4-6:12-2001** (Immunity to conducted disturbances, induced by radio-frequency fields)

**EN 61000-4-11:02-2005** (Voltage dips, short interruptions and voltage variations immunity tests)

Offenburg, 2007-07-31

Quality Representative N. Herzog

Managing Director T. Räßiger

BMG LABTECH GmbH  
Hanns-Martin-Schleyer-Str. 10  
77656 Offenburg (GERMANY)