

CE**EU Declaration of Conformity**

Product:

FLUOstar *

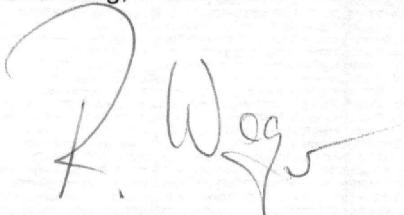
BMG LabTechnologies GmbH declares the conformity of FLUOstar (type 0403) to the following standards or normative documents:

Security of product:

**DIN EN 61010 Part 1/03.94
Classification VDE 0411 Part 1/03.94
IEC 1010-1/09.90 and Amendment 1/09.92
EU Low Voltage (73/23/EWG)**

EMC:

EN 61326-1/1997 and A1/06.1998 (El. equipment for measurement, control and laboratory use...)
EN 55011/1998 (Industrial, scientific and medical (ISM) radio-frequency equipment...)
EN 61000-3-2/1995 and A12/1995 (Limits for harmonic current emissions)
EN 61000-3-3/1995 (Limitation of voltage fluctuations and flicker in low-voltage s. systems...<=16A)
EN 61000-4-2/1995 (Electrostatic discharge immunity test)
EN 61000-4-3/1995 (Radiated, radio-frequency, electromagnetic field immunity test)
EN 61000-4-4/1995 (Electrical fast transient/burst immunity test)
EN 61000-4-5/1995 (Surge immunity test)
EN 61000-4-6/1996 (Immunity to conducted disturbances, induced by radio-frequency fields)
EN 61000-4-11/1994 (Voltage dips, short interruptions and voltage variations immunity tests)

Offenburg, November 24th 1999

Quality Manager René Wagner



Managing Director T. Räbiger

BMG Labtechnologies GmbH
Hanns-Martin-Schleyer-Str. 10
D – 77656 Offenburg (GERMANY)