



EU Declaration of Conformity

Product:

STACKER (0900)

BMG LABTECH GmbH declares the conformity of the product STACKER (type 0900) to the following standards or normative documents:

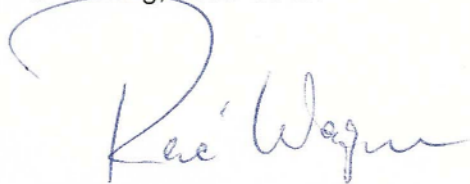
Safety:

EN 61010-1:2001 (Safety requirements for electrical equipment for measurement, control and laboratory use: General requirements), **Classification VDE 0411 Part 1 + A1:2002**
EU Low Voltage (73/23/EWG) 1993

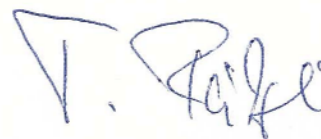
EMC:

DIN EN 61326-1/1997 + DIN EN 61326-A1:1998 + DIN EN 61326-A2:2001 + DIN EN 61326-A3:2003
(Electrical equipment for measurement, control and laboratory use...)
DIN EN 55011/2003 (Industrial, scientific and medical (ISM) radio-frequency equipment...)
DIN EN 61000-3-2/2001 (Limits for harmonic current emissions)
DIN EN 61000-3-3/2002 (Limitation of voltage fluctuations and flicker in low-voltage systems...<=16A)
DIN EN 61000-4-2/2001 (Electrostatic discharge immunity test)
DIN EN 61000-4-3/2003 (Radiated, radio-frequency, electromagnetic field immunity test)
DIN EN 61000-4-4/2002 (Electrical fast transient/burst immunity test)
DIN EN 61000-4-5/2001 (Surge immunity test)
DIN EN 61000-4-6/2001 (Immunity to conducted disturbances, induced by radio-frequency fields)
DIN EN 61000-4-11/2001 (Voltage dips, short interruptions and voltage variations immunity tests)

Offenburg, 2005-03-24

A handwritten signature in blue ink, appearing to read "R. Wagner".

Quality Assurance Manager R. Wagner

A handwritten signature in blue ink, appearing to read "T. Rübiger".

Managing Director T. Rübiger

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