

EC-Declaration

We herewith confirm that the construction of the following indicated products / systems / units is brought into circulation to comply with the relevant EC regulations listed.

Description: **Concentration- and Moisture-Measuring Systems
Micro-Polar 2, Micro-Polar 2 ++,
Micro-Polar Moist, and Micro-Polar Moist ++**

Type: **LB 567-XX and LB 568-XX**

	Richtlinie und Änderungen	angewendete Normen	
EMC	2004/108/EC	EN 55011	1998 +A1:1999 +A2:2002
		EN 61326-1	2006-05
		EN 61000-4-2	1995 +A1:1998 +A2:2001
		EN 61000-4-3	2006-12
		EN 61000-4-4	2004
		EN 61000-4-5	1995 +A1:2001
		EN 61000-4-6	1996 +A1:2001
		EN 61000-4-11	1994-08 +A1:2001-02
		Namur NE21	2004
LVD	2006/95/EC	EN 61010 part 1	2002-08

This declaration is issued by the manufacturer

BERTHOLD TECHNOLOGIES GmbH & Co. KG
Calmbacher Str. 22
75323 Bad Wildbad, Germany

released by



Dr. Wilfried Reuter – Technical Director
Bad Wildbad, 28th of April 2010

Registergericht / Court of Registration
Persönlich haftende Gesellschafterin / Fully liable Associates
Registergericht / Court of Registration
Geschäftsführung / Management
USt.-Id-Nr. / VAT Reg. No.
Deutsche Steuernummer / German Tax No.
WEE-Reg. No.

Stuttgart HRA 330991
BERTHOLD TECHNOLOGIES Verwaltungs-GmbH
Stuttgart HRB 331520
Hans J. Oberhofer, Dr. Wilfried Reuter
DE813050511
49038/08038
DE99468690

Sparkasse PF-CW	75323 Bad Wildbad	Konto/Account No. 8 045 003 (BLZ 666 500 85)	SWIFT-BIC PZHSDE66	IBAN: DE37 6665 0085 0008 0450 03
Volksbank	75119 Pforzheim	Konto/Account No. 957 004 (BLZ 666 900 00)	SWIFT-BIC VBPFDE66	IBAN: DE85 6669 0000 0000 9570 04
Commerzbank	75105 Pforzheim	Konto/Account No. 6 511 120 (BLZ 666 800 13)	SWIFT-BIC DRESDEFF 666	IBAN: DE05 6668 0013 0651 1120 00