



IECEx Certificate of Conformity

INTERNATIONAL ELECTROTECHNICAL COMMISSION IEC Certification System for Explosive Atmospheres

for rules and details of the IECEx Scheme visit www.iecex.com

Certificate No.: **IECEx EPS 19.0073**

Page 1 of 4

Certificate history:

Status: **Current**

Issue No: 1

Issue 0 (2020-01-23)

Date of Issue: 2021-03-15

Applicant: **Berthold Technologies GmbH & Co. KG**
Calmbacher Str. 22
75323 Bad Wildbad
Germany

Equipment: **Evaluation unit, type LB 47x-xx-xx**

Optional accessory:

Type of Protection: **ib**

Marking: [Ex ib Gb] IIB

[Ex ib Db] IIIC

[Ex ib Gb] IIC

Approved for issue on behalf of the IECEx
Certification Body:

Position:

Signature:
(for printed version)

Date:



1. This certificate and schedule may only be reproduced in full.
2. This certificate is not transferable and remains the property of the issuing body.
3. The Status and authenticity of this certificate may be verified by visiting www.iecex.com or use of this QR Code.



Certificate issued by:

Bureau Veritas Consumer Products Services Germany GmbH
Businesspark A96
86842 Türkheim
Germany





IECEx Certificate of Conformity

Certificate No.: **IECEx EPS 19.0073**

Page 2 of 4

Date of issue: 2021-03-15

Issue No: 1

Manufacturer: **Berthold Technologies GmbH & Co. KG**
Calmbacher Str. 22
75323 Bad Wildbad
Germany

Additional
manufacturing
locations:

This certificate is issued as verification that a sample(s), representative of production, was assessed and tested and found to comply with the IEC Standard list below and that the manufacturer's quality system, relating to the Ex products covered by this certificate, was assessed and found to comply with the IECEx Quality system requirements. This certificate is granted subject to the conditions as set out in IECEx Scheme Rules, IECEx 02 and Operational Documents as amended

STANDARDS :

The equipment and any acceptable variations to it specified in the schedule of this certificate and the identified documents, was found to comply with the following standards

IEC 60079-0:2017 Explosive atmospheres - Part 0: Equipment - General requirements
Edition:7.0

IEC 60079-11:2011 Explosive atmospheres - Part 11: Equipment protection by intrinsic safety "i"
Edition:6.0

This Certificate **does not** indicate compliance with safety and performance requirements other than those expressly included in the Standards listed above.

TEST & ASSESSMENT REPORTS:

A sample(s) of the equipment listed has successfully met the examination and test requirements as recorded in:

Test Report:

[DE/EPS/ExTR19.0071/01](#)

Quality Assessment Report:

[DE/PTB/QAR06.0011/06](#)



IECEx Certificate of Conformity

Certificate No.: **IECEx EPS 19.0073**

Page 3 of 4

Date of issue: 2021-03-15

Issue No: 1

EQUIPMENT:

Equipment and systems covered by this Certificate are as follows:

The LB 47x-xx-xx (associated apparatus) is used for the evaluation and display of signals as well as for power supply and data transmission to the detector (LB 44xx or LB 4700) by means of FSK modulation. The LB 47x-xx-xx is installed in the safe area and provides an intrinsically safe interface to the detector, which may be installed in zone 1 (gas) or zone 21 (dust).

The evaluation unit type LB 47x-xx-xx needs a partition wall between the intrinsically safe circuit and the non-is circuits to ensure the required 50mm distances between them.

SPECIFIC CONDITIONS OF USE: NO



IECEx Certificate of Conformity

Certificate No.: **IECEx EPS 19.0073**

Page 4 of 4

Date of issue: 2021-03-15

Issue No: 1

DETAILS OF CERTIFICATE CHANGES (for issues 1 and above)

Additional inclusion of a new variant for IIC (see below).

Annex:

[IECEx EPS 19.0073 - Annex_01.pdf](#)



Annex to Certificate
IECEx EPS 19.0073 Issue No.: 1



Description of equipment:

The LB 47x-xx-xx (associated apparatus) is used for the evaluation and display of signals as well as for power supply and data transmission to the detector (LB 44xx or LB 4700) by means of FSK modulation. The LB 47x-xx-xx is installed in the safe area and provides an intrinsically safe interface to the detector, which may be installed in zone 1 (gas) or zone 21 (dust).

Type key:

LB 47x	-	-	-	-	-	Description
x = 0,2..9	0					non Ex
	1					Ex-ib IIB / IIIC
	2					Ex-ib IIC / IIIC
		1				24 V _{DC}
		2				100...240 V _{AC} (50/60 Hz)
				M		Master
					x	one or more digits, not relevant for approval

Electrical data:

Power supply (ST1-30A,30C, PE: ST1-32A,32C)

U_m = 100 – 240 VAC

U_n = 32 VDC (Option 1)

U_n = 100 – 240 VAC (Option 2)

FSK interface

(Pins ST1-2a, ST1-2c)

in type of protection Intrinsic Safety Ex ib IIIC, IIB, IIC only for the connection of certified intrinsically safe circuits.

Max. values: (output)

at IIC:

U_o = 16.23 V

I_o = 60 mA

P_o = 1.0 W

L_o = 0.37 mH

C_o = 0.16 µF

C_i = 22 nF

L_i = 13 µH



Annex to Certificate
IECEx EPS 19.0073 Issue No.: 1



at IIB/ IIIC:

$U_o = 16.23 \text{ V}$
 $I_o = 60 \text{ mA}$
 $P_o = 1.0 \text{ W}$
 $L_o = 22.0 \text{ mH}$
 $C_o = 0.6 \text{ }\mu\text{F}$

$C_i = 22 \text{ nF}$
 $L_i = 13 \text{ }\mu\text{H}$

Ambient temperature:

19'' Rack: $-20 \text{ }^\circ\text{C} \leq T_a \leq +50 \text{ }^\circ\text{C}$

Wall mounted housing: $-20 \text{ }^\circ\text{C} \leq T_a \leq +45 \text{ }^\circ\text{C}$