



SITIiAS
Worldwide Access

防爆合格证

证 号: GYJ23.1352X

制 造 商 Berthold Technologies GmbH & Co. KG

(地址: Calmbacher Str. 22 75323 Bad Wildbad Germany)

产 品 名 称 评估单元

型 号 规 格 LB 47x-xx-xx

防 爆 标 志 [Ex ib Gb] IIB
[Ex ib Db] IIIC
[Ex ib Gb] IIC

产 品 标 准 /

图 样 编 号 57724、65101、68339、65098

经图样及技术文件的审查和样品检验, 确认上述产品符合下列标准:

GB/T 3836.1-2021, GB/T 3836.4-2021

特颁发此证。

本证书有效期: 2023 年 07 月 19 日 至 2028 年 07 月 18 日

- 备 注
1. 安全使用注意事项见本证书附件。
 2. 证书编号后缀“X”表明产品具有安全使用特殊条件, 内容见本证书附件。
 3. 型号规格说明见本证书附件。
 4. 本安电气参数见本证书附件。



批 准

上海仪器仪表自控系统检验测试所有限公司
国家级仪器仪表防爆安全监督检验站
颁发日期二〇二三年七月十九日

本证书仅对与认可文件和样品一致的产品有效。

地址: 上海市漕宝路103号
邮编: 200233

网址: www.nepsi.org.cn
Email: info@nepsi.org.cn

电话: +86 21 64368180
传真: +86 21 64844580



SitiAs
Worldwide Access

EXPLOSION PROTECTION CERTIFICATE OF CONFORMITY

Cert No. GYJ23.1352X

Manufacturer	Berthold Technologies GmbH & Co. KG (Address: Calmbacher Str. 22 75323 Bad Wildbad Germany)
Product	Evaluation unit
Model	LB 47x-xx-xx
Ex marking	[Ex ib Gb] IIB [Ex ib Db] IIIC [Ex ib Gb] IIC
Product standard	/
Drawing number	57724、65101、68339、65098

The product was found to comply with the following standard(s):

GB/T 3836.1-2021,GB/T 3836.4-2021

Valid until: 2028.07.18

Remarks

- 1.Conditions for safe use are specified in the attachment(s) to this certificate.
- 2.Symbol "X" placed after the certification number denotes specific conditions of use, which are specified in the attachment(s) to this certificate.
- 3.Model designation is specified in the attachment(s) to this certificate.
- 4.Intrinsic safety parameters specified in the attachment(s) to this certificate.



Approval

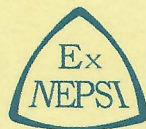
Shanghai Inspection and Testing Institute of
Instruments and Automation Systems Co., Ltd.
National Supervision and Inspection Center for
Explosion Protection and Safety of Instrumentation
Date of issue 2023.07.19

This Certificate is valid for products compatible with the documents and samples approved by NEPSI.

103 Cao Bao Road
Shanghai 200233, China

<http://www.nepsi.org.cn>
Email: info@nepsi.org.cn

Tel: +86 21 64368180
Fax: +86 21 64844580



(GYJ23.1352X)

(Attachment I)

GYJ23.1352X防爆合格证附件 I

由Berthold Technologies GmbH & Co. KG生产的LB 47x-xx-xx评估单元，经国家级仪器仪表防爆安全监督检验站(NEPSI)检验，符合下列标准：

GB/T 3836.1-2021 爆炸性环境 第1部分：设备 通用要求

GB/T 3836.4-2021 爆炸性环境 第4部分：由本质安全型“i”保护的设备

产品防爆标志：[Ex ib Gb] IIB、[Ex ib Db] IIIC、[Ex ib Gb] IIC，防爆合格证号为GYJ23.1352X。

本证书认可的产品具体型号为：

LB 47x-xx-xx:

LB47 *a-b c-d e*

a 代表产品选型代码：0、2、3、4、5、6、7、8、9；

b 代表防爆选型代码：0（非防爆型）、1（防爆标志为[Ex ib Gb] IIB、[Ex ib Db] IIIC）、2（防爆标志为[Ex ib Gb] IIC、[Ex ib Db] IIIC）；

c 代表电气参数选型代码：1（24 VDC）、2（100~240 VAC 50/60 Hz）；

d 代表产品描述代码：可为空白、或M（Master）；

e 代表批次选型，一个或多个数字，与防爆性能无关。

一、产品安全使用特殊条件

产品防爆合格证号后缀“X”代表产品安全使用有特殊条件：

产品的使用环境温度范围，19”机架：-20℃~+50℃，壁挂式外壳：-20℃~+45℃。

二、产品使用注意事项

1. 电气安全参数：

电源（ST1-30A，30C，PE：ST1-32A，32C）

$U_m = 100 \sim 240 \text{ VAC}$

$U_n = 32 \text{ VDC (Option 1)}$

$U_n = 100 \sim 240 \text{ VDC (Option 2)}$

FSK接口（引脚ST1-2a、ST1-2c）：本质安全型 Ex ib IIIC、IIB、IIC 仅用于连接经认证的本质安全电路。

本安参数及内部最大等效参数（输出值）如下：

选型IIC：

最高输出 电压 U_o (V)	最大输出 电流 I_o (mA)	最大输出 功率 P_o (W)	最大内部等效参数			
			C_o (μ F)	C_i (nF)	L_o (mH)	L_i (μ H)
16.23	60	1.0	0.16	22	0.37	13

选型IIB/ IIIC：

最高输出 电压 U_o (V)	最大输出 电流 I_o (mA)	最大输出 功率 P_o (W)	最大内部等效参数			
			C_o (μ F)	C_i (nF)	L_o (mH)	L_i (μ H)
16.23	60	1.0	0.6	22	22.0	13

2. 该产品必须与已通过防爆认证的关联设备配套共同组成本安防爆系统方可使用于爆炸性气体环境。其系统接线必须同时遵守本产品 and 所配关联设备的使用说明书要求，接线端子不得接错。

3. 用户不得自行更换该产品的零部件，应会同产品制造商共同解决运行中出现的故障，以杜绝损坏现象的发生。

4. 产品的安装、使用和维护应同时遵守产品使用说明书、GB/T3836.13-2021“爆炸性环境第13部分：设备的修理、检修、修复和改造”、GB/T 3836.15-2017“爆炸性环境 第15部分：电气装置的设计、选型和安装”、GB/T 3836.16-2022“爆炸性环境 第16部分：电气装置的检查与维护”、GB/T3836.18-2017 爆炸性环境 第18部分：本质安全电气系统和GB50257-2014“电气装置安装工程爆炸和火灾危险环境电力装置施工及验收规范”的有关规定。

三、制造厂责任

1. 制造厂必须将上述产品安全使用特殊条件及使用注意事项纳入该产品使用说明书。
2. 制造厂必须严格按照NEPSI认可的文件资料生产。

上海仪器仪表自控系统检验测试所有限公司

国家级仪器仪表防爆安全监督检验站

二〇二三年七月十九日



(GYJ23.1352X)

(Attachment I)

Attachment I to GYJ23.1352X

1. Description

Type LB 47x-xx-xx Evaluation unit, manufactured by Berthold Technologies GmbH & Co. KG have been certified by National Supervision and Inspection Center for Explosion Protection and Safety of Instrumentation (NEPSI), according with following standards:

GB/T 3836.1-2021 Explosive atmospheres - Part 1: Equipment - General requirements

GB/T 3836.4-2021 Explosive atmospheres - Part 4: Equipment protection by intrinsic safety “i”

Type	Ex marking
LB 47x-xx-xx	[Ex ib Gb] IIB 、 [Ex ib Db] IIIC 、 [Ex ib Gb] IIC

Certificate number is GYJ23.1352X.

Type LB 47 series

LB47 *a-b c-d e*

a representative product selection code: 0、2、3、4、5、6、7、8、9;

b representative explosion-proof selection code: 0 (Non explosion-proof type) 、 1 (The Ex mark [Ex ib Gb] IIB、 [Ex ib Db] IIIC) 、 2 (The Ex mark [Ex ib Gb] IIC、 [Ex ib Db] IIIC) ;

c representative electrical parameter selection code: 1 (24 VDC) 、 2 (100~240 VAC 50/60 Hz) ;

d representative product description code: Can be blank, or M (Master);

e representative batch selection, one or more numbers, independent of explosion-proof performance.

2. Special Condition for Safe Use

Ambient temperature range is: $-20^{\circ}\text{C} \leq T_a \leq +50^{\circ}\text{C}$ (19" Rack)、 $-20^{\circ}\text{C} \leq T_a \leq +45^{\circ}\text{C}$ (Wall mounted housing) .

3. Conditions for Safe Use

3.1 Electrical safety parameters:

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

$U_m = 100 \sim 240 \text{ VAC}$

$U_n = 32 \text{ VDC}$ (Option 1)

$U_n = 100 \sim 240 \text{ VDC}$ (Option 2)

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

PINS: ST1-2a、ST1-2c,Max values:(output):

IIC:

Uo (V)	Io (mA)	Po (W)	Maximum internal equivalent parameter			
			Co(μ F)	Ci(nF)	Lo(mH)	Li(μ H)
16.23	60	1.0	0.16	22	0.37	13

IIB/ IIIC:

Uo (V)	Io (mA)	Po (W)	Maximum internal equivalent parameter			
			Co(μ F)	Ci(nF)	Lo(mH)	Li(μ H)
16.23	60	1.0	0.6	22	22.0	13

3.2 This product must be paired with associated equipment that has passed explosion-proof certification to form an intrinsic safety explosion-proof system before it can be used in explosive gas environments. The system wiring must comply with the user manual requirements of both this product and the associated equipment, and the wiring terminals must not be connected incorrectly.

3.3 Users are not allowed to replace the components of the product on their own. They should work together with the product manufacturer to solve any faults that occur during operation to prevent damage.

3.4 When installation, use and maintenance of this product, observe following standards:

GB/T3836.13-2021 Explosive atmospheres - Part 13: Equipment repair, overhaul and reclamation

GB/T3836.15-2017 Explosive atmospheres - Part 15: Electrical installations design, selection and erection

GB/T3836.16-2022 Explosive atmospheres - Part 16: Electrical installations inspection and maintenance

GB/T3836.18-2017 Explosive atmospheres - Part 18: Intrinsically safe electrical systems

GB50257-2014 Code for construction and acceptance of electric device for explosion atmospheres and fire hazard electrical equipment installation engineering

4. Manufacturer's Responsibility

4.1 Special condition for safe use specified above should be included in the instruction manual.

4.2 Manufacturing should be done according to the documentation approved by NEPSI.

Shanghai Inspection and Testing Institute of
Instruments and Automation Systems Co., Ltd.
National Supervision and Inspection Center for
Explosion Protection and Safety of Instrumentation
July 19, 2023