

Human Bcl-2 ELISA Kit ab119506

21 References 1 图像

概述

产品名称

检测方法

精确度

样品类型

检测类型

灵敏度

范围

回收率

实验步骤

种属反应性

产品概述

人Bcl-2 ELISA试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall				8.6%

批次间

样品	n	Mean	SD	CV%
Overall				12%

Serum, Plasma, Cell culture extracts

Sandwich (quantitative)

0.5 ng/ml

0.5 ng/ml - 32 ng/ml

71 %

Multiple steps standard assay

与反应: Human

Human Bcl-2 in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit (ab119506) is designed for accurate quantitative measurement of Human Bcl-2 concentrations in cell lysates, serum and plasma.

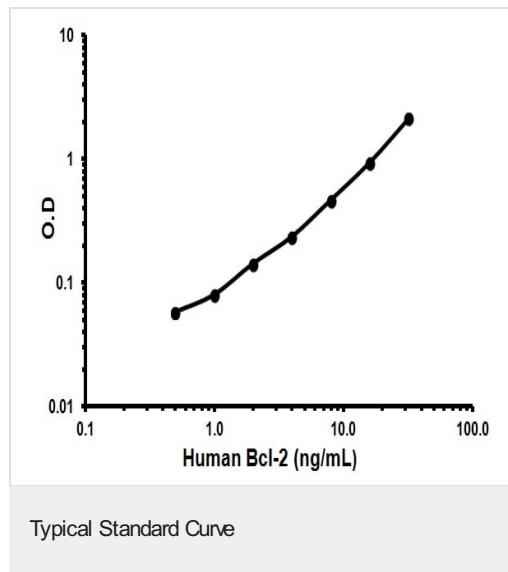
平台	Microplate
性能	
存放说明	Store at +4°C. Please refer to protocols.
组 件	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Biotin-Conjugate anti-human Bcl-2 monoclonal antibody	1 x 70µl
Human Bcl-2 Standard lyophilized (64 ng/mL upon reconstitution)	2 vials
Lysis Buffer (10x)	1 x 15ml
Microplate coated with monoclonal antibody to Bcl-2 (12 x 8 wells)	1 unit
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
Streptavidin-HRP	1 x 150µl
TMB Substrate Solution	1 x 15ml
功 能	Suppresses apoptosis in a variety of cell systems including factor-dependent lymphohematopoietic and neural cells. Regulates cell death by controlling the mitochondrial membrane permeability. Appears to function in a feedback loop system with caspases. Inhibits caspase activity either by preventing the release of cytochrome c from the mitochondria and/or by binding to the apoptosis-activating factor (APAF-1). May attenuate inflammation by impairing NLRP1-inflammasome activation, hence CASP1 activation and IL1B release (PubMed:17418785).
组织特异性	Expressed in a variety of tissues.
疾病相关	A chromosomal aberration involving BCL2 has been found in chronic lymphatic leukemia. Translocation t(14;18)(q32;q21) with immunoglobulin gene regions. BCL2 mutations found in non-Hodgkin lymphomas carrying the chromosomal translocation could be attributed to the Ig somatic hypermutation mechanism resulting in nucleotide transitions.
序列相似性	Belongs to the Bcl-2 family.
结构域	BH1 and BH2 domains are required for the interaction with BAX and for anti-apoptotic activity. The BH4 motif is required for anti-apoptotic activity and for interaction with RAF1 and EGLN3. The loop between motifs BH4 and BH3 is required for the interaction with NLRP1.
翻译后修饰	Phosphorylation/dephosphorylation on Ser-70 regulates anti-apoptotic activity. Growth factor-stimulated phosphorylation on Ser-70 by PKC is required for the anti-apoptosis activity and occurs during the G2/M phase of the cell cycle. In the absence of growth factors, BCL2 appears to be phosphorylated by other protein kinases such as ERKs and stress-activated kinases.

Phosphorylated by MAPK8/JNK1 at Thr-69, Ser-70 and Ser-87, which stimulates starvation-induced autophagy. Dephosphorylated by protein phosphatase 2A (PP2A). Proteolytically cleaved by caspases during apoptosis. The cleaved protein, lacking the BH4 motif, has pro-apoptotic activity, causes the release of cytochrome c into the cytosol promoting further caspase activity. Monoubiquitinated by PARK2, leading to increase its stability. Ubiquitinated by SCF(FBXO10), leading to its degradation by the proteasome.

细胞定位

Mitochondrion outer membrane. Nucleus membrane. Endoplasmic reticulum membrane.

图片



Representative Standard Curve using ab119506.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors