

Human Bcl-2 ELISA Kit ab119506

21 References    画像数 1

製品の概要

製品名	Human Bcl-2 ELISA Kit			
検出方法	Colorimetric			
再現性	Intra-Assay (同時再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				8.6%
	Inter-Assay (日差再現性)			
	サンプル	N	平均値	SD
	Overall			CV%
				12%
サンプルの種類	Serum, Plasma, Cell culture extracts			
アッセイタイプ	Sandwich (quantitative)			
検出感度	0.5 ng/ml			
検出範囲	0.5 ng/ml - 32 ng/ml			
添加回収試験	71 %			
ステップ	Multiple steps standard assay			
種交差性	交差種: Human			
製品の概要	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.			

Bcl-2 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells along with a biotin-conjugated Bcl-2 detection antibody then incubated at room temperature. Following washing, a Streptavidin-HRP conjugate is added to each well, incubated at room temperature and washed. TMB is added and then catalyzed by HRP to produce a blue product that changes to yellow after the addition of acidic stop solution. The density of yellow coloration is directly proportional to the amount of Bcl-2 captured on the plate.

Get results in 90 minutes with Human Bcl-2 ELISA Kit ([ab202411](#)) from our SimpleStep ELISA® range.

## 製品の特性

## 保存方法

Store at +4°C. Please refer to protocols.

内容	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

## 機能

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).

## 組織特異性

Expressed in a variety of tissues.

## 関連疾患

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.

## 配列類似性

Belongs to the Bcl-2 family.

## ドメイン

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.

## 翻訳後修飾

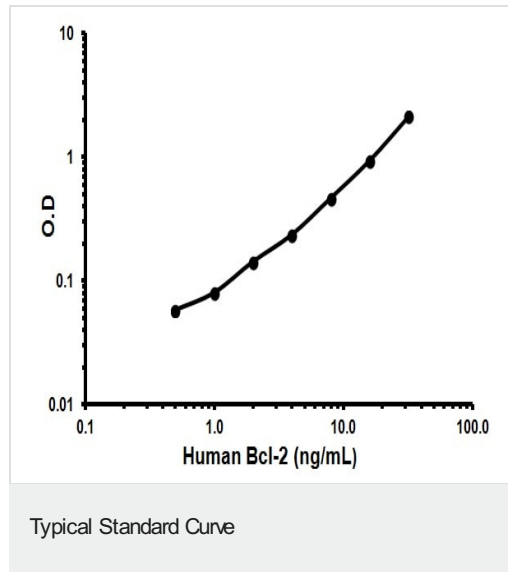
Phosphorylation/dephosphorylation on Ser-70 regulates anti-apoptotic activity. Growth factor-stimulated phosphorylation on Ser-70 by PKC is required for the anti-apoptosis activity and occurs during the G2/M phase of the cell cycle. In the absence of growth factors, BCL2 appears to be phosphorylated 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.

#### 細胞内局在

Mitochondrion outer membrane. Nucleus membrane. Endoplasmic reticulum membrane.

#### 画像



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.co.jp/abpromise> or contact our technical team.

#### Terms and conditions

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