

TREVIGEN[®] Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Anti-Bcl-X_L YTH-2H12 Monoclonal Antibody

Catalog #: 2300-MC-100

Size: 100 µg

Description: The Bcl-2 family of proteins plays a crucial role in the regulation of cell death in many eukaryotic systems. Bcl-X_L has been shown to promote cell survival through the regulation of mitochondrial homeostasis.

Physical State: This antibody is provided as purified immunoglobulin from mouse ascites in 1X PBS containing 0.01% sodium azide.

Immunogen: A synthetic peptide corresponding to amino acids 3 to 14 of the human Bcl-X_L sequence.

Ig Class: IgG_{2a}

Specificity: Anti-Bcl-X_L is specific for human, mouse and rat species.

Storage: Store at 4°C.

Applications: Western analysis, and immunocytochemistry. For western blots, an antibody dilution of 1:1000 is recommended.

Cell Lysates for Western Blotting:

To prepare total cell lysates, cells are solubilized in 1X SDS gel sample buffer (63 mM Tris, pH 6.8, 10% glycerol, 2% SDS, 2.5% β-mercaptoethanol, and 0.0025% bromophenol blue) at 5×10^5 - 1×10^6 cells per ml. The extracts are heated in a boiling water bath for 5 minutes. Electrophoresis on 4-20% Tris-Glycine SDS-PAGE gels.

Procedure for Immunoblotting using Peroxidase Detection:

Blotting buffer: 12 mM Tris base, 96 mM Glycine, and 20% MeOH.

Blocking solution: 5% (w/v) nonfat dry milk in TBS, 0.05% Tween.

Antibody solution: 5% (w/v) nonfat dry milk, 0.05% Tween in TBS.

Transfer the electrophoresed proteins to a nitrocellulose or PVDF membrane and incubate the membrane for 1 hour at room temperature in blocking solution. Incubate the membrane overnight at 4°C in antibody solution containing a 1:1000 dilution of anti-Bcl-X_L YTH-2H12 antibody. Empirical determination of primary antibody concentration will be required for optimal results.

Wash the membrane at room temperature for 15 minutes with 3 changes of 0.05% Tween in TBS. Changing the membrane containers often reduces background.

Incubate the membrane at room temperature for 1 hour in antibody solution containing anti-rabbit conjugated to horseradish peroxidase. Empirical determination of the secondary antibody concentration will be required for optimal results.

Wash the membrane for 15 minutes with 3 changes of 0.05% Tween in TBS.

TREVIGEN[®]

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

Develop peroxidase reaction using e.g. chemiluminescence (Trevigen's PeroxyGlow A, cat# 4855-20-13, and PeroxyGlow B, cat# 4855-20-14).

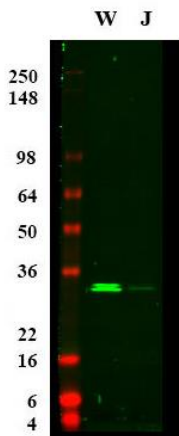


Fig 1. Western blot analysis of mouse Wehi (W) and human Jurkat (J) cell lines. Cells were lysed in Tris-Glycine SDS sample buffer at 10^7 cells/ml and 10 μ l of lysates were loaded per well of a 4-20% Tris-Glycine gel. Proteins were transferred onto an Immobilon FL membrane and mouse and human Bcl-X_L proteins were detected with anti-Bcl-X_L (YTH-2H12) antibody (cat# 2300- MC-100) followed by IR800-conjugated secondary antibody. The membrane was scanned using an Odyssey Infrared Imaging System (Licor).

anti-Bcl-X_L (YTH-2H12) MC (1:1,000)
+
Goat anti-Mouse IR800 (1:5,000)

- References:**
1. Hsu, Y.-T., K.G. Wolter, and R.J. Youle. 1997. Cytosol-to-membrane redistribution of Bax and Bcl-X_L during apoptosis. *Proc. Natl. Acad. Sci. USA* **94**: 3668-3672.
 2. Hsu, Y.-T., and R.J. Youle. 1997. Nonionic detergents induce dimerization among members of the Bcl-2 family. *J. Biol. Chem.* **272**: 13829-13834.
 3. Neuchushtan, A., C.L. Smith, Y.-T. Hsu, and R.J. Youle. 1999. Conformation of the Bax C terminus regulates subcellular location and cell death. *EMBO J.* **18**: 2330-2341

Related Products:

Catalog #	Description	Size
4411-PC-100	Anti-Phosphorylated Histone H2AX polyclonal	100 μ l
6370-MC-100	Anti-human/murine-Cytochrome C	100 μ g
6380-MC-100	Anti-human/murine-Holocytochrome C	100 μ g
2291-MC-100	Anti-human-Bcl-2 mAb (clone YTH-8C8)	100 μ g
2290-MC-100	Anti-Bcl-2 mAb (clone YTH 10C4)	100 μ g
2281-MC-100	Anti-human-Bax mAb (clone YTH-6A7)	100 μ g
6361-PC-100	Anti-human/mouse-PBR polyclonal	100 μ l
4335-MC-100	Anti-PAR polymer mAb (10HA)	100 μ l
4336-BPC-100	Anti- PAR polymer polyclonal	100 μ l
4338-MC-50	Anti-human/murine-PARP mAb (clone C2-10)	50 μ g

Anti-Bcl-X_L YTH-2H12
Monoclonal Antibody
Catalog #: 2300-MC-100
Storage: 4 °C
TREVIGEN®
1-800-873-8443