

PRODUCT INFORMATION



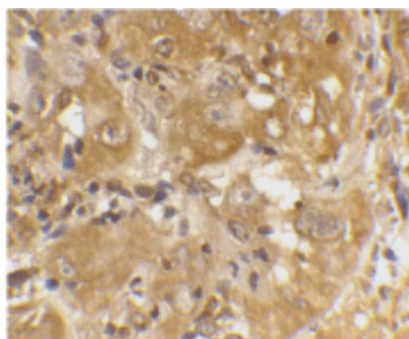
Bim/BOD (IN) Polyclonal Antibody

Item No. 10011385

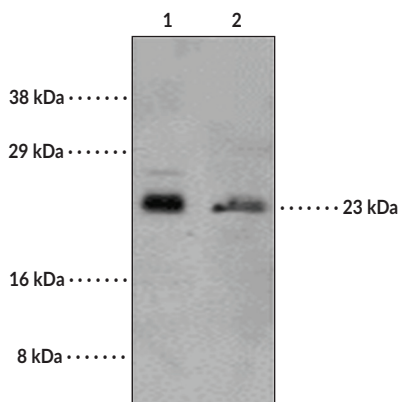
Overview and Properties

Contents:	This vial contains 25 or 100 µg protein A-affinity purified polyclonal antibody.
Immunogen:	internal central human Bim amino acids
Species Reactivity:	(+) Human, mouse, and rat Bim/BOD. Detects a 23 kDa protein corresponding to the molecular mass of Bim on SDS PAGE immunoblots.
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS, pH 7.4, with 50% glycerol and 0.09% sodium azide
Concentration:	1 mg/ml
Host:	Rabbit
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100 for IHC and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunohistochemical staining of human BIM in skin cancer cells using Bim/BOD (IN) Polyclonal Antibody at a dilution of 1:100.



Lane 1: Human K562 whole cell lysates
Lane 2: Human A549 whole cell lysates

WB of Bim/BOD (IN) Polyclonal Antibody at 1:1,000 dilution.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

5

Copyright Cayman Chemical Company, 04/25/2023

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897

[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM

WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Members in the Bcl-2 family are critical regulators of apoptosis by either inhibiting or promoting cell death. Bim/BOD is a group of three splice variants, BimEL, BimL and BimS, with apparent molecular masses of ~23, 16, and 13 kDa, respectively. Bcl-2 homology 3 (BH3) domain is a potent death domain. BH3 domain containing pro-apoptotic proteins, including Bad, Bax, Bid, Bik, and Hrk, form a growing subclass of the Bcl-2 family. A novel BH3 domain containing protein was recently identified and designated Bim or BOD in human, mouse, and rat.^{1,2} Bim/BOD interacts with diverse members in the pro-survival Bcl-2 sub-family including Bcl-2, Bcl-xL and Bcl-w. Bim/BOD induces apoptosis. The messenger RNA of Bim is ubiquitously expressed in multiple tissues and cell lines.^{1,2} Cayman's Bim/BOD (IN) Polyclonal Antibody can be used for Immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. O'Connor, L., Strasser, A., O'Reilly, L.A., *et al.* Bim: A novel member of the Bcl-2 family that promotes apoptosis. *EMBO J.* **17**, 384-395 (1998).
2. Hsu, S.Y., Lin, P., and Hsueh, A.J.W. BOD (Bcl-2-related ovarian death gene) is an ovarian BH3 domain-containing proapoptotic Bcl-2 protein capable of dimerization with diverse antiapoptotic Bcl-2 members. *Mol. Endocrinol.* **12**, 1432-1440 (1998).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM