

# Product Information



## Bim/BOD (IN) Polyclonal Antibody

Item No. 10011385

<b>Contents:</b>	This vial contains immunoaffinity chromatography purified IgG at 1 mg/ml in PBS containing 0.02% sodium azide.
<b>Antigen:</b>	Internal central human Bim amino acids
<b>Host:</b>	Rabbit
<b>Cross Reactivity:</b>	(+) human, mouse, and rat Bim/BOD. Detects a 23 kDa protein corresponding to the molecular mass of Bim on SDS PAGE immunoblots.
<b>Stability:</b>	≥1 year at -20°C
<b>Applications:</b>	The recommended starting dilution for western blot is 1:1,000 and 1:50 for immunohistochemistry. Other applications were not attempted and therefore optimal working dilutions should be determined empirically.

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.<sup>1,2</sup> 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.<sup>1,2</sup>

### 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).

### Related Products

For a list of related products please visit: [www.caymanchem.com/catalog/10011385](http://www.caymanchem.com/catalog/10011385)

**WARNING: THIS PRODUCT IS NOT FOR HUMAN OR ANIMAL DISEASE DIAGNOSIS OR THERAPEUTIC DRUG USE.**

#### MATERIAL SAFETY DATA

This material should be considered hazardous until information to the contrary becomes available. Do not ingest, swallow, or inhale. Do not get in eyes, on skin, or on clothing. Wash thoroughly after handling. This information contains some, but not all, of the information required for the safe and proper use of this material. Before use, the user must review the complete Material Safety Data Sheet, which has been sent under separate cover to the MSDS supervisor at your institution.

#### WARRANTY AND LIMITATION OF REMEDY

Cayman Chemical Company makes **no warranty or guarantee** of any kind, whether written or oral, expressed or implied, including without limitation, any warranty of fitness for a particular purpose, suitability and merchantability, which extends beyond the description of the chemicals hereof. Cayman **warrants only** to the original customer that the material will **meet our specifications at the time of delivery**.

Cayman will carry out its delivery obligations with due care and skill. Thus, in no event will Cayman have **any obligation or liability**, whether in tort (including negligence) or in contract, for any direct, indirect, incidental or consequential damages, even if Cayman is informed about their possible existence.

This limitation of liability does not apply in the case of intentional acts or negligence of Cayman, its directors or its employees. Buyer's **exclusive remedy** and Cayman's sole liability hereunder shall be limited to a **refund** of the purchase price, or at Cayman's option, the **replacement**, at no cost to Buyer, of all material that does not meet our specifications.

Said refund or replacement is conditioned on Buyer giving written notice to Cayman within thirty (30) days after arrival of the material at its destination. Failure of Buyer to give said notice within thirty (30) days shall constitute a waiver by Buyer of all claims hereunder with respect to said material.

For further details, please refer to our **Warranty and Limitation of Remedy located on our website and in our catalog.**

Copyright Cayman Chemical Company, 12/20/2011

### Cayman Chemical

#### Mailing address

1180 E. Ellsworth Road  
Ann Arbor, MI  
48108 USA

#### Phone

(800) 364-9897  
(734) 971-3335

#### Fax

(734) 971-3640

#### E-Mail

[custserv@caymanchem.com](mailto:custserv@caymanchem.com)

#### Web

[www.caymanchem.com](http://www.caymanchem.com)