



## Anti-OMI (Htr2A) (CT)

(HtrA2, High temperature requirement protein A2)

CATALOG NO.: 54171

### BACKGROUND:

Inhibitor of apoptosis proteins (IAPs) were initially identified in baculoviruses as proteins that inhibit apoptosis of the host cells to allow time for viral replication (1). Cellular homologues containing at least one baculoviral IAP repeat (BIR) motif essential for their anti-apoptosis activity have been identified in yeasts and higher organisms, often act by binding and inhibiting processed caspases (reviewed in 2). The activity of these proteins can be modulated by the expression of proteins such as Smac/DIABLO and XAF-1, which displace or prevent the binding of caspases by IAPs (reviewed in 3). Recently, a mitochondrial serine protease termed OMI/HtrA2 has been found to bind IAPs (4). Similar to Smac, OMI possesses a conserved IAP-binding motif, but acts to cleave IAPs to irreversibly inactivate IAPs and promote apoptosis (5).

### SOURCE & REACTIVITY:

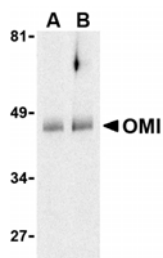
Rabbit anti-OMI polyclonal antibody was raised against a peptide corresponding to 15 amino acids near the C-terminus of human Omi (GenBank accession no. AAB94569). Anti-OMI reacts with Omi at the molecular weight of 45 kDa on western blot. Species reactivity includes human, while others are not tested.

### APPLICATION:

The following concentration ranges are recommended starting points for this product.

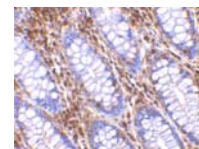
WB: 0.5 to 1 µg/ml.

### IHC



Western blot analysis of OMI in human colon tissue lysate with anti-OMI at (A) 0.5 and (B) 1 µg/ml.

Immunohistochemistry of OMI in human colon tissue with anti-OMI at 10 µg/ml.



*This product is for in vitro research purposes only.*

### RELATED PRODUCTS:

Anti-OMI (C2), Catalog No. **54126**  
Anti-Smac (CT), Catalog No. **54074**  
Anti-XIAP (CT), Catalog No. **54174**  
Anti-XAF-1 (CT), Catalog No. **54153**

### STORAGE:

Anti-OMI is supplied as ion exchange chromatography purified IgG, in 1X PBS containing 0.02% sodium azide. Store at 2-8 °C for up to 1 year. Avoid repeated freeze thaw cycles.

### REFERENCES:

1. Crook, NE. et al. *J. Virol.* **67**, 2168 (1993).
2. Liston, P. et al. *Oncogene* **22**, 8568 (2003).
3. Vaux, DL. et al. *Biochem. Biophys. Res. Comm.* **304**, 499 (2003).
4. Suzuki, Y. et al. *Mol. Cell* **8**, 613 (2001).
5. Yang, QH. et al. *Genes Dev.* **17**, 1487 (2003).

AnaSpec, Inc. • 2149 O'Toole Ave. • San Jose, CA 95131

408-452-5055 • Fax 408-452-5059 • service@anaspec.com • www.anaspec.com