

XIAP Protein

Full length recombinant protein expressed in Sf9 cells

Catalog # X08-30G

Lot # M357-1

Product Description

Recombinant full-length human XIAP was expressed by baculovirus in Sf9 insect cells using a C-terminal GST tag. The gene accession number is [NM_001167](#).

Gene Aliases

API3, ILP1, MIHA, XLP2, BIRC4

Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

Storage and Stability

Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

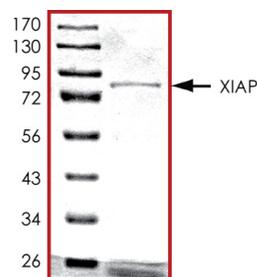
Scientific Background

XIAP is a member of a family of proteins which inhibit apoptosis and act as mammalian cell-death suppressors. XIAP inhibits at least 2 members of the caspase family of cell-death proteases, caspase-3 and caspase-7 (1). In addition, XIAP inhibits apoptosis induced by menadione, a potent inducer of free radicals, and ICE. Furthermore, XIAP can mediate protection of cells from apoptosis by utilizing both a JNK1 activation pathway that involves ILPIP and a caspase inhibition pathway that is independent of ILPIP (2). Disruption of the XIAP gene in human colon cancer cells can enhance sensitivity of these cells to exogenously added TRAIL suggesting that XIAP is a nonredundant modulator of TRAIL-mediated apoptosis.

References

1. Deveraux Q L, et al: X-linked X-linked IAP is a direct inhibitor of cell-death proteases. Nature 388: 300-304, 1997.
2. Tang G, et al: Inhibition of JNK activation through NF- κ B target genes. Nature 414: 313-317, 2001.

Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **84kDa**.

XIAP Protein

Full length recombinant protein expressed in Sf9 cells

| | |
|---------------------|--------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------------|
| Catalog Number | X08-30G |
| Specific Lot Number | M357-1 |
| Purity | >70% |
| Concentration | 0.2 µg/µl |
| Stability | 1yr at -70°C from date of shipment |
| Storage & Shipping | Store product at -70°C . For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice. |

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: orders@signalchem.com
www.signalchem.com

FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.