

Data Sheet

XIAP (Bir2, Bir3)

Human, recombinant, N-terminal His-tag
Catalog #: 75002
Lot#: 120918 **Conc.:** 0.45 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, and 20% glycerol.

Stability: >6 months at -80 °C

References:

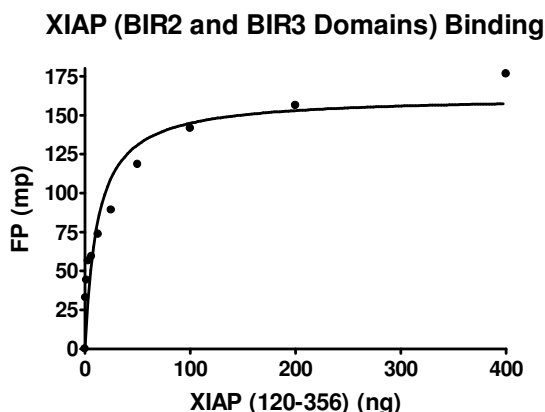
1. Nikolovska-Coleska, Z., *et al.*, *Anal Biochem.* 2008 Mar 1;**374(1)**:87-98.
2. Verhagen, A.M., *et al.*, *Cell Death Differ.* 2007 Feb;**14(2)**:348-57.
3. Fulda, S., Vucic., D., *Nat Rev Drug Discov.* 2012 Feb 1;**11(2)**:109-24.

Description: Human x-linked inhibitor of apoptosis (XIAP), also known as IAP-like protein, API3, X-linked IAP, and BIRC4, amino-acids 120 – 356 containing Bir2 and Bir3 domains (GenBank Accession No. NM_001167) with N-terminal His-tag, MW = 28.3 kDa, expressed in an *E. coli* expression system.

Assay Conditions: 50 mM Tris-HCl, pH 7.4, 25 mM NaCl, 0.05% Tween 20, 0.01% BSA, 0.5 nM FAM-SMAC peptide. Incubate 30 minutes at room temperature. Fluorescence polarization is measured at λ_{exc} 475 nm, λ_{em} 528 nm.

Application: Useful in binding and inhibition studies of Caspases 3, 7, and 9, screening inhibitors, and selectivity profiling.

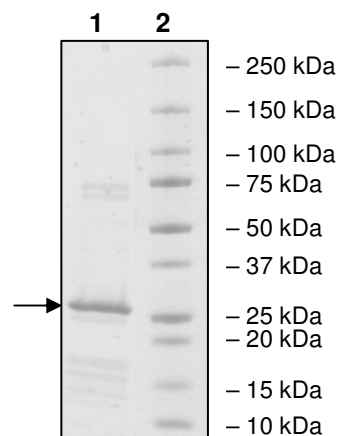
Quality Assurance



4-20% SDS-PAGE
Coomassie staining

Lane 1:
1.8 µg XIAP
Lane 2:
Protein Marker

MW: 28.3 kDa
Purity: ≥60%



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
 To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
 Or you can Email us at: info@bpsbioscience.com
 Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence:

MHHHHHHYLGSRDHFALDRPSETHADYLLRRTGQVVDISDTIYPRNPAMYSEEARLKSF
QNWPDYAHLTPRELASAGLYYTGIGDQVQCFCGGLKNWEPDRAWSEHRRHFPN
CFFVLGRNLNIRSEDAVSSDRNFPNSTNLPRNPSMADYEARIFTFGTWIYSVNKEQLA
RAGFYALGEGDKVKCFHCGGLTDWKPSDPWEQHAKWYPGCKYLLEQKGQEYINNI
HLTHSLEECLVRTT

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.
To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com