

## Data Sheet

### **XIAP (Bir3)**

Human, recombinant, N-terminal His-tag  
**Catalog #:** 75001  
**Lot#:** 120718      **Conc.:** 2.42 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, 3 mM DTT and 20% glycerol.

**Stability:** >6 months at -80 °C

**References:**

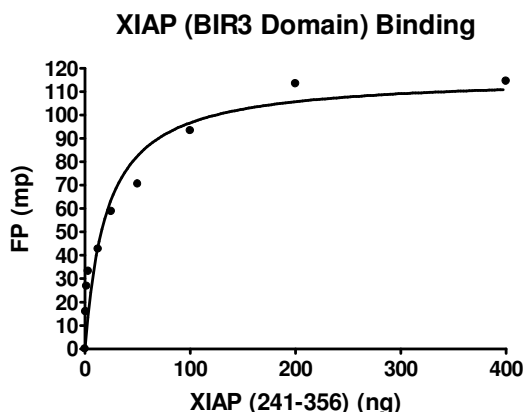
1. Nikolovska-Coleska, Z., *et al.*, *Anal Biochem.* 2008 Mar 1;**374(1)**:87-98.
2. Verhagen, A.M., *et al.*, *Cell Death Differ.* 2007 Feb;**14(2)**:348-57.
3. Fulda, S., Vucic., D., *Nat Rev Drug Discov.* 2012 Feb 1;**11(2)**:109-24.

**Description:** Human x-linked inhibitor of apoptosis (XIAP) Bir3 domain, also known as IAP-like protein, API3, X-linked IAP, and BIRC4, amino-acids 241 – 356 (GenBank Accession No. NM\_001167) with N-terminal His-tag, MW = 14.2 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  $\lambda_{exc}$  475 nm,  $\lambda_{em}$  528 nm.

**Application:** Useful in binding and inhibition studies of Caspase 9, screening inhibitors, and selectivity profiling.

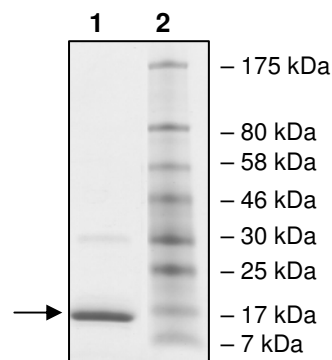
### Quality Assurance



**4-20% SDS-PAGE  
 Coomassie staining**

**Lane 1:**  
 7 µg XIAP  
**Lane 2:**  
 Protein Marker

**MW:** 14.2 kDa  
**Purity:** ≥92%



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](mailto:info@bpsbioscience.com)  
 Please visit our website at: [www.bpsbioscience.com](http://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](mailto:info@bpsbioscience.com)

**Sequence:**

MHHHHHSDAVSSDRNFPNSTNLPRNPSMADYEARIFTFGTWIYSVNKEQLARAGFYA  
LGEGDKVKCFHCGGLTDWKPSDPWEQHAKWYPGCKYLLEQKGQEYINNIHLTHSL  
EECLVRTT

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](mailto:info@bpsbioscience.com)  
Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)