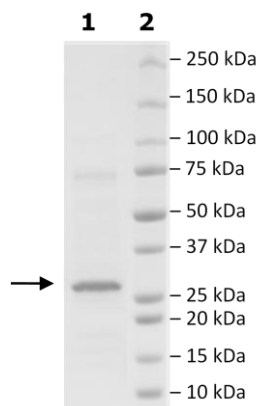


Product Information

Description:	Recombinant human XIAP (x-linked inhibitor of apoptosis protein), encompassing amino acids 120-356, corresponding to the Bir2-Bir3 (baculovirus IAP repeat) domains. This construct contains an N-terminal His-tag (6xHis). The protein was affinity purified.
Background:	The XIAP (X-linked inhibitor of apoptosis protein), also known as IAP3 (inhibitor of apoptosis protein 3) and BIRC4 (baculoviral IAP repeat-containing protein 4), is a member of the IAP family of proteins. XIAP directly regulates caspases and suppresses apoptotic cell death pathways. XIAP has three Bir (baculovirus IAP repeat) domains, involved in binding to caspases. Bir2 inhibits caspase 3 and 7, while Bir3 is involved in caspase 9 inhibition. An increased expression level of XIAP has been observed in many cancer types, such as prostate cancer. Dysfunction of XIAP has been linked to inflammatory bowel disease, cancer, neurodegenerative and autoimmunity disorders.
Species:	Human
Construct:	XIAP (His-120-356)
Concentration:	0.98 mg/ml
Expression System:	<i>E. coli</i>
Purity:	80%
Format:	Aqueous buffer solution
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole, 20% glycerol, and 3 mM DTT
MW:	28 kDa
Genbank Accession:	NM_001167
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	The assay was performed in 50 mM Tris-HCl, pH 7.4, 25 mM NaCl, 0.05% Tween-20, 0.01% BSA, with 0.5 nM FAM-SMAC peptide (Anaspec #SA-64758), incubated 30 minutes at room temperature. Fluorescence polarization was measured at $\lambda_{exc}/\lambda_{em}$ = 475 nm/528 nm.
Applications:	Useful for binding studies.

Quality Control Data

4-20% SDS-Page Coomassie Staining



XIAP Bir2-Bir3 Domains Binding Activity

