

**XIAP BIR2 domain (His tag)**

**CATALOG NO.:** APT-11-491

**LOT NO.:**

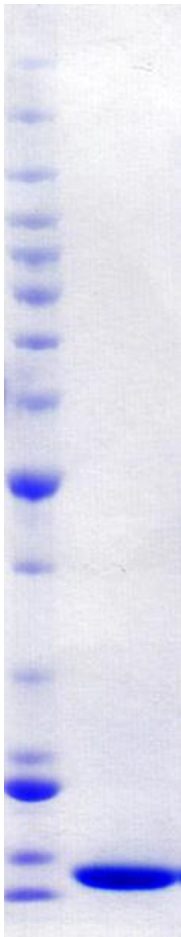
**DESCRIPTION:** Human recombinant BIR2 domain of XIAP (residues 152-236; NCBI Reference Sequence NM\_001167.3; MW = 13.59 kDa) expressed in *E. coli* with an N-terminal 6xHis-tag and C-terminal Strep-tag.

**PURITY:** >90% by SDS-PAGE

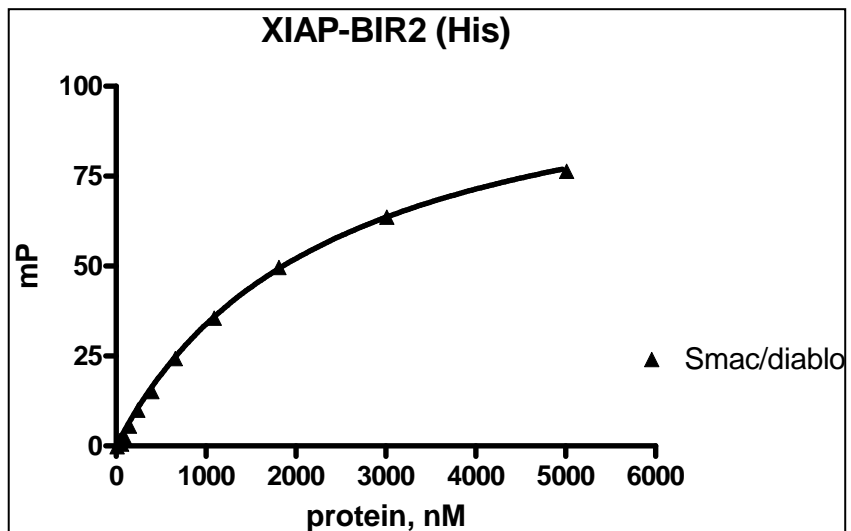
**ASSAY CONDITIONS:** RBC's XIAP-BIR2 (His) domain displays binding affinity for Smac/DIABLO substrate peptide. Reactions were performed with 25 nM FAM-labelled peptide at room temperature in 100mM Potassium Phosphate, pH 7.5, 0.1mg/mL BSA. The resulting polarization values (Ex. 480nm/Em. 535nm) was read following a 60 minute incubation. (See figure, below.)

**SUPPLIED AS:** \_\_\_ µg/µL in 50 mM Tris pH 7.5, 500 mM NaCl, 10% glycerol, 1 mM TCEP as determined by OD<sub>280</sub>.

**STORAGE:** -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted enzyme should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC XIAP-BIR2 (His).** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**XIAP-BIR2 (His) Binding Assay.** FAM-Smac/DIABLO substrate peptide binding was detected using fluorescence polarization. The 20µL reaction contained 25nM peptide and variable concentration of XIAP-BIR2 (His). Polarization values (mP) were measured at an excitation wavelength at 480nm and an emission wavelength at 535nm using a ClarioStar reader (BMG Labtech) following a 60 minute incubation.

This product is NOT intended for therapeutic or diagnostic use in animals or in humans.