

cIAP1-BIR3 (GST)

CATALOG NO.: APT-11-371

LOT NO.:

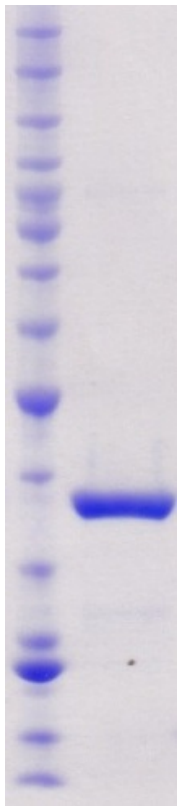
DESCRIPTION: Human recombinant cIAP-BIR3 (residues 258-363; Genbank Accession # NM_001166.4; MW = 40.3 kDa) expressed with an N-terminal GST tag and C-terminal StrepII tag in *E.coli*.

PURITY: >95% by SDS-PAGE

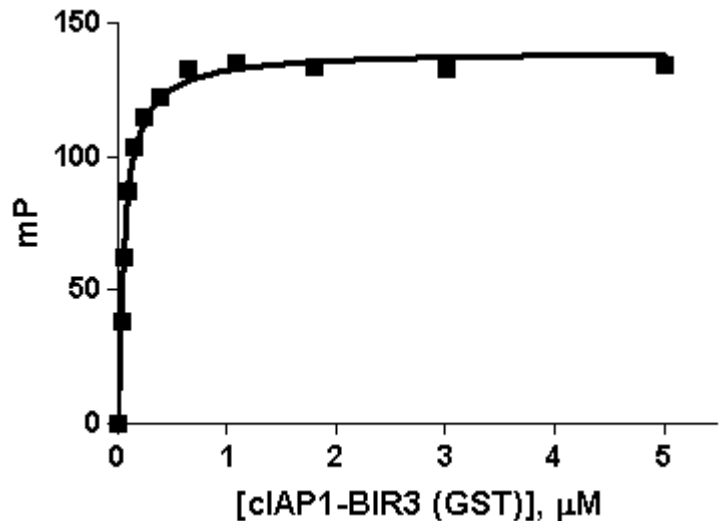
ASSAY CONDITIONS: RBC's cIAP1-BIR3 domain displays binding affinity for Smac/DIABLO peptide. Reactions were performed with 20nM FAM-labelled peptide at room temperature in 50mM HEPES, pH 7.5, 200mM NaCl, 0.01% Triton X100. The resulting polarization values (Ex. 480 nm/Em. 535 nm) was read following a 30 minute incubation. (See figure, below.)

SUPPLIED AS: _ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein 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 (12% acrylamide) of 4µg of RBC cIAP1-BIR3 (nGST, cStrepII). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



cIAP1-BIR3 Binding Assay. FAM-Smac/DIABLO peptide binding was detected using fluorescence polarization. The 20µL reaction contained 20nM peptide and variable concentration of cIAP1-BIR3. Polarization values (mP) were measured at an excitation wavelength at 480 nm and an emission wavelength at 535 nm using an EnVision reader (Perkin Elmer) following a 30 minute incubation.

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