

Bcl-XL (GST)

CATALOG NO.: APT-11-442

LOT NO.:

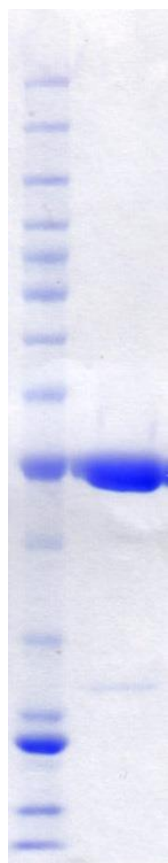
DESCRIPTION: Apoptosis regulator Bcl-XL (residues 1-209; Genbank Accession # NM_138578; MW = 49.8 kDa) expressed with a C-terminal GST tag in *E.coli*.

PURITY: >90% by SDS-PAGE

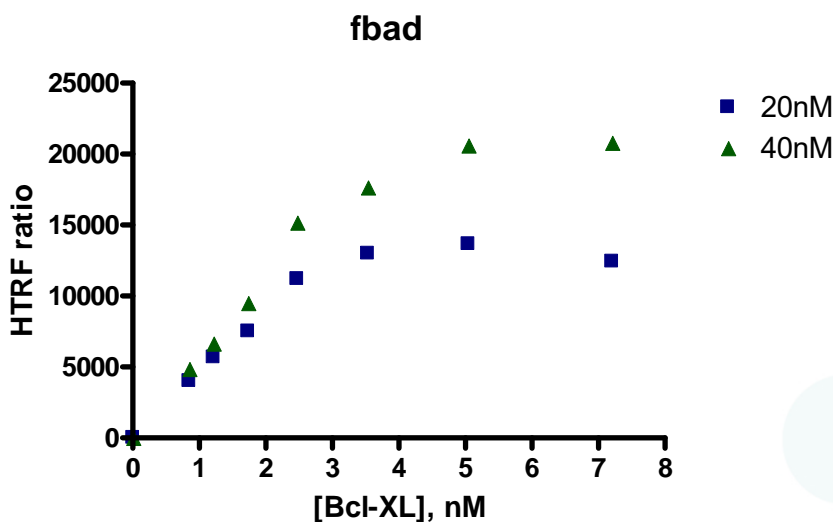
ASSAY CONDITIONS: RBC's Bcl-XL domain displays binding affinity for fbad peptide in TR-FRET based assay. Reactions were performed with the indicated FAM-labeled peptide at room temperature in 50mM potassium phosphate, pH 7.5, 50mM NaCl, 1mM EDTA, 0.01% NP-40 and 0.5% DMSO. The resulting fluorescence emission (520 and 490 nm) was read following 60 minute incubation with detection mixture. (See figure, below.)

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris-HCl pH 8, 500 mM NaCl, 10% glycerol, 1 mM TCEP as determined by OD₂₈₀

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 (4-12% acrylamide) of 4 μg of RBC Bcl-XL (GST). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Bcl-XL/peptide interaction assay. Bcl-XL binding to FAM-labeled bad peptide was assessed using Cisbio HTRF detection (cat # 61GSTTLA). The 15 μL reaction contained variable protein, fixed peptide concentrations as indicated above and detection mix. Fluorescence emission (520 and 490nm) was read using Envision reader (Perkin Elmer) following 1h incubation.

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