

**Bcl-2 (GST)**

**CATALOG NO.:** APT-11-441

**LOT NO.:**

**DESCRIPTION:** Apoptosis regulator Bcl-2 (residues 1-207; Genbank Accession # NM\_000633; MW = 49.2 kDa) expressed with a C-terminal GST tag in *E.coli*.

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

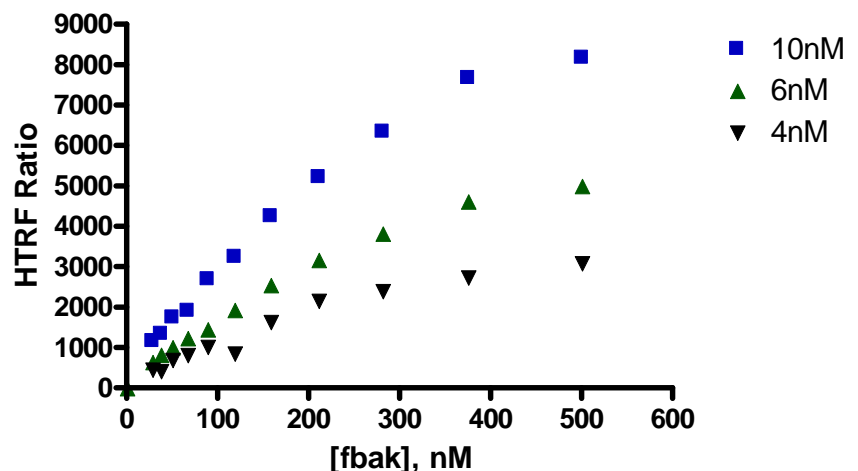
**ASSAY CONDITIONS:** RBC's Bcl-2 domain displays binding affinity for fbak peptide in TR-FRET based assay. Reactions were performed with variable concentration of FAM-labeled peptide at room temperature in 20mM potassium phosphate, pH 7.5, 50mM NaCl, 1mM EDTA, and 0.01% NP-40. The resulting fluorescence emission (520 and 490 nm) was read following 60 minute incubation with detection mix (see figure, below).

**SUPPLIED AS:** \_ µg/µL in 50 mM Tris-HCl 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 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-2 (GST). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



**Bcl-2/peptide interaction assay.** Bcl-2 binding to FAM-labeled bak peptide was assessed using Cisbio HTRF detection (cat # 61GSTTLA). The 15µL reaction contained the above indicated fixed protein concentration, variable peptide concentration 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.