

**Bcl-2 (G101V) (GST)**
**CATALOG NO.:** APT-11-633

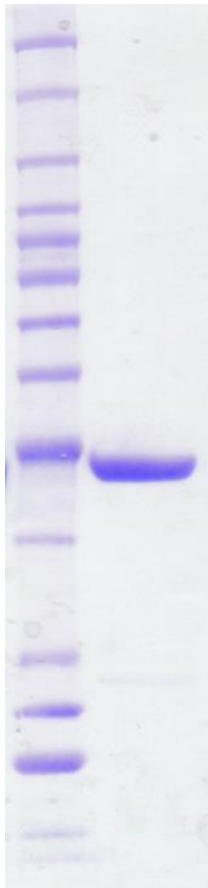
**LOT NO.:**

**DESCRIPTION:** Mutant apoptosis regulator Bcl-2 (G101V) with valine (V) substituted for glycine-101 (G) (otherwise contains wildtype residues 1-207; Genbank Accession # NM\_000633.3; MW = 49.3 kDa) expressed with a C-terminal GST fusion protein in *E. coli*.

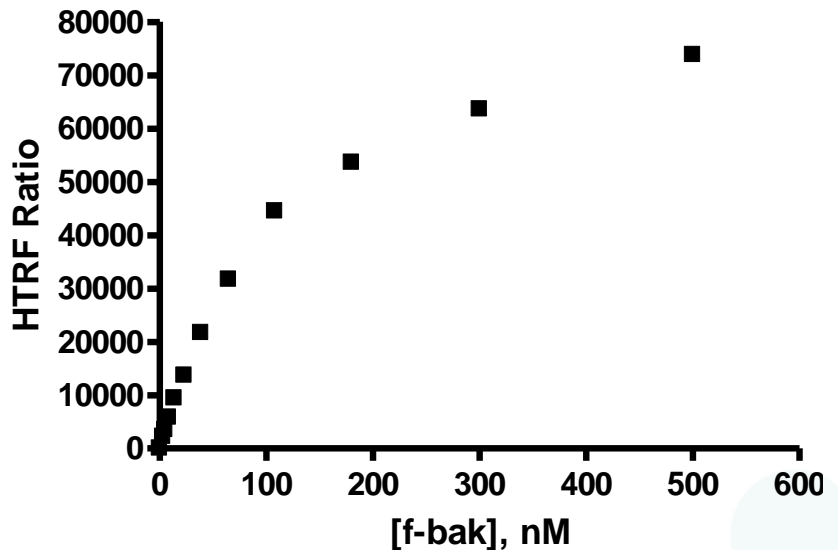
**PURITY:** > 95% by SDS-PAGE.

**SUPPLIED AS:** \_\_\_  $\mu\text{g}/\mu\text{L}$  in 50 mM Tris, pH 7.5, 500 mM NaCl, 10% glycerol, 0.25 mM TCEP as determined by  $\text{OD}_{280}$ .

**STORAGE:**  $-70^{\circ}\text{C}$ . 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  $\mu\text{l}$ ) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 2  $\mu\text{g}$  of RBC Bcl-2 (G101V) (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 15  $\mu\text{L}$  reaction contained 4 nM protein, variable peptide concentration and detection mix. Fluorescence emission (520 and 490 nm) was read using PHERAstar reader (BMG Labtech) following 1h incubation.

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

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com