

BCL2

(Apoptosis regulator Bcl-2)

- **CAT#:** BCL2
- **Enzyme:** Human BCL2
- **Substrate:** Bak-BH3 peptide, 5 nM
- **Reaction:**
Competition binding assay with Bak-BH3 peptide.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
ABT-737	5.1
ABT-263	11.6

- **Scientific Information:** P10415 (<http://www.expasy.org/uniprot/P10415>)

Bak-BH3 peptide displacement IC50 for BCL2

