

BCLXL

(Bcl-2-like protein 1, Bcl2-L-1, Apoptosis regulator Bcl-X, BCL2L1, BCL2L, BCLX, Bcl-X(L))

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

Compounds	IC50 (nM)
ABT-737	2.8
ABT-263	6.0

- **Scientific Information:** Q07817 (<http://www.expasy.org/uniprot/Q07817>)

Bak-BH3 peptide displacement IC50 for BCLXL

