

HSP90b

(Heat shock protein HSP 90-beta, Heat shock 84 kDa, HSP90AB1, HSP90B, HSPC2, HSPCB, HS90B)

- **CAT#:** HSP90b
- **Enzyme:** Human HSP90b
- **Probe:** Geldanamycin, 5 nM
- **Reaction:**
Competition binding assay with fluorescently labeled Geldanamycin.
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Geldanamycin	50.3
Radicicol	89.0

- **Scientific Information:** P08238 (<http://www.expasy.org/uniprot/P08238>)

Control Compound IC50 for HSP90b

