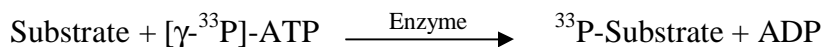


AKT2 (Δ PH, S474D)

(RAC-beta serine/threonine-protein kinase, RAC-PK-beta, Protein kinase Akt-2, Protein kinase B beta, PKB β)

- **CAT#:** AKT2 (Δ PH, S474D)
- **Enzyme:** Human AKT2 (Δ PH, S474D)
- **Substrate:** Peptide substrate, [KGSGSGRPRTSSFAEG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.8
Ro 31-8220	73
GW 5074	7,200
H-89	280
Rottlerin	162,000

- **Scientific Information:** P31751 (<http://www.expasy.org/uniprot/P31751>)

Reference compound IC50 for Akt2

