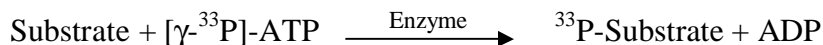


AKT3 (Δ PH, S472D)

(RAC-gamma serine/threonine-protein kinase, RAC-PK-gamma, Protein kinase Akt-3, Protein kinase B gamma, PKB gamma, STK-2)

- **CAT#:** AKT3 (Δ PH, S472D)
- **Enzyme:** Human AKT3 (Δ PH, S472D)
- **Substrate:** Peptide substrate, [KGSGRPRTSSFAEG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.2
Ro 31-8220	12.6
GW 5074	10,400
H-89	525
Rottlerin	304,000

- **Scientific Information:** Q9Y243 (<http://www.expasy.org/uniprot/Q9Y243>)

Reference compound IC50 for AKT3

