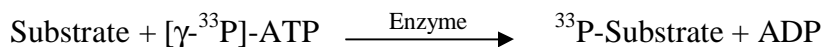


AKT1 (Δ PH, S473D)

(RAC-alpha serine/threonine-protein kinase,
RAC-PK-alpha, Protein kinase B α , C-AKT)

- **CAT#:** AKT1 (Δ PH, S473D)
- **Enzyme:** Human AKT1 (Δ PH, S473D)
- **Substrate:** Peptide substrate, [KGS γ SGSRPRTSSFAEG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.6
Ro 31-8220	208
GW 5074	16,000
H-89	290
H-89	290

- **Scientific Information:** P31749 (<http://www.expasy.org/uniprot/P31749>)

Reference compound IC50 for AKT1

