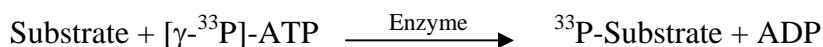


## GSK3b

(Glycogen synthase kinase-3 beta)

- **CAT#:** GSK3b
- **Enzyme:** Human GSK3b
- **Substrate:** Peptide substrate, [YRRAAVPPSPSLSRHSSPHQ(pS)EDEEE], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	11.5
Ro 31-8220	2.5
GW 5074	970
Rottlerin	6,700

- **Scientific Information:** P49841 (<http://www.expasy.org/uniprot/P49841>)

### Reference compound IC50 for GSK3b

