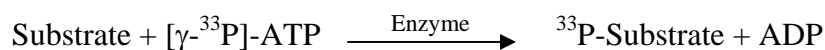


## IKKb/IKBKB

(Inhibitor of nuclear factor kappa-B kinase subunit beta, I-kappa-B-kinase beta, IkbKB, IKK-beta, IKK-B, I-kappa-B kinase 2, IKK2, Nuclear factor NF-kappa-B inhibitor kinase beta, NFKB1KB, IKBKB, IKKB)

- **CAT#:** IKKb/IKBKB
- **Enzyme:** Human IKKb/IKBKB
- **Substrate:** Peptide substrate, [KKKKERLLDDRHDSGLDSMKDEE], 20 μM  
ATP 10 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6,600

- **Scientific Information:** O14920 (<http://www.expasy.org/uniprot/O14920>)

### Reference compound IC50 for IKKb/IKBKB

