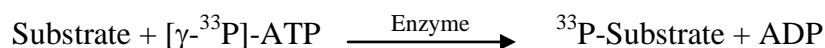


EIF2AK2

(Interferon-induced, double-stranded RNA-activated protein kinase, Interferon-inducible RNA-dependent, protein kinase, Eukaryotic translation initiation factor 2-alpha kinase 2, eIF-2A protein kinase 2, Protein, kinase RNA-activated, PKR, p68 kinase, P1/eIF-2A protein kinase, PKR, PRKR, E2AK2)

- **CAT#:** EIF2AK2
- **Enzyme:** Human EIF2AK2
- **Substrate:** Peptide substrate, [H-GRSRSRSRSR-OH], 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	39
Ro 31-8220	5,500
GW 5074	3,900

- **Scientific Information:** P19525 (<http://www.expasy.org/uniprot/P19525>)

Reference compound IC50 for EIF2AK2/RStide

