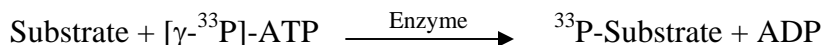


## EPHA7

(Ephrin type-A receptor 7 [Precursor], Tyrosine-protein kinase receptor EHK-3, EPH homology kinase 3, Receptor protein-tyrosine kinase HEK11, EHK3, HEK11)

- **CAT#:** EPHA7
- **Enzyme:** Human EPHA7
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	17.8
PP2	2,500
AG1478	481
D-64406	77,500

- **Scientific Information:** Q15375 (<http://www.expasy.org/uniprot/Q15375>)

### Reference compound IC50 for EPHA7

