

## EPHA4

(Ephrin type-A receptor 4 [Precursor], Tyrosine-protein kinase receptor SEK, Receptor protein-tyrosine kinase HEK8)

- **CAT#:** EPHA4
- **Enzyme:** Human EPHA4
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 1  $\mu$ M
- **Reaction:**  

$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	9.4
PP2	44
AG1478	924
PDGFR I-III	1430,000
D-64406	6,500

- **Scientific Information:** P54764 (<http://www.expasy.org/uniprot/P54764>)

### Reference compound IC50 for EPHA4

