

EPHA5

(Ephrin type-A receptor 5 [Precursor], Tyrosine-protein kinase receptor EHK-1, EPH homology kinase 1, Receptor protein-tyrosine kinase HEK7)

- **CAT#:** EPHA5
- **Enzyme:** Human EPHA5
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ 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	14.5
PP2	19
AG1478	612
D-64406	9,000

- **Scientific Information:** P54756 (<http://www.expasy.org/uniprot/P54756>)

Reference compound IC50 for EPHA5

