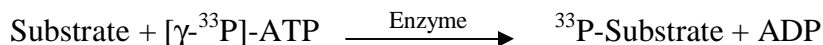


EPHA3

(Ephrin type-A receptor 3 [Precursor], Tyrosine-protein kinase receptor ETK1, HEK, HEK4)

- **CAT#:** EPHA3
- **Enzyme:** Human EPHA3
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	18
PP2	368
AG1478	3,400
PDGFR I-III	72,000
D-64406	16,300

- **Scientific Information:** P29320 (<http://www.expasy.org/uniprot/P29320>)

Reference compound IC50 for EPHA3

