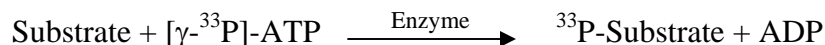


EPHA8

(Ephrin type-A receptor 8 [Precursor], Tyrosine-protein kinase receptor EEK
EPH- and ELK-related kinase, HEK3, EEK, KIAA1459)

- **CAT#:** EPHA8
- **Enzyme:** Human EPHA8
- **Substrate:** Peptide substrate, , poly[Glu:Tyr] (4:1), 0.1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	32
PP2	31.4
AG1478	1,600
D-64406	43,300

- **Scientific Information:** P29322 (<http://www.expasy.org/uniprot/P29322>)

Reference compound IC50 for EPHA8

