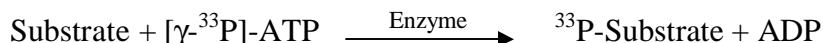


AXL

(Tyrosine-protein kinase receptor UFO [Precursor], AXL oncogene, UFO)

- **CAT#:** AXL
- **Enzyme:** Human AXL
- **Substrate:** Peptide substrate, [KKSARGDYMTMQIG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	90
AG1478	25,000
D-64406	17,600

- **Scientific Information:** P30530 (<http://www.expasy.org/uniprot/P30530>)

Reference Compound IC50 for AXL

