

AXL

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

- **CAT#:** AXL
- **Enzyme:** Human AXL
- **Substrate:** Peptide substrate, [EAIYAAPFAKKK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.2
PP2	54,000
AG1478	8,800
PDGFR I-III	60,000
D-64406	13,000

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

Reference compound IC50 for AXL

