

ALK (R1275Q)

(ALK tyrosine kinase receptor [Precursor], Anaplastic lymphoma kinase)

- **CAT#:** ALK (R1275Q)
- **Enzyme:** Human ALK (R1275Q)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.5
PP2	40,000
AG1478	6,000
PDGFR I-III	>100,000
D-64406	24,000

- **Scientific Information:** Q9UM73 (<http://www.expasy.org/uniprot/Q9UM73>)

Reference compound IC50 for ALK (R1275Q)

