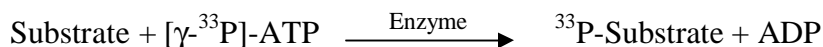


## ALK

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

- **CAT#:** ALK
- **Enzyme:** Human ALK
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	16
PP2	90,000
AG1478	9,300

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

