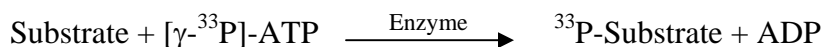


## JAK1

(Tyrosine-protein kinase JAK1, Janus kinase 1, JAK-1, JAK1A, JAK1B)

- **CAT#:** JAK1
- **Enzyme:** Human JAK1
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	< 0.1
AG1478	14,200
D-64406	11,800

- **Scientific Information:** P23458 (<http://www.expasy.org/uniprot/P23458>)

### Reference compound IC50 for JAK1

