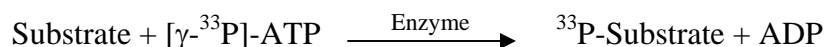


## ZAP70

(Tyrosine-protein kinase ZAP-70, 70 kDa zeta-associated protein, Syk-related tyrosine kinase, SRK)

- **CAT#:** ZAP70
- **Enzyme:** Human ZAP70
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.9

- **Scientific Information:** P43403 (<http://www.expasy.org/uniprot/P43403>)

### Reference compound IC50 for ZAP70

