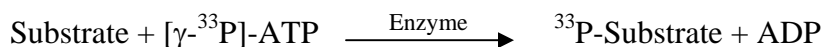


BLK

(Tyrosine-protein kinase BLK, B lymphocyte kinase, p55-BLK)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	32.7
PP2	379
AG1478	12,000
D-64406	4,100

- **Scientific Information:** P51451 (<http://www.expasy.org/uniprot/P51451>)

Reference Compound IC50 for BLK

