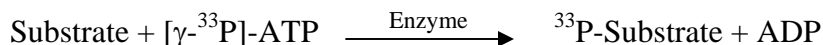


BTK (E41K)

(Tyrosine-protein kinase BTK, Bruton tyrosine kinase, Agammaglobulinaemia tyrosine kinase, ATK, B cell progenitor kinase, BPK)

- **CAT#:** BTK (E41K)
- **Enzyme:** Human BTK (E41K)
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	9.0
PP2	233
AG1478	24,400
D-64406	22,000

- **Scientific Information:** Q06187 (<http://www.expasy.org/uniprot/Q06187>)

Reference Compound IC50 for BTK (E41K)

