

## BTK

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

- **CAT#:** BTK
- **Enzyme:** Human BTK
- **Substrate:** Peptide substrate, [KVEKIGEGTYGVVYK], 20  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	16
PP2	11
AG1478	1,600
D-64406	4,900

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

### Reference compound IC50 for BTK

