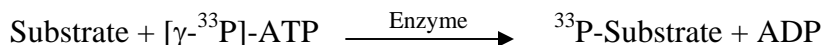


## BRK

(Tyrosine-protein kinase 6, Breast tumor kinase  
Tyrosine-protein kinase BRK, PTK6)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	270
PP2	6.8
AG1478	103
D-64406	12,500

- **Scientific Information:** Q13882 (<http://www.expasy.org/uniprot/Q13882>)

### Reference compound IC50 for BRK

