

BMX/ETK

(Cytoplasmic tyrosine-protein kinase BMX, Bone marrow tyrosine kinase gene in chromosome X protein, Epithelial and endothelial tyrosine kinase, ETK, NTK38)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	10
PP2	62
AG1478	7,400
D-64406	14,000

- **Scientific Information:** P51813 (<http://www.expasy.org/uniprot/P51813>)

Reference compound IC50 for BMX/ETK

