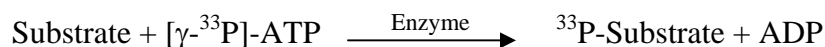


BMPR2

(Bone morphogenetic protein receptor type-2, PPH1)

- **CAT#:** BMPR2
- **Enzyme:** Human BMPR2
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	670
Ro 31-8220	2,000
GW5074	2,000
H-89	>50,000
Rottlerin	>50,000

- **Scientific Information:** Q13873 (<http://www.expasy.org/uniprot/Q13873>)

Reference compound IC50 for BMPR2

