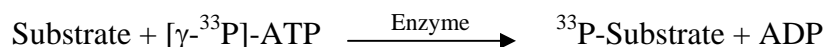


TGFβR2

(TGF-beta receptor type-2 [Precursor], Transforming growth factor-beta receptor type II, TGF-beta receptor type II, TGF-beta type II receptor, TbetaR-II, TGFR-2, TGFBR2, TGFR2)

- **CAT#:** TGFβR2
- **Enzyme:** Human TGFβR2
- **Substrate:** MBP (Myelin Basic Protein), 10 μM
ATP 1 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	6,600
Ro 31-8220	>10,000
GW 5074	3,500
H-89	>10,000
Rottlerin	43,000

- **Scientific Information:** P37173 (<http://www.expasy.org/uniprot/P37173>)

Reference Compound IC50 for TGFβR2

