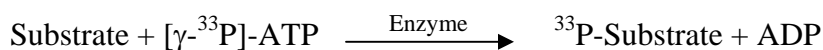


ALK5/TGFBR1

(TGF-beta receptor type-1 [Precursor], TGFR-1, TGF-beta type I receptor, Transforming growth factor-beta receptor type I, TbetaR-I, Serine/threonine-protein kinase receptor R4, SKR4, Activin receptor-like kinase 5, TGFR1)

- **CAT#:** ALK5
- **Enzyme:** Human ALK5
- **Substrate:** Casein, 1 mg/ml
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1,900
Ro 31-8220	10,000
GW 5074	6,800
H-89	93,400

- **Scientific Information:** P36897 (<http://www.expasy.org/uniprot/P36897>)

Reference Compound IC50 for ALK5/TGFB-R1

