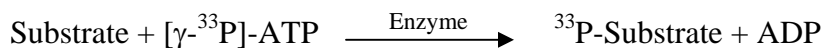


ALK2/ACVR1

(Activin receptor type-1 [Precursor], Activin receptor type I, ACTR-I, Serine/threonine-protein kinase receptor R1, SKR1, Activin receptor-like kinase 2, ALK-2 TGF-B superfamily receptor type I, TSR-I, ACVRLK2)

- **CAT#:** ALK2/ACVR1
- **Enzyme:** Human ALK2/ACVR1
- **Substrate:** Casein, 1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4,800
GW 5074	26,000

- **Scientific Information:** Q04771 (<http://www.expasy.org/uniprot/Q04771>)

Reference Compound IC50 for ALK2/ACVR1

