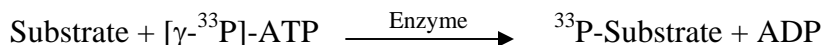


ALK4/ACVR1B

(Activin receptor type 1B [Precursor], ACTR-IB, Serine/threonine-protein kinase receptor R2 SKR2, Activin receptor-like kinase 4, ALK-4, ACVR1B)

- **CAT#:** ALK4/ACVR1B
- **Enzyme:** Human ALK4/ACVR1B
- **Substrate:** Casein, 1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3,700
GW 5074	4,300

- **Scientific Information:** P36896 (<http://www.expasy.org/uniprot/P36896>)

Reference compound IC50 for ALK4/ACVR1B

