

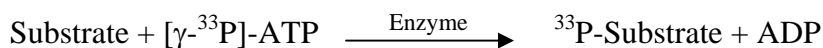
TAK1

(TAK1: Mitogen-activated protein kinase kinase kinase 7, Transforming growth factor-beta-activated kinase 1, TGF-beta-activated kinase 1, M3K7, MAP3K7)

TAB1: Mitogen-activated protein kinase kinase kinase 7-interacting protein 1, TGF-beta-activated kinase 1-binding protein 1, TAK1-binding protein 1, MAP3K7IP1)

- **CAT#:** TAK1
- **Enzyme:** Human TAK1
- **Substrate:** Casein, 1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	36
Ro 31-8220	16,700
GW 5074	14,000

- **Scientific Information:** O43318 (<http://www.expasy.org/uniprot/O43318>)

Reference compound IC50 for TAK1

