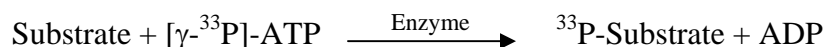


LATS1

(Serine/threonine-protein kinase LATS1, Large tumor suppressor homolog 1, WARTS protein kinase)

- **CAT#:** LATS1
- **Enzyme:** Human LATS1
- **Substrate:** Peptide substrate, [KKRNRRLSVA], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.5
Ro 31-8220	10
GW 5074	14,000
H-89	500
Rottlerin	>50,000

- **Scientific Information:** O95835 (<http://www.expasy.org/uniprot/Q95835>)

Reference compound IC50 for LATS1

