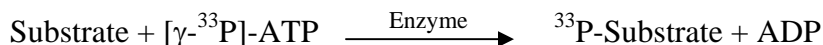


LATS2

(Serine/threonine-protein kinase LATS2, Kinase phosphorylated during mitosis protein, Large tumor suppressor homolog 2, Serine/threonine-protein kinase kpm, Warts-like kinase, KPM)

- **CAT#:** LATS2
- **Enzyme:** Human LATS2
- **Substrate:** Peptide substrate, [KKRNRRLSVA], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.6
Ro 31-8220	15
GW 5074	1,700
H-89	600
Rottlerin	>50,000

- **Scientific Information:** Q9NRM7 (<http://www.expasy.org/uniprot/Q9NRM7>)

Reference compound IC50 for LATS2

