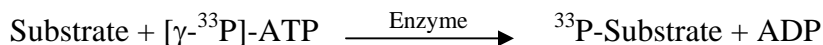


LIMK1

(LIM domain kinase 1)

- **CAT#:** LIMK1
- **Enzyme:** Human LIMK1
- **Substrate:** Cofilin 1, 3 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.9
Ro 31-8220	17,000
GW 5074	1,500
H-89	2,400
Rottlerin	3,900

- **Scientific Information:** P53667 (<http://www.expasy.org/uniprot/P53667>)

Reference compound IC50 for LIMK1

