

MSSK1

(Serine/threonine-protein kinase SRPK3, Serine/arginine-rich protein specific kinase 3, Serine/threonine-protein kinase 23, Muscle-specific serine kinase 1, SRPK3, STK23)

- **CAT#:** MSSK1
- **Enzyme:** Human MSSK1
- **Substrate:** Peptide substrate, [GRSRSRSRSR], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1,750
Ro 31-8220	540
GW 5074	5,800
Rottlerin	11,000

- **Scientific Information:** Q9UPE1 (<http://www.expasy.org/uniprot/Q9UPE1>)

Reference Compound IC50 for MSSK1

