

VRK1

(Serine/threonine-protein kinase VRK1, Vaccinia-related kinase 1)

- **CAT#:** VRK1
- **Enzyme:** Human VRK1
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	8,400
Ro 31-8220	2,300
GW 5074	57,000
H-89	61,000

- **Scientific Information:** Q99986 (<http://www.expasy.org/uniprot/Q99986>)

Reference Compound IC50 for VRK1

