

GRK7

(G protein-coupled receptor kinase 7 [Precursor], G protein-coupled receptor kinase GRK7, GPRK7)

- **CAT#:** GRK7
- **Enzyme:** Human GRK7
- **Substrate:** Casein, 1 mg/ml
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.5
Ro 31-8220	1,500
GW5074	40,900
H-89	15,000

- **Scientific Information:** Q8WTQ7 (<http://www.expasy.org/uniprot/Q8WTQ7>)

Reference compound IC50 for GRK7

