

ROS/ROS1

(Proto-oncogene tyrosine-protein kinase ROS [Precursor], c-ros-1)

- **CAT#:** ROS/ROS1
- **Enzyme:** Human ROS/ROS1
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 20 μ M
ATP 10 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	<0.1
PP2	3,100
AG1478	2,600
PDGFR I-III	>20,000
D-64406	>20,000

- **Scientific Information:** P08922 (<http://www.expasy.org/uniprot/P08922>)

Reference compound IC50 for ROS/ROS1

