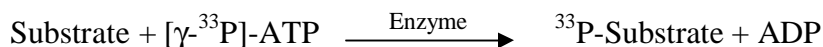


ROS

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

- **CAT#:** ROS
- **Enzyme:** Human ROS
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.9
PP2	5,200
AG1478	13,600
PDGFR I-III	48,000
D-64406	365,000

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

Reference Compound IC50 for ROS

