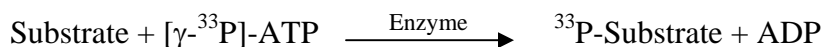


## SRMS

(Tyrosine-protein kinase Srms, C20orf148)

- **CAT#:** SRMS
- **Enzyme:** Human SRMS
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.2 mg/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	15,000
PP2	16,000

- **Scientific Information:** Q9H3Y6 (<http://www.expasy.org/uniprot/Q9H3Y6>)

### Reference Compound IC50 for SRMS

