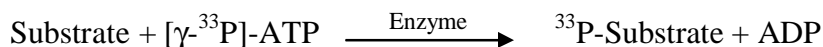


## FES

(Proto-oncogene tyrosine-protein kinase Fes/Fps, C-Fes)

- **CAT#:** FES
- **Enzyme:** Human FES
- **Substrate:** Peptide substrate, poly[Glu:Tyr] (4:1), 0.1 mg/ml  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.0
PP2	29,000
AG1478	605,000
PDGFR I-III	39,000
D-64406	8,300

- **Scientific Information:** P07332 (<http://www.expasy.org/uniprot/P07332>)

### Reference Compound IC50 for FES

