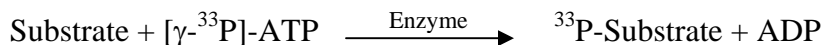


FES/FPS

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

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

- **Reaction:**

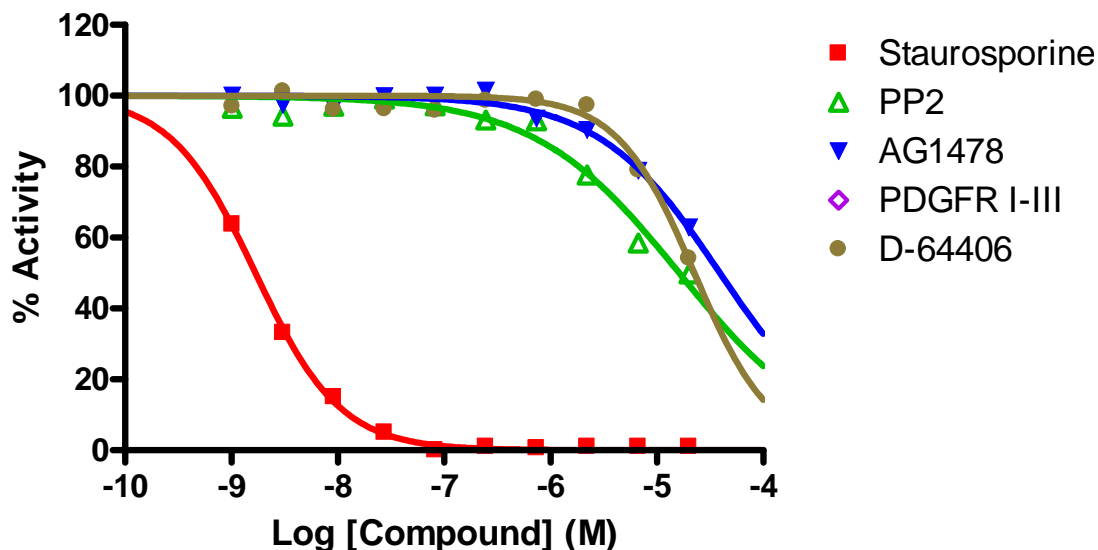


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	1.7
PP2	16,000
AG1478	39,000
D-64406	23,000

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

Reference compound IC50 for FES/FPS



Service performed
using kinases
under license from  **invitrogen**[™]
part of *life* technologies[™]