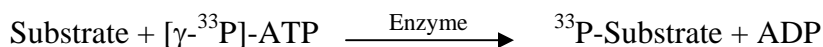


FGR

(Proto-oncogene tyrosine-protein kinase FGR, P55-FGR, C-FGR)

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

- **Reaction:**

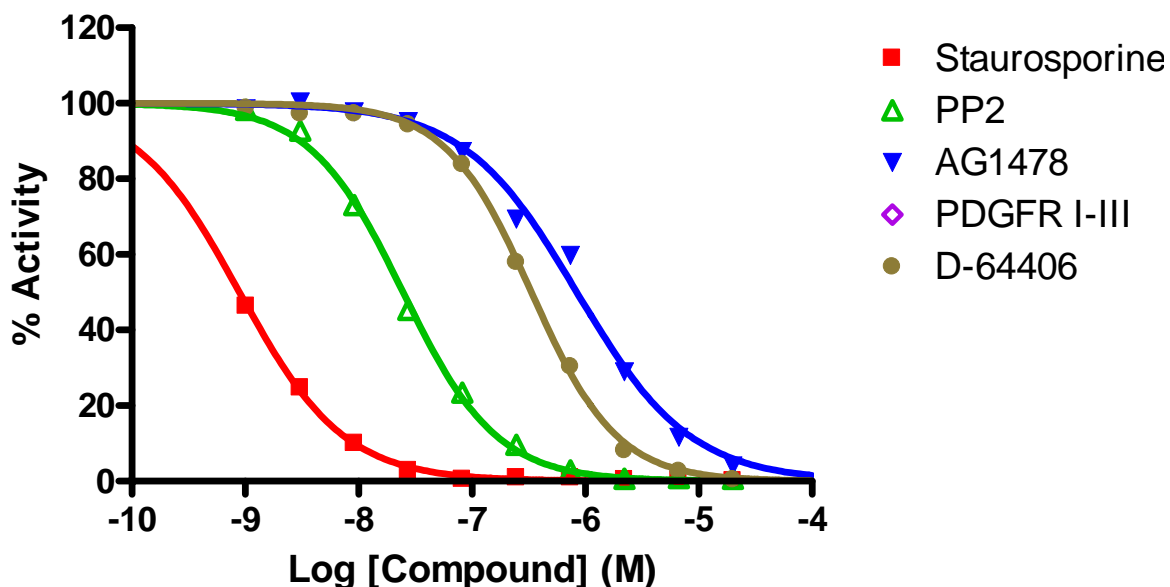


- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.9
PP2	25
AG1478	840
D-64406	330

- **Scientific Information:** P09769 (<http://www.expasy.org/uniprot/P09769>)

Reference compound IC50 for FGR



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