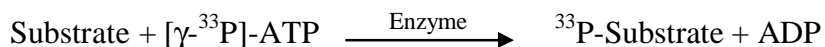


IR

(Insulin receptor [Precursor], Insulin receptor subunit beta, INSR)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.7
PP2	22,400
AG1478	14,200
PDGFR I-III	101,000
D-64406	201,000

- **Scientific Information:** P06213 (<http://www.expasy.org/uniprot/P06213>)

Reference compound IC50 for IR

