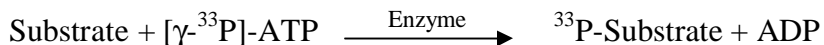


IRR/INSRR

(Insulin receptor-related protein [Precursor], IR-related receptor)

- **CAT#:** IRR
- **Enzyme:** Human IRR
- **Substrate:** MBP (Myelin Basic Protein), 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.6
PP2	5,600
AG1478	7,300
PDGFR I-III	40,000
D-64406	1,800

- **Scientific Information:** P14616 (<http://www.expasy.org/uniprot/P14616>)

Reference Compound IC50 for IRR

