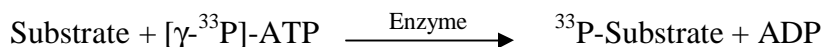


IGF1R

(Insulin-like growth factor 1 receptor [Precursor], IGF-I receptor)

- **CAT#:** IGF1R
- **Enzyme:** Human IGF1R
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 397 |
| PDGFR I-III | 132,000 |
| D-64406 | 106,000 |

- **Scientific Information:** P08069 (<http://www.expasy.org/uniprot/P08069>)

Reference Compound IC50 for IGF1R

