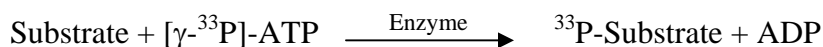


## IGF1R

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

- **CAT#:** IGF1R
- **Enzyme:** Human IGF1R
- **Substrate:** Peptide substrate, [KKKSPGEYVNIEFG], 20  $\mu$ M  
ATP 10  $\mu$ M
- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 9.0       |
| PP2           | >20,000   |

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

### Reference compound IC50 for IGF1R

