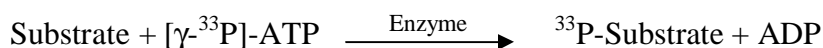


DDR2

(Discoidin domain-containing receptor 2 [Precursor], Discoidin domain receptor 2, Receptor protein-tyrosine kinase TKT, Tyrosine-protein kinase TYRO 10, Neurotrophic tyrosine kinase receptor-related 3, NTRKR3)

- **CAT#:** DDR2
- **Enzyme:** Human DDR2
- **Substrate:** Peptide substrate, [KKSARGDYMTMQIG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	18.3
PP2	629
AG1478	83,000
D-64406	20,300

- **Scientific Information:** Q16832 (<http://www.expasy.org/uniprot/Q16832>)

Reference Compound IC50 for DDR2

