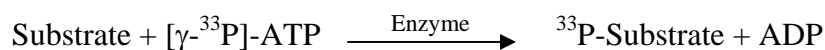


DDR2

(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], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	16
PP2	260
AG1478	9,500
D-64406	30,000

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

Reference compound IC50 for DDR2

