

DYRK2

(Dual specificity tyrosine-phosphorylation-regulated kinase 2)

- **CAT#:** DYRK2
- **Enzyme:** Human DYRK2
- **Substrate:** Peptide substrate, [RRRFRPASPLRGPPK], 20 μ M
ATP 10 μ M
- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	280
Ro 31-8220	5,000
GW 5074	1,500
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q92630 (<http://www.expasy.org/uniprot/Q92630>)

Reference compound IC50 for DYRK2

