

DYRK4

(Dual specificity tyrosine-phosphorylation-regulated kinase 4)

- **CAT#:** DYRK4
- **Enzyme:** Human DYRK4
- **Substrate:** Peptide substrate, [RRRFRPASPLRGPPK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	18,500
Ro 31-8220	>20,000
GW5074	11,500
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q9NR20 (<http://www.expasy.org/uniprot/Q9NR20>)

Reference compound IC50 for DYRK4

