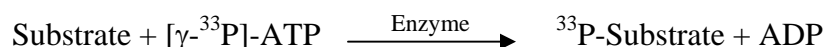


DYRK1/ DYRK1A

(Dual specificity tyrosine-phosphorylation-regulated kinase 1A, Protein kinase minibrain homolog, MNBH, hMNB, HP86, Dual specificity YAK1-related kinase, DYRK, MNB)

- **CAT#:** DYRK1/DYRK1A
- **Enzyme:** Human DYRK1/DYRK1A
- **Substrate:** Peptide substrate, [RRRFRPASPLRGPPK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	3.8
Ro 31-8220	23
GW5074	420
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q13627 (<http://www.expasy.org/uniprot/Q13627>)

Reference compound IC50 for DYRK1/DYRK1A

