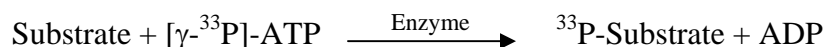


DYRK1B

(Dual specificity tyrosine-phosphorylation-regulated kinase 1B, Mirk protein kinase
Minibrain-related kinase, MIRK)

- **CAT#:** DYRK1B
- **Enzyme:** Human DYRK1B
- **Substrate:** Peptide substrate, [RRRFRPASPLRGPPK], 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.0
Ro 31-8220	68
GW 5074	470
H-89	>20,000
Rottlerin	>20,000

- **Scientific Information:** Q9Y463 (<http://www.expasy.org/uniprot/Q9Y463>)

Reference compound IC50 for DYRK1B

