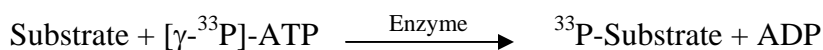


Aurora B

(Serine/threonine-protein kinase 12, Aurora-B, Aurora- and Ip11-like midbody-associated protein 1, AIM-1, Aurora/IPL1-related kinase 2, Aurora-related kinase 2, STK-1, AURKB, AIK2, AIM1, ARK2, STK12)

- **CAT#:** Aurora B
- **Enzyme:** Human Aurora B
- **Substrate:** Peptide substrate, [H-LRRASLG], 100 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4.6
Ro 31-8220	4,000
GW 5074	86
H-89	4,000
Rottlerin	5,300

- **Scientific Information:** Q96GD4 (<http://www.expasy.org/uniprot/Q96GD4>)

Reference compound IC50 for Aurora B

