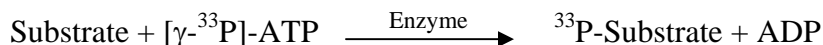


Aurora C

(Serine/threonine-protein kinase 13, AURKC, Aurora/Ipl1/Eg2 protein 2, Aurora/Ipl1-related kinase 3, Aurora-C, AIE2, AIK3, STK13)

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

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	2.6
Ro 31-8220	12,200
GW 5074	6,200
H-89	40,300
Rottlerin	55,700

- **Scientific Information:** Q9UQB9 (<http://www.expasy.org/uniprot/Q9UQB9>)

Reference Compound IC50 for Aurora C

