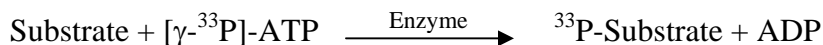


## CHK2 (I157T)

(Serine/threonine-protein kinase Chk2, Cds1, CHEK2)

- **CAT#:** CHK2 (I157T)
- **Enzyme:** Human CHK2 (I157T)
- **Substrate:** Peptide substrate, [KKKVSRSGLYRSPSPENLNRPR], 5  $\mu$ M  
ATP 1  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

| Compounds     | IC50 (nM) |
|---------------|-----------|
| Staurosporine | 3.6       |
| Ro 31-8220    | 957       |
| GW 5074       | 2,200     |
| H-89          | 2,900     |
| Rottlerin     | 6,200     |

- **Scientific Information:** O96017 (<http://www.expasy.org/uniprot/O96017>)

### Reference Compound IC50 for CHK2 (I157T)

