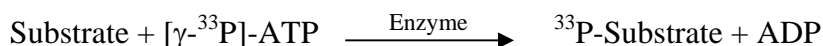


## DNA-PK

(DNA-dependent protein kinase catalytic subunit, DNA-PK catalytic subunit, DNA-PKcs, DNPk1, p460, PRKDC, HYRC, HYRC1)

- **CAT#:** DNA-PK
- **Enzyme:** Human DNA-PK
- **Substrate:** Peptide substrate, [EPPLSQEAFADLWKK], 20  $\mu$ M  
DNA, 10  $\mu$ g/ml  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
LY294002	450
Wortmannin	43
PI3K I	56
PI-103	12

- **Scientific Information:** P78527 (<http://www.expasy.org/uniprot/P78527>)

### Reference Compound IC50 for DNA-PK

