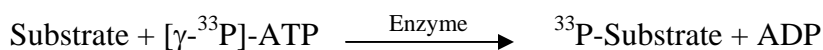


HIPK3

(Homeodomain-interacting protein kinase 3, Homolog of protein kinase YAK1,
Fas-interacting serine/threonine-protein kinase, FIST,
Androgen receptor-interacting nuclear protein kinase, ANPK, DYRK6, FIST3, PKY)

- **CAT#:** HIPK3
- **Enzyme:** Human HIPK3
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	4,600
Ro 31-8220	2,800
GW 5074	3,200
Rottlerin	14,000

- **Scientific Information:** Q9H422 (<http://www.expasy.org/uniprot/Q9H422>)

Reference compound IC50 for HIPK3

