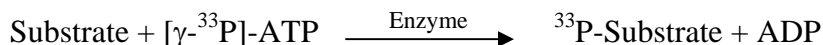


Haspin/GSG2

(Serine/threonine-protein kinase haspin, Haploid germ cell-specific nuclear protein kinase, H-haspin, Germ cell-specific gene 2 protein, HASP)

- **CAT#:** Haspin/GSG2
- **Enzyme:** Human Haspin/GSG2
- **Substrate:** Histone H3, 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	12.3
Ro 31-8220	8,700
GW 5074	25.8
H-89	301
Rottlerin	30,000

- **Scientific Information:** Q8TF76 (<http://www.expasy.org/uniprot/Q8TF76>)

Reference compound IC50 for Haspin/GSG2

