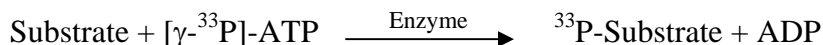


PASK

(PAS domain-containing serine/threonine-protein kinase, PAS-kinase, PASKIN, hPASK)

- **CAT#:** PASK
- **Enzyme:** Human PASK
- **Substrate:** Peptide substrate, [KKLNRTL^SF^AE^PG], 20 μM
ATP 10 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	8.5
Ro 31-8220	178
GW 5074	400
H-89	18,000

- **Scientific Information:** Q96RG2 (<http://www.expasy.org/uniprot/Q96RG2>)

Reference compound IC50 for PASK

