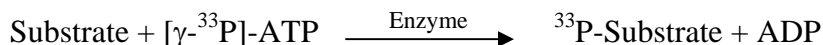


DRAK1/STK17A

(Serine/threonine-protein kinase 17A, DAP kinase-related apoptosis-inducing protein kinase 1, ST17A, STK17A)

- **CAT#:** DRAK1
- **Enzyme:** Human DRAK1
- **Substrate:** Peptide substrate, [KKLNRTL^SF^AE^PG], 20 μM
ATP 10 μM

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	17
Ro 31-8220	5,000
GW 5074	1,400

- **Scientific Information:** Q9UEE5 (<http://www.expasy.org/uniprot/Q9UEE5>)

Reference compound IC50 for DRAK1/STK17A

