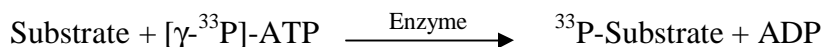


LOK/STK10

(Serine/threonine-protein kinase 10, Lymphocyte-oriented kinase)

- **CAT#:** LOK
- **Enzyme:** Human LOK
- **Substrate:** Peptide substrate, [KKSARGDYMTMQIG], 5 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	65
Ro 31-8220	41
GW 5074	1,100
H-89	10,700
Rottlerin	33,700

- **Scientific Information:** O94804 (<http://www.expasy.org/uniprot/O94804>)

Reference Compound IC50 for LOK

