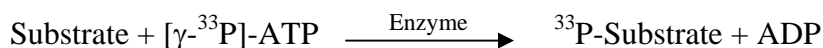


## TAOK3/JIK

(Serine/threonine-protein kinase TAO3, Thousand and one amino acid protein 3, Jun kinase-inhibitory kinase, JNK/SAPK-inhibitory kinase, Dendritic cell-derived protein kinase, Cutaneous T-cell lymphoma tumor antigen HD-CL-09, CTCL tumor antigen HD-CL-09 Kinase from chicken homolog A, hKFC-A, DPK, JIK, KDS, MAP3K18)

- **CAT#:** TAOK3
- **Enzyme:** Human TAOK3
- **Substrate:** MBP (Myelin Basic Protein), 10  $\mu$ M  
ATP 10  $\mu$ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.5
Ro 31-8220	200
GW 5074	1,040
H-89	21,900

- **Scientific Information:** Q9H2K8 (<http://www.expasy.org/uniprot/Q9H2K8>)

Reference compound IC50s for TAOK3/JIK

