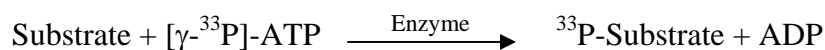


PBK

(Lymphokine-activated killer T-cell-originated protein kinase, T-LAK cell-originated protein kinase, PDZ-binding kinase, Spermatogenesis-related protein kinase, SPK, MAPKK-like protein kinase, Nori-3, Cancer/testis antigen 84, CT84, TOPK)

- **CAT#:** PBK
- **Enzyme:** Human PBK
- **Substrate:** MBP (Myelin Basic Protein), 10 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	141
Ro 31-8220	22,900
GW5074	153,000
Rottlerin	12,700

- **Scientific Information:** Q96KB5 (<http://www.expasy.org/uniprot/Q96KB5>)

