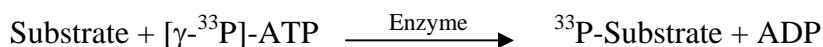


NLK

(Serine/threonine kinase NLK, Nemo-like kinase, Protein LAK1)

- **CAT#:** NLK
- **Enzyme:** Human NLK
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 10 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	86
Ro 31-8220	4,600
GW 5074	490
H-89	23,000
Rottlerin	14,000

- **Scientific Information:** Q9UBE8 (<http://www.expasy.org/uniprot/Q9UBE8>)

Reference compound IC50 for NLK

