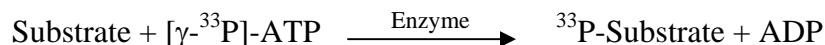


NIK/MAP3K14

(Mitogen-activated protein kinase kinase kinase 14, NF-kappa beta-inducing kinase
Serine/threonine-protein kinase NIK, HsNIK, M3K14)

- **CAT#:** NIK/MAP3K14
- **Enzyme:** Human NIK/MAP3K14
- **Substrate:** MBP (Myelin Basic Protein), 20 μ M
ATP 1 μ M

- **Reaction:**



- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	13.4
Ro 31-8220	27,400
GW 5074	2,100

- **Scientific Information:** Q99558 (<http://www.expasy.org/uniprot/Q99558>)

Reference Compound IC50 for NIK

