

MLK1

(Mitogen-activated protein kinase kinase kinase 9, Mixed lineage kinase 1, PRKE1, MAP3K9, M3K9)

- **CAT#:** MLK1
- **Enzyme:** Human MLK1
- **Substrate:** Casein, 1 mg/ml
ATP 1 μ M
- **Reaction:**
$$\text{Substrate} + [\gamma\text{-}^{33}\text{P}]\text{-ATP} \xrightarrow{\text{Enzyme}} {}^{33}\text{P}\text{-Substrate} + \text{ADP}$$
- **Reference Compound IC50s:**

Compounds	IC50 (nM)
Staurosporine	0.5
Ro 31-8220	146
GW 5074	153
H-89	3,600
Rottlerin	808,000

- **Scientific Information:** P80192 (<http://www.expasy.org/uniprot/P80192>)

Reference Compound IC50 for MLK1



