

---

---

## MLK1(MAP3K9)

Product code: 09-115

Human MLK1, catalytic domain [110-422 amino acids of accession number NP\_149132.2] was expressed as N-terminal GST-fusion protein (62kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	MAP2K1 Cascade Assay *
ATP ( $\mu$ M)	:	1000
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at 1 mM ATP (nM)	:	11

\* MAP2K1/Erk2/Modified Erktide

---

---