
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 (μ M)	: 1000
Metal	: Mg
Reference compound	: Staurosporine
IC50 at 1 mM ATP (nM)	: 11

* MAP2K1/Erk2/Modified Erktide
