
MLK2(MAP3K10)

Product code: 09-116

Human MLK2, catalytic domain and leucine-zipper domain [75-462 amino acids of accession number NP_002437.2] was expressed as N-terminal GST-fusion protein (71kDa) 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)	:	45

* MAP2K1/Erk2/Modified Erktide
