
DLK(MAP3K12)

Product code: 09-111

Human DLK, catalytic domain [1-520 amino acid of accession number NP_006292.3] was expressed as N-terminal GST-fusion protein (86 kDa) using baculovirus expression system.

Assay platform	:	MSA
Substrate	:	MAP2K4/MAP2K7 Cascade Assay *
ATP (μ M)	:	1000
Metal	:	Mg
Reference compound	:	Staurosporine
IC50 at 1 mM ATP (nM)	:	460

* (MAP2K4/MAP2K7)/JNK2/Modified Erktide
