
MNK1(MKKNK1)

Product code: 02-145

Full-length human MNK1 [1-424(end) amino acids and T344D of accession number NP_001129025.1] was expressed as N-terminal GST-fusion protein (74 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: RS peptide
ATP (μ M) Km app / Bin	: 460 / 450
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 21
IC50 at 1 mM ATP (nM)	: n.a.

