
MNK2(MKMK2)

Product code: 02-146

Full-length human MNK2 [1-465(end) amino acids and T379D of accession number NP_951009.1] was expressed as N-terminal GST-fusion protein (79 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: RS peptide
ATP (μ M) Km app / Bin	: 110 / 100
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 7.5
IC50 at 1 mM ATP (nM)	: 44
