
MNK1(MKNK1)

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. GST-MNK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

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.
