

---

## MNK2(MKNK2)

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. GST-MNK2 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : RS peptide

ATP ( $\mu$ 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

---