
MARK4

Product code 02-123

Full-length human MARK4 [1-688(end) amino acids of accession number NP_113605.2] was expressed as N-terminal GST-fusion protein (103 kDa) using baculovirus expression system. GST-MARK4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay

Substrate : CHKtide

ATP (μ M) Km app / Bin : 12 / 10

Metal : Mg

Reference compound : Staurosporine

IC50 at ATP Bin (nM) : 0.12

IC50 at 1 mM ATP (nM) : 5.6
