
NEK4

Product code 05-128

Full-length human NEK4 [1-841(end) amino acids of accession number NP_003148.2] was expressed as N-terminal GST-fusion protein (122 kDa) using baculovirus expression system. GST-NEK4 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : GS peptide

ATP (μ M) Km app / Bin : 51 / 50
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 120
IC50 at 1 mM ATP (nM) : n.a.
