
NEK6

Product code 05-130

Full-length human NEK6 [1-313(end) amino acids of accession number NP_055212.2] was expressed as N-terminal GST-fusion protein (63 kDa) using baculovirus expression system. GST-NEK6 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : CDK7 peptide

ATP (μ M) Km app / Bin : 69 / 75
Metal : Mg
Reference compound : PKR Inhibitor
IC50 at ATP Bin (nM) : 19000
IC50 at 1 mM ATP (nM) : >10000
