
NEK1

Product code 05-123

Human NEK1, catalytic domain [1-505 amino acids of accession number NP_036356.1] was expressed as N-terminal GST-fusion protein (85 kDa) using baculovirus expression system. GST-NEK1 was purified by using glutathione sepharose chromatography.

Assay platform : Mobility Shift Assay
Substrate : CDK7 peptide

ATP (μ M) Km app / Bin : 64 / 75
Metal : Mg
Reference compound : Staurosporine
IC50 at ATP Bin (nM) : 51
IC50 at 1 mM ATP (nM) : 650
