
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.

Assay platform	: MSA
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

