
NEK7

Product code: 05-131

Full-length human NEK7 [1-302(end) amino acids of accession number NP_598001.1] was expressed as N-terminal GST-fusion protein (62 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CDK7 peptide
ATP (μ M) Km app / Bin	: 40 / 50
Metal	: Mg
Reference compound	: PKR Inhibitor
IC50 at ATP Bin (nM)	: 8500
IC50 at 1 mM ATP (nM)	: >10000
