
NEK9

Product code: 05-133

Truncated human NEK9 [1-346, 733-979(end) amino acids of accession number NP_149107.4] was expressed as N-terminal GST-fusion protein (93 kDa) using baculovirus expression system.

Assay platform	: MSA
Substrate	: CDK7 peptide
ATP (μ M) Km app / Bin	: 190 / 200
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 150
IC50 at 1 mM ATP (nM)	: 400
