
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
