
NEK2

Product code: 05-226

Full-length human NEK2 [1-445(end) amino acids of accession number NP_002488.1] was expressed as N-terminal His-tagged protein (55 kDa) using baculovirus expression system. His-tagged NEK2 was purified by using Ni-NTA affinity chromatography. Purified His-NEK2 was digested by recombinant His-TEV protease, and His-tag free NEK2 (ca. 52 kDa) was collected as flow-through fraction from Ni-NTA affinity chromatography.

Assay platform	: MSA
Substrate	: CDK7 peptide
ATP (μ M) Km app / Bin	: 65 / 75
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 3700
IC50 at 1 mM ATP (nM)	: >10000