
NIM1K(MGC42105)

Product code 02-175

Full-length human NIM1K [1-436(end) amino acids of accession number NP_699192.1] was expressed as N-terminal GST-fusion protein (76 kDa) using baculovirus expression system. GST-NIM1K was purified by using glutathione sepharose chromatography and anion exchange chromatography.

Assay platform : Mobility Shift Assay

Substrate : CHKtide

ATP (μ M) K_{mapp} / B_{in} : 21 / 25

Metal : Mg

Reference compound : Staurosporine

IC₅₀ at ATP B_{in} (nM) : 300

IC₅₀ at 1 mM ATP (nM) : n.a.
