
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.

Assay platform	: MSA
Substrate	: CHKtide
ATP (μ M) Km app / Bin	: 21 / 25
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 300
IC50 at 1 mM ATP (nM)	: n.a.

