
NDR1(STK38)

Product code: 01-125

Full-length human NDR1[1-465(end) amino acids of accession number NP_009202.1] was co-expressed as N-terminal GST-fusion protein (81kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP_775739.1] using baculovirus expression system. GST-NDR1 was purified by using glutathione sepharose chromatography.

Assay platform	: MSA
Substrate	: SGKtide
ATP (μ M) Km app / Bin	: 12 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 1.8
IC50 at 1 mM ATP (nM)	: n.a.