

Product Information

NDR1(STK38)

Product Number : **01-125**

Product description

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.

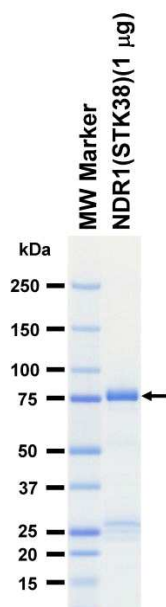
Storage buffer:

50 mM Tris-HCl, 150 mM NaCl, 0.05% Brij35,
1 mM DTT, 10% glycerol, pH7.5

Storage and Handling:

Store at -80C.
Avoid repeating freeze-thaws.

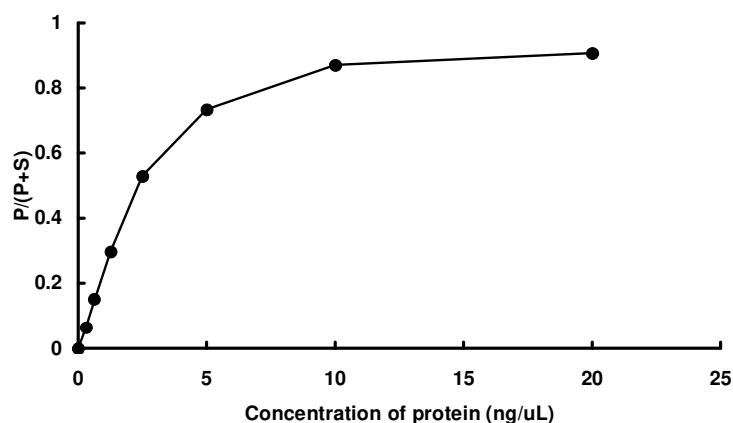
SDS-PAGE



Purity: 72 %

The purity was assessed by SDS-PAGE/CBB staining.

Activity data



The activity was measured by off-chip mobility shift assay(MSA). The enzyme was incubated with fluorescence-labeled substrate and Mg(or Mn)/ATP. The phosphorylated and unphosphorylated substrates were separated and detected by MSA device.

Substrate : SGKtide

ATP : 100 µM