

---

## **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 ( $\mu$ 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.

---