

---

---

## **NDR2(STK38L)**

Product code: 01-126

Full-length human NDR2 [1-464(end) amino acids of accession number NP\_055815.1] was co-expressed as N-terminal GST-fusion protein (81 kDa) with human His-tagged MOBKL1A [1-216(end) amino acids of accession number NP\_775739.1] using baculovirus expression system.

Assay platform	: MSA
Substrate	: SGKtide
ATP ( $\mu$ M) Km app / Bin	: 7.6 / 10
Metal	: Mg
Reference compound	: Staurosporine
IC50 at ATP Bin (nM)	: 2.1
IC50 at 1 mM ATP (nM)	: n.a.

