

## NDR, Unactive

Full-length recombinant protein expressed in Sf9 cells

**Catalog # S49-14G**

Lot # M071-1

### Product Description

Recombinant full-length human NDR was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_007271](#).

### Gene Aliases

NDR1, STK38

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 150mM NaCl, 10mM glutathione, 0.1mM EDTA, 0.25mM DTT, 0.1mM PMSF, 25% glycerol.

### Storage and Stability

Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles.

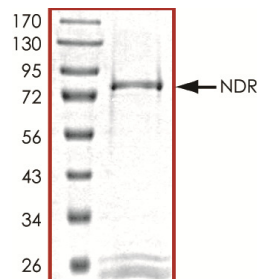
### Scientific Background

NDR or Nuclear Dbf2-related protein is a serine/threonine kinase in the subgroup of the AGC kinase family that is regulated by Mob proteins (1). Fas and TNFα receptor stimulation activates human NDR by promoting phosphorylation at the hydrophobic motif. NDR acts as a novel pro-apoptotic kinase and key member of the RASSF1A/MST1 signaling cascade (2). NDR is crucial for the fidelity of mitotic chromosome alignment in mammalian cells. NDR is incorporated into the HIV-1 particles and cleaved by the HIV-1 protease. NDR-driven centrosome duplication requires Cdk2 activity, and Cdk2-induced centrosome amplification is affected upon reduction of NDR activity.

### References

1. Millward, T. et.al: Molecular cloning and characterization of a conserved nuclear serine(threonine) protein kinase. Proc. Nat. Acad. Sci. 92: 5022-5026, 1995.
2. Vichalkovski A. et al: NDR kinase is activated by RASSF1A/MST1 in response to Fas receptor stimulation and promotes apoptosis. Curr Biol. 2008 Dec 9;18(23):1889-95.

### Purity



The purity of NDR was determined to be **>80%** by densitometry. Approx. MW **81kDa**.

## NDR, Unactive

Full-length recombinant protein expressed in Sf9 cells

Catalog Number **S49-14G**

Specific Lot Number **M071-2**

Purity	>80%
Concentration	0.1 µg/µl
Stability	1yr at -70°C from date of shipment
Storage & Shipping	Store product at -70°C. For optimal storage, aliquot target into smaller quantities after centrifugation and store at recommended temperature. For most favorable performance, avoid repeated handling and multiple freeze/thaw cycles. Product shipped on dry ice.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**