

## STYK1(NOK) Protein

Recombinant human protein expressed in Sf9 cells

**Catalog # S53-35G**

Lot # 1159-2

### Product Description

Recombinant human STYK1(NOK) (58-end) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_018423](#).

### Gene Aliases

NOK; STYK1; DKFZp761P1010; SuRTK106

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM 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.

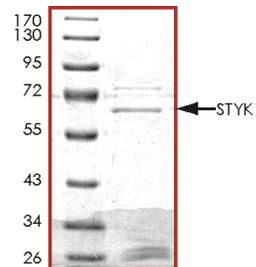
### Scientific Background

STYK1 belongs to serine/threonine/tyrosine kinase 1 family which play important roles in diverse cellular and developmental processes, such as cell proliferation, differentiation, and survival (1). STYK1 concomitantly activate both MAP kinase and PI3K pathways in stably transfected mouse pre-B cells and STYK1 expression is elevated in 92.3% cancerous tissues in comparison with normal tissues (2).

### References

1. Liu, L. et.al: A novel protein tyrosine kinase NOK that shares homology with platelet-derived growth factor/fibroblast growth factor receptors induces tumorigenesis and metastasis in nude mice. *Cancer Res.* 64: 3491-3499, 2004.
2. Amachika, T. et.al: Diagnostic relevance of overexpressed mRNA of novel oncogene with kinase-domain (NOK) in lung cancers. *Lung Cancer* 56: 337-340, 2007.

### Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **67 kDa**.

## STYK1(NOK) Protein

Recombinant human protein expressed in Sf9 cells

Catalog Number	S53-35G
Specific Lot Number	1159-2
Purity	>70%
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.**