

## POMK (SGK196) Protein

Full-length recombinant human protein expressed in Sf9 cells

**Catalog # S55-34G**

Lot # A1423-5

### Product Description

Full-length recombinant human POMK (SGK196) was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. This gene accession number is [NM\\_032237](#).

### Gene Aliases

MDDGA12; SGK196

### Formulation

Recombinant protein stored in 50mM Tris-HCl, pH 7.5, 50mM NaCl, 10mM glutathione, 0.1 mM 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.

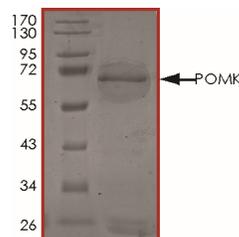
### Scientific Background

POMK or protein-O-mannose kinase encodes a protein that may be involved in the presentation of the laminin-binding O-linked carbohydrate chain of alpha-dystroglycan (α-DG), which forms transmembrane linkages between the extracellular matrix and the exoskeleton. POMK is involved in cell surface presentation of laminin-binding carbohydrates on alpha-dystroglycan (1). Mutations in GTDC2, B3GALNT2, and SGK196 disrupt dystroglycan receptor function and lead to congenital muscular dystrophy (2).

### References

1. Jae, L. T.et.al: Deciphering the glycosylome of dystroglycanopathies using haploid screens for Lassa virus entry. *Science* 340: 479-483, 2013.
2. Yoshida-Moriguchi, T. et.al: SGK196 is a glycosylation-specific O-mannose kinase required for dystroglycan function. *Science* 341: 896-899, 2013.

### Purity



The purity of POMK (SGK196) protein was determined to be **>80%** by densitometry. Approx. MW **68 kDa**.

## POMK (SGK196) Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	S55-34G
Lot #	A1423-5
Purity	>80%
Concentration	0.05 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.**