

## MOS Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # M57-34G**

Lot # N006-4

### Product Description

Recombinant full-length human MOS was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number is [NM\\_005372](#).

### Gene Aliases

MGC119962; MGC119963; MSV

### 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  $-70^{\circ}\text{C}$ . 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.

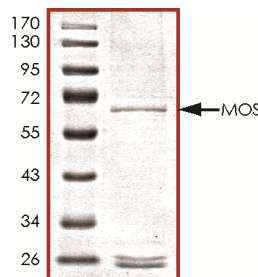
### Scientific Background

MOS or v-mos Moloney murine sarcoma viral oncogene homolog is a serine/threonine kinase that activate the MAP kinase cascade through direct phosphorylation of the MAP kinase activator MEK (1). The early site-specific phosphorylation of CPEB, which was catalyzed by Eg2 is essential for the polyadenylation of c-mos mRNA, subsequent translation of c-mos, and oocyte maturation (2). The interaction with and phosphorylation of MYOD by MOS enhances muscle differentiation. MOS is an essential component of the cytostatic factor that acts almost exclusively in the second meiotic metaphase arrest.

### References

1. Prasad, C. K. et.al: Mos 3-prime UTR regulatory differences underlie species-specific temporal patterns of Mos mRNA cytoplasmic polyadenylation and translational recruitment during oocyte maturation. *Molec. Reprod. Dev.* 75: 1258-1268, 2008.
2. Mendez, R. et.al: Phosphorylation of CPE binding factor by Eg2 regulates translation of c-mos mRNA. *Nature* 404: 302-307, 2000.

### Purity



The purity was determined to be **>70%** by densitometry. Approx. MW **66 kDa**.

## MOS Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	M57-34G
Specific Lot Number	N006-4
Purity	>70%
Concentration	0.1 µg/µl
Stability	1yr At $-70^{\circ}\text{C}$ from date of shipment
Storage & Shipping	Store product at $-70^{\circ}\text{C}$ . 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.**