

## MASTL Protein

Full-length recombinant human protein expressed in Sf9 cells

**Catalog # M72-30G**

Lot # Q2405-4

### Product Description

Full-length recombinant human MASTL was expressed by baculovirus in Sf9 insect cells using an N-terminal GST tag. The gene accession number for MASTL is [NM\\_001172303](#).

### Gene Aliases

GREATWALL; GW; GWL; hGWL; RP11-85G18.2; THC2

### Formulation

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

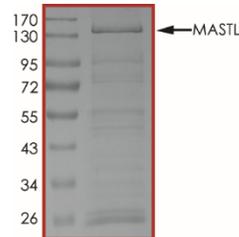
### Scientific Background

Microtubule associated serine/threonine kinase-like (MASTL) belongs to the MAST family, AGC Ser/Thr protein kinase group. MASTL promotes mitotic progression and cell cycle reentry after DNA damage. Loss of MASTL leads to defects in chromosome condensation, separation, and other aspects of mitotic progression, which is primarily mediated by inhibiting PP2A/B55δ. Overexpression of MASTL is correlated with cancer progression, poor patient survival, and tumor recurrence in various cancers.

### References

1. Wang L, et al. Mastl kinase, a promising therapeutic target, promotes cancer recurrence. *Oncotarget*. 5:11479-89, 2014.
2. Zhao Y, et al. Roles of Greatwall kinase in the regulation of cdc25 phosphatase. *Mol Biol Cell*. 19:1317-27, 2008.

### Purity



The purity of MASTL was determined to be **~70%** by densitometry, MASTL approx. MW **150kDa**.

## MASTL Protein

Full-length recombinant human protein expressed in Sf9 cells

Catalog #	M72-30G
Lot #	Q2405-4
Purity	~70%
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.**