

## MGMT Protein

Full length recombinant protein expressed in Sf9 cells

**Catalog # M354-30H**

Lot # F660-2

### Product Description

Recombinant full length human MGMT was expressed by baculovirus in Sf9 insect cells using an N-terminal His tag. The gene accession number is [BC000824](#).

### Gene Aliases

(None)

### Formulation

Recombinant protein stored in 50mM sodium phosphate, pH 7.0, 300mM NaCl, 150mM imidazole, 0.1mM PMSF, 0.25mM DTT, 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.

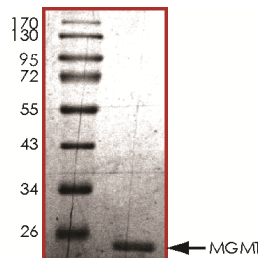
### Scientific Background

MGMT or O-6-methylguanine-DNA methyltransferase is the major mutagenic and carcinogenic lesion in DNA induced by simple alkylating mutagens. The methylation of the MGMT promoter in gliomas is a useful predictor of the responsiveness of the tumors to alkylating agents (1). The clinical trial substantiates the predictive value of O-6-methylguanine-DNA methyltransferase promoter methylation in glioblastoma patients treated with temozolomide (2). Methylation of the MGMT promoter in tumors is associated with longer survival in glioblastoma patients.

### References

1. Esteller, M. et.al: Inactivation of the DNA-repair gene MGMT and the clinical response of gliomas to alkylating agents. New Eng. J. Med. 343: 1350-1354, 2000.
2. Hegi, M. E. et.al: Clinical trial substantiates the predictive value of O-6-methylguanine-DNA methyltransferase promoter methylation in glioblastoma patients treated with temozolomide. Clin. Cancer Res. 10: 1871-1874, 2004.

### Purity



The purity of MGMT was determined to be **>95%** by densitometry. Approx. MW **25kDa**.

## MGMT Protein

Full length recombinant protein expressed in Sf9 cells

Catalog Number	M354-30H
Specific Lot Number	F660-2
Purity	>95%
Concentration	0.1µg/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.**