
MGMT Polyclonal Antibody

(Catalog # A-1010)

Background

O6-methylguanine methyltransferase (MGMT) is a ubiquitous DNA repair protein that removes O6 -alkyl-guanine lesions, primarily O6 -methyl-guanine from damaged DNA. It is a major contributor to cellular protection from the mutagenic, carcinogenic, and cytotoxic effects of DNA alkylation. Also, MGMT is frequently silenced due to abnormal methylation during cancer development. Consequently, the loss of MGMT activity that made the cell more likely to become a cancer cell in the first place also makes tumor cells more sensitive to radiation therapy and certain alkylating drugs.

Description

MGMT Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH7.3.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P16455

Purification

Affinity Purification

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 1-207 of human MGMT (XP_005252739.1).

Storage

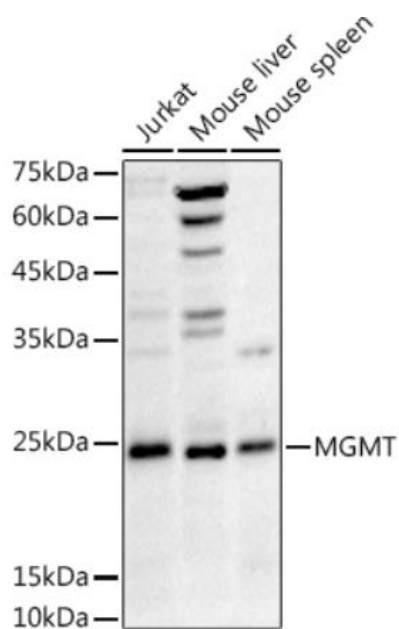
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

6 O methylguanine DNA methyltransferase antibody, Agat antibody, AGT antibody, AI267024 antibody, EC 2.1.1.63 antibody, Methylated DNA protein cysteine methyltransferase antibody, Methylguanine DNA methyltransferase antibody, MGC107020 antibody, MGMT antibody, O 6 methylguanine DNA alkyltransferase antibody, O 6 methylguanine DNA methyltransferase antibody

Application

WB; Recommended dilution: WB: 1:500 - 1:2000



Western blot analysis of extracts of various cell lines, using MGMT Polyclonal Antibody at 1:1000 dilution.
Secondary antibody: HRP Goat Anti-Rabbit IgG (H+L) (AS014) at 1:10000 dilution.
Lysates/proteins: 25ug per lane.
Blocking buffer: 3% nonfat dry milk in TBST.
Exposure time: 180s.