

## MGMT antibody (pAb)

**Catalog Nos:** 39951, 39952

**RRID:** AB\_2793405

**Isotype:** IgG

**Application(s):** IP, WB

**Reactivity:** Human

**Volumes:** 100 µl, 10 µl

**Purification:** Affinity Purified

**Host:** Rabbit

**Molecular Weight:** 28 kDa

**Background:** MGMT (O-6 methylguanine DNA methyltransferase) is an enzyme involved in DNA damage repair. MGMT transfers methyl groups from O-6 methylguanine methylated DNA, and other methylated adducts, to a cysteine residue in itself, thus reversing DNA damage caused by alkylation. This is a suicide reaction as the enzyme is irreversibly inactivated. MGMT helps prevent DNA alkylation events from causing mutations. It also contributes to resistance to chemotherapy in cancer patients being treated with alkylating drugs.

**Immunogen:** This MGMT antibody was raised against a peptide derived from the C-terminal region of human MGMT.

**Buffer:** Purified IgG in 70 mM Tris (pH 8), 105 mM NaCl, 31 mM glycine, 0.07 mM EDTA, 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

### Application Notes:

Applications Validated by Active Motif:

IP: 10 µl per IP

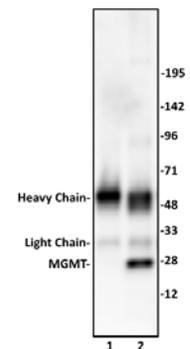
WB\*: 1:500 - 1:1,000 dilution

The addition of 0.05% Tween 20 in the blocking buffer and primary antibody incubation buffer is recommended to aid in detection by Western blot. Individual optimization may be required.

\*Note: many chromatin-bound proteins are not soluble in a low salt nuclear extract and fractionate to the pellet. Therefore, we recommend a High Salt / Sonication Protocol when preparing nuclear extracts for Western blot.

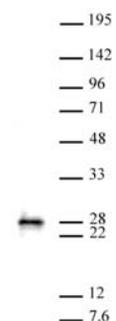
**Storage and Guarantee:** Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.



### MGMT antibody (pAb) tested by Immunoprecipitation.

10 µl of MGMT antibody was used to immunoprecipitate MGMT from 250 µg of MCF-7 nuclear cell extract (lane 2). 10 µl of rabbit IgG was used as a negative control (lane 1). The immunoprecipitated protein was detected by Western blotting using the MGMT antibody at a dilution of 1:500.



### MGMT antibody (pAb) tested by Western blot.

Nuclear extract of MCF-7 cells (20 µg) probed with MGMT antibody (pAb) at a dilution of 1:500.