

---

## GAPDH Monoclonal Antibody [RMC232A]

(Catalog # A68405)

---

### Background

Has both glyceraldehyde-3-phosphate dehydrogenase and nitrosylase activities, thereby playing a role in glycolysis and nuclear functions, respectively. Participates in nuclear events including transcription, RNA transport, DNA replication and apoptosis. Nuclear functions are probably due to the nitrosylase activity that mediates cysteine S-nitrosylation of nuclear target proteins such as SIRT1, HDAC2 and PRKDC. Modulates the organization and assembly of the cytoskeleton. Facilitates the CHP1-dependent microtubule and membrane associations through its ability to stimulate the binding of CHP1 to microtubules (By similarity). Glyceraldehyde-3-phosphate dehydrogenase is a key enzyme in glycolysis that catalyzes the first step of the pathway by converting D-glyceraldehyde 3-phosphate (G3P) into 3-phospho-D-glyceroyl phosphate. Component of the GAIT (gamma interferon-activated inhibitor of translation) complex which mediates interferon-gamma-induced transcript-selective translation inhibition in inflammation processes. Upon interferon-gamma treatment assembles into the GAIT complex which binds to stem loop-containing GAIT elements in the 3'-UTR of diverse inflammatory mRNAs (such as ceruplasmin) and suppresses their translation.

### Description

GAPDH Monoclonal Antibody [RMC232A]. Unconjugated. Raised in: Rabbit.

### Formulation

Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.

### Specificity

Human, Mouse

### Isotype

IgG

### Uniprot ID

P04406

### Purification

Affinity Purified

### Immunogen

A synthesized peptide

### Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

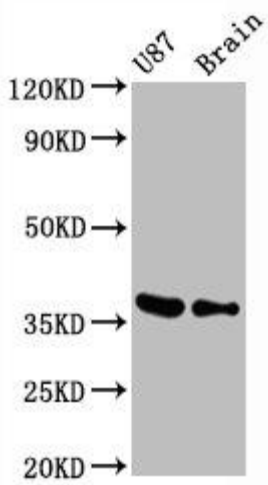
### Alternative Names

Glyceraldehyde-3-phosphate dehydrogenase, GAPDH, Peptidyl-cysteine S-nitrosylase GAPDH, GAPDH, GAPD, CDABP0047, OK/SW-cl.12

### Application

ELISA, WB; Recommended dilution: WB:1:3000-1:10000

---



Western Blot

Positive WB detected in: U87 whole cell lysate, Mouse brain tissue

All lanes: GAPDH Monoclonal Antibody [RMC232A] at 0.31 ug/ml

Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 36 KDa

Observed band size: 36 KDa