

PRODUCT INFORMATION

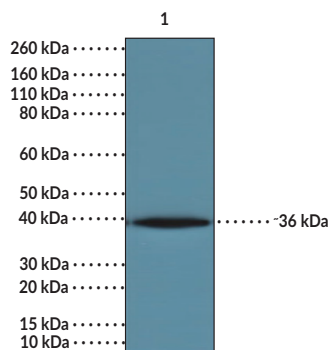
GAPDH (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM114)

Item No. 32378

Overview and Properties

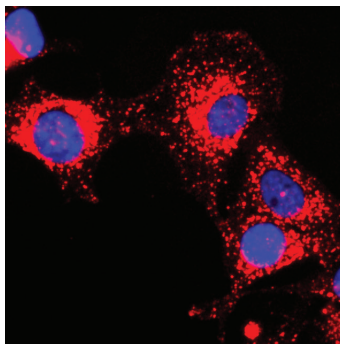
| | |
|----------------------------|--|
| Contents: | This vial contains 50 µl of protein A-affinity purified monoclonal antibody. |
| Synonyms: | G3PD Protein, Human, GAPD Protein, Human, Glyceraldehyde-3-phosphate dehydrogenase, HEL-S-162eP Protein, Peptidyl-cysteine S-nitrosylase |
| Immunogen: | Peptide from the C-terminal region of GAPDH |
| Cross Reactivity: | (+) GAPDH |
| Species Reactivity: | Species independent |
| Form: | Liquid |
| Storage: | -20°C (as supplied) |
| Stability: | ≥1 year |
| Storage Buffer: | PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide |
| Clone: | RM114 |
| Host: | Rabbit |
| Isotype: | IgG |
| Applications: | Chromatin IP (ChIP), Immunocytochemistry (ICC), Immunohistochemistry (IHC), Immunoprecipitation (IP), and Western blot (WB); the recommended starting dilution is 1:200 for ChIP, ICC, IHC, and IP, and 1:1,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically. |

Images



Lane 1: A431 cells

WB of A431 cells using GAPDH (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM114) at a dilution of 1:1,000.



Immunocytochemical staining of HeLa cells using GAPDH (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM114) at a dilution of 1:200 (red). Nuclei have been labeled with DAPI (blue).

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 03/13/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Glyceraldehyde-3-phosphate dehydrogenase (GAPDH) is an enzyme that catalyzes the conversion of glyceraldehyde-3-phosphate (Item No. 17865) to 1,3-bisphosphoglycerate during glycolysis and is involved in numerous additional cellular processes, including intracellular trafficking, receptor-mediated signaling, apoptosis, DNA repair, and the oxidative stress response.^{1,2} It exists as a tetramer and is composed of an N-terminal domain, which contains binding sites for NAD⁺, phosphatidylserine, RNA, and glutathione, and a C-terminal catalytic domain.³ GAPDH is widely expressed and primarily localizes to the cytosol, where it has roles in glycolysis and intracellular trafficking.^{1,4} It also localizes to the nucleus, mediating DNA integrity, gene transcription, and apoptosis, as well as to cellular membranes, where it has roles in membrane fusion and iron transport.⁴ GAPDH expression is increased by insulin, hypoxia-inducible factor-1 (HIF-1), p53, and nitric oxide (NO) and decreased by acetylated histones.^{1,5} Aberrant mRNA and protein levels of GAPDH have been found in tumor biopsies from patients with a variety of cancers, including lung, renal cell, colorectal, or breast cancer.⁶ Cayman's GAPDH (C-Term) Rabbit Monoclonal Antibody - Biotinylated (Clone RM114) can be used for immunocytochemistry (ICC), immunohistochemistry (IHC), immunoprecipitation (IP), chromatin immunoprecipitation (ChIP), and Western blot (WB) applications. The antibody recognizes the C-terminal region of GAPDH at approximately 36 kDa.

References

1. Tristan, C., Shahani, N., Sedlak, T.W., *et al.* The diverse functions of GAPDH: Views from different subcellular compartments. *Cell. Signal.* **23(2)**, 317-323 (2011).
2. Sirover, M.A. On the functional diversity of glyceraldehyde-3-phosphate dehydrogenase: Biochemical mechanisms and regulatory control. *Biochim. Biophys. Acta* **1810(8)**, 741-751 (2011).
3. Sirover, M.A. Structural analysis of glyceraldehyde-3-phosphate dehydrogenase functional diversity. *Int. J. Biochem. Cell Biol.* **57**, 20-26 (2014).
4. Sirover, M.A. Pleiotropic effects of moonlighting glyceraldehyde-3-phosphate dehydrogenase (GAPDH) in cancer progression, invasiveness, and metastases. *Cancer Metastasis Rev.* **37(4)**, 665-676 (2018).
5. Zhang, J.-Y., Zhang, F., Hong, C.-Q., *et al.* Critical protein GAPDH and its regulatory mechanisms in cancer cells. *Cancer Biol. Med.* **12(1)**, 10-22 (2015).
6. Guo, C., Liu, S., and Sun, M.-Z. Novel insight into the role of GAPDH playing in tumor. *Clin. Transl. Oncol.* **15(3)**, 167-172 (2013).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM