

PRODUCT INFORMATION



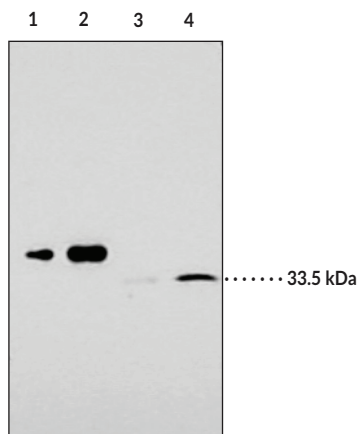
Monoacylglycerol Lipase (FL) Polyclonal Antibody

Item No. 10212

Overview and Properties

Contents:	This vial contains 500 µg protein A-purified polyclonal antibody.
Synonyms:	HU-K5, Lysophospholipase Homology, MAGL, MGL, MGLL
Immunogen:	Purified recombinant human MAGL
Species Reactivity:	(+) Human and rat; other species not tested
Uniprot No.:	Q99685
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥3 years
Storage Buffer:	PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host:	Rabbit
Applications:	Immunocytochemistry (ICC), Immunohistochemistry (IHC), and Western blot (WB); the recommended starting dilution for ICC, IHC, and WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: Human recombinant MAGL (His-tagged) (0.01 µg)
Lane 2: Human recombinant MAGL (His-tagged) (0.02 µg)
Lane 3: Rat adipose tissue homogenate (10 µg)
Lane 4: Rat adipose tissue homogenate (20 µg)

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, 11/03/2023

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

Endocannabinoids, such as arachidonoyl ethanolamide (AEA) and 2-arachidonoyl glycerol (2-AG), function as short-range modulators of cell and synaptic activity. Monoacylglycerol lipase (MAGL) hydrolyzes 2-AG to terminate its biological actions and works consecutively with hormone-sensitive lipase to mobilize fatty acids from the triglyceride stores of adipocytes.^{1,2} MAGL has a molecular weight of ~33 kDa and exhibits a high degree of homology among human, murine, and rat at the amino acid level.¹⁻⁴ The enzyme is expressed in a variety of tissues with most abundant expression in skeletal muscle and adipose tissue, suggesting a role of MAGL in monoglyceride hydrolysis in diverse tissues. Cayman's MAGL (FL) Polyclonal Antibody can be used for Western blot, immunocytochemistry, and immunohistochemistry analysis for MAGL.

References

1. Dinh, T.P., Carpenter, D., Leslie, F.M., *et al.* Brain monoglyceride lipase participating in endocannabinoid inactivation. *Proc.Natl.Acad.Sci. USA* **99(16)**, 10819-10824 (2002).
2. Karlsson, M., Reue, K., Xia, Y.-R., *et al.* Exon-intron organization and chromosomal localization of the mouse monoglyceride lipase gene. *Gene* **272**, 11-18 (2001).
3. Karlsson, M., Contreras, J.A., Hellman, U., *et al.* cDNA cloning, tissue distribution, and identification of the catalytic triad of monoglyceride lipase. Evolutionary relationship to esterases, lysophospholipases, and haloperoxidases. *J.Biol.Chem.* **272**, 27218-27223 (1997).
4. Dinh, T.P., Freund, T.F., and Piomelli, D. A role for monoglyceride lipase in 2-arachidonoyl glycerol inactivation. *Chem.Phys.Lipids* **121**, 149-158 (2002).

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