

PRODUCT INFORMATION



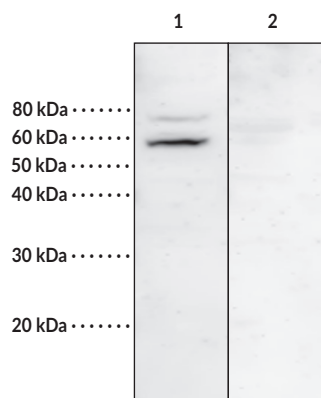
MBOAT1 Polyclonal Antibody

Item No. 14698

Overview and Properties

Contents: This vial contains 500 μ l of peptide affinity-purified polyclonal antibody.
Synonym: Membrane-bound O-Acyltransferase Domain-containing Protein 1
Immunogen: Human MBOAT1, C-terminal region
Species Reactivity: (+) Human; other species not tested
Uniprot No.: Q6ZNC8
Form: Liquid
Storage: -20°C (as supplied)
Stability: \geq 3 years
Storage Buffer: PBS, pH 7.2, with 50% glycerol and 0.02% sodium azide
Host: Rabbit
Applications: Western blot (WB); the recommended starting dilution for WB is 1:200. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Image



Lane 1: RT4 Cell Lysates (50 μ g)
Lane 2: RT4 Cell Lysates (50 μ g) +
10 μ g/ml immunizing peptide

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, 08/11/2025

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

Membrane bound O-acyl transferases (MBOATs) are a group of conserved, multiple transmembrane spanning enzymes involved in many biological functions including lipid biosynthesis, embryogenesis, nutrient sensing, and membrane lipid remodeling. This group of proteins is associated with pathologies such as diabetes, obesity, atherosclerosis, and Alzheimer's disease.¹ MBOAT1 functions as an acyltransferase which mediates the conversion of lysophosphatidylserine (1-acyl-2-hydroxy-*sn*-glycero-3-phospho-L-serine or LPS) into phosphatidylserine (1,2-diacyl-*sn*-glycero-3-phospho-L-serine or PS) (LPSAT activity), and prefers oleoyl-CoA as the acyl donor. Lysophospholipid acyltransferases catalyze the reacylation step of the phospholipid remodeling pathway also known as the Lands cycle.² The predicted size of MBOAT1 is 56.6 kDa, and Cayman's MBOAT1 Polyclonal Antibody detects a band at approximately 60 kDa by Western blot.

References

1. Chang, C.C.Y., Sun, J., and Chang, T.Y. Membrane-bound O-acyltransferases (MBOATs). *Front. Biol.* **6(3)**, 177-182 (2011).
2. Gijón, M.A., Riekhof, W.R., Zarini, S., *et al.* Lysophospholipid acyltransferases and arachidonate recycling in human neutrophils. *J. Biol. Chem.* **283(44)**, 30235-30245 (2008).

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