

PRODUCT INFORMATION



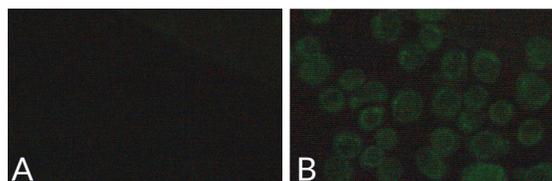
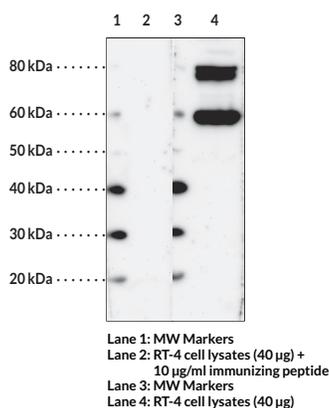
HHATL Polyclonal Antibody

Item No. 15648

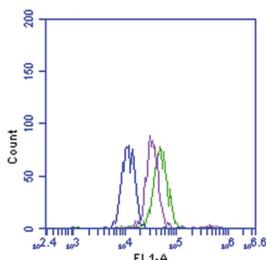
Overview and Properties

Contents: This vial contains 500 μ l of peptide affinity-purified polyclonal antibody.
Synonyms: C3orf3, Glycerol Uptake Transporter Homolog, GUP1, Hedgehog Acyltransferase-like Protein, KIAA1173, MBOAT3, MSTP002, OACT3, Protein-Cysteine N-Palmitoyltransferase HHAT-like Protein
Immunogen: Synthetic peptide from the C-terminal region of human HHAT
Species Reactivity: (+) Human HHATL; other species not tested
Uniprot No.: Q9HCP6
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: Flow cytometry (FC), immunofluorescence (IF), and Western blot (WB); the recommended starting dilution for FC and IF is 1:100 and 1:200 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



Immunofluorescent staining of RT-4 cells. RT-4 cells were fixed with cytospin, washed with 95% ethanol, and blocked with 1% fetal bovine serum. Cells were probed with the indicated antibodies, washed between steps, and images were captured using a Leica DMIL inverted fluorescence microscope (40X objective). Panel A: Control secondary antibody alone Panel B: HHATL Polyclonal Antibody (1:50)



Blue: Goat Anti-Rabbit IgG FITC (Item No. 10006588) (1:200)
Purple: HHATL Polyclonal Antibody (1:100)
Green: HHATL Polyclonal Antibody (1:50)
RT-4 cells were fixed with 4% formaldehyde and permeabilized with Cayman permeabilization buffer, followed by blocking with 1% fetal bovine serum. Cells were probed with indicated antibodies, washed between steps, and fluorescence was detected with a BD Accuri C6 flow cytometer.

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, 11/14/2023

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

Hedgehog acyltransferase-like (HHATL) is the mammalian homolog of yeast Gup1 that is encoded by the *HHATL* gene in humans.¹ HHATL is a membrane-bound O-acyltransferase (MBOAT) superfamily member and contains multiple transmembrane domains common to this family but lacks acyltransferase function due to a single amino acid substitution of histidine for leucine at position 447.^{1,2} It is localized to the endoplasmic reticulum (ER) and is expressed in the heart, skeletal muscle, and brain.^{1,3,4} HHATL colocalizes with sonic hedgehog (Shh) and is a negative regulator of Shh N-terminal palmitoylation, a post-translational modification that is critical for Shh signaling in neural development and embryogenesis.³ *Hhntl*^{-/-} neonatal mice fail to develop proper suckling ability, leading to malnutrition and death by postnatal day 14.⁵ *HHATL* expression is decreased in six nasopharyngeal carcinoma cell lines, as well as in tissue isolated from patients with nasopharyngeal carcinoma or skin squamous cell carcinoma.^{6,7} Cayman's HHATL Polyclonal Antibody can be used for flow cytometry (FC), immunofluorescence (IF), and Western blot applications. The antibody recognizes HHATL at ~60 kDa from human samples.

References

1. Lucas, C., Ferreira, C., Cazzanelli, G., *et al.* Yeast Gup1(2) proteins are homologues of the hedgehog morphogens acyltransferases HHAT(L): Facts and implications. *J. Dev. Biol.* **4(4)**, 33 (2016).
2. Chang, C.C.Y., Sun, J., and Chang, T.Y. Membrane-bound O-acyltransferases (MBOATs). *Front. Biol.* **6(3)**, 177-182 (2011).
3. Abe, Y., Kita, Y., and Niikura, T. Mammalian Gup1, a homolog of *Saccharomyces cerevisiae* glycerol uptake/transporter 1, acts as a negative regulator for N-terminal palmitoylation of Sonic hedgehog. *FEBS J.* **275(2)**, 318-331 (2008).
4. Bleve, G., Zacheo, G., Cappello, M.S., *et al.* Subcellular localization and functional expression of the glycerol uptake protein 1 (GUP1) of *Saccharomyces cerevisiae* tagged with green fluorescent protein. *Biochem. J.* **390(Pt. 1)**, 145-155 (2005).
5. Van, B., Nishi, M., Komazaki, S., *et al.* Mitsugumin 56 (hedgehog acyltransferase-like) is a sarcoplasmic reticulum-resident protein essential for postnatal muscle maturation. *FEBS Lett.* **589(10)**, 1095-1104 (2015).
6. Zhang, S.-Q., Peng, H., Song, L.-Y., *et al.* Detection of KIAA1173 gene expression in nasopharyngeal carcinoma tissues and cell lines on tissue microarray. *Ai Zheng* **24(11)**, 1322-1326 (2005).
7. Zhang, S.-Q., Tian, X., Luo, Y.-W., *et al.* Expression, clinical and pathological significance of KIAA1173 gene in skin squamous cell carcinoma. *Zhonghua Yi Xue Za Zhi* **90(18)**, 1243-1246 (2010).

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