

LAPTM5 Polyclonal Antibody

(Catalog # A64814)

Background

May have a special functional role during embryogenesis and in adult hematopoietic cells.

Description

LAPTM5 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

Q13571

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Lysosomal-associated transmembrane protein 5 protein (205-262AA)

Storage

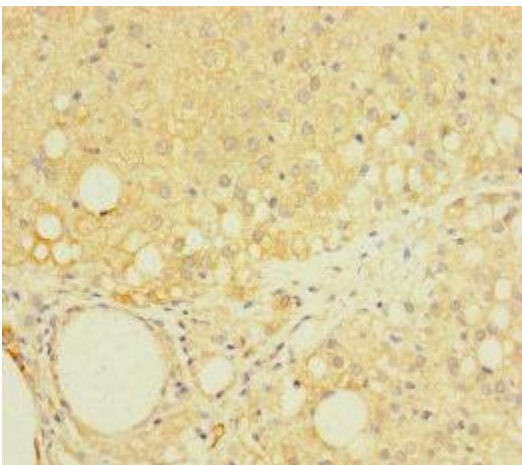
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

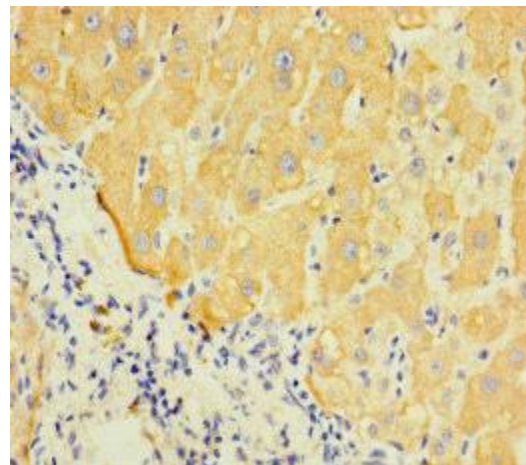
Lysosomal-associated transmembrane protein 5, Lysosomal-associated multitransmembrane protein 5, Retinoic acid-inducible E3 protein, LAPTM5, KIAA0085

Application

ELISA, IHC; Recommended dilution: IHC:1:20-1:200



Immunohistochemistry of paraffin-embedded human adrenal gland tissue using LAPTM5 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human liver cancer using LAPTM5 Polyclonal Antibody at dilution of 1:100