
ANGPTL3 Polyclonal Antibody

(Catalog #A72615)

Background

This gene encodes a member of a family of secreted proteins that function in angiogenesis. The encoded protein, which is expressed predominantly in the liver, is further processed into an N-terminal coiled-coil domain-containing chain and a C-terminal fibrinogen chain. The N-terminal chain is important for lipid metabolism, while the C-terminal chain may be involved in angiogenesis. Mutations in this gene cause familial hypobetalipoproteinemia type 2.

Description

ANGPTL3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Buffer: PBS with 0.02% sodium azide, 50% glycerol, pH 7.3.

Specificity

Mouse

Isotype

IgG

Uniprot ID

Q9Y5C1

Purification

Affinity Purified

Immunogen

Recombinant fusion protein containing a sequence corresponding to amino acids 231-460 of human ANGPTL3 (NP_055310.1).

Storage

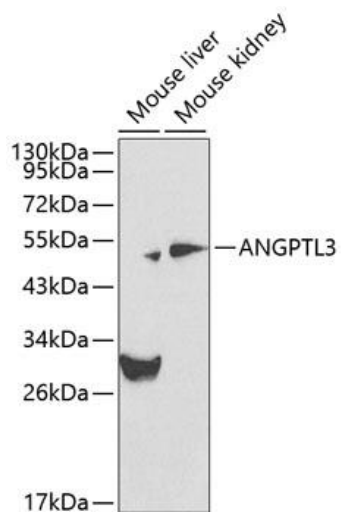
Shipped at 4°C. Upon receipt, store at -20°C. Avoid freeze / thaw cycles

Alternative Names

ANGPTL3; ANG-5; ANGPT5; ANL3; FHBL2; angiopoietin like 3

Application

ELISA, WB; Recommended dilution: WB 1:500 – 1:2000



Western blot analysis of extracts of various cell lines, using ANGPTL3 Polyclonal Antibody at 1:1000 dilution.

Secondary Antibody: HRP Goat Anti-Rabbit IgG (H+L) at 1:10000 dilution.

Lysates/proteins: 25ug per lane.

Blocking buffer: 3% nonfat dry milk in TBST.