

SAMD7 Polyclonal Antibody

(Catalog # A65970)

Description

SAMD7 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

Q7Z3H4

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Sterile alpha motif domain-containing protein 7 protein (1-300AA)

Storage

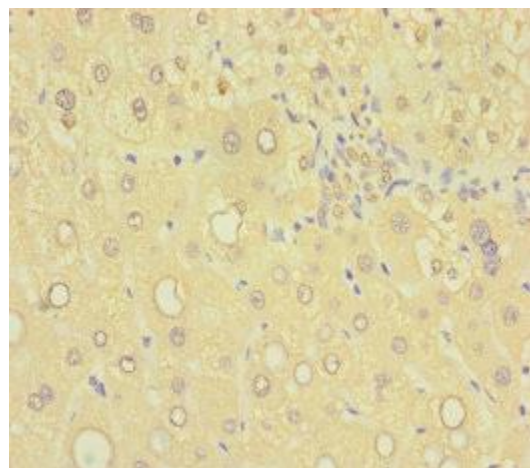
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

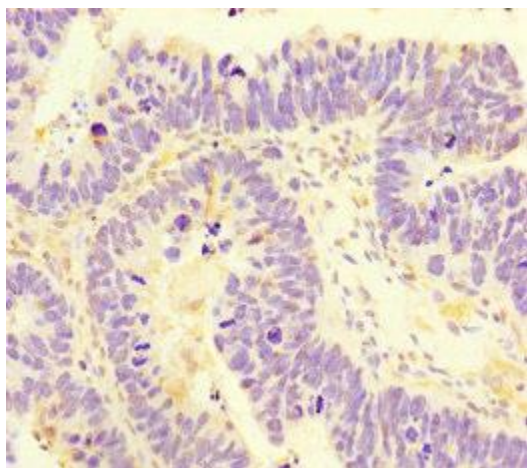
Sterile alpha motif domain-containing protein 7, SAM domain-containing protein 7, SAMD7

Application

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



Immunohistochemistry of paraffin-embedded human liver tissue using SAMD7 Polyclonal Antibody at dilution of 1:100



Immunohistochemistry of paraffin-embedded human ovarian cancer using SAMD7 Polyclonal Antibody at dilution of 1:100