

TM4SF5 Polyclonal Antibody

(Catalog # A63568)

Background

integral component of plasma membrane

Description

TM4SF5 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

O14894

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Transmembrane 4 L6 family member 5 protein (112-157AA)

Storage

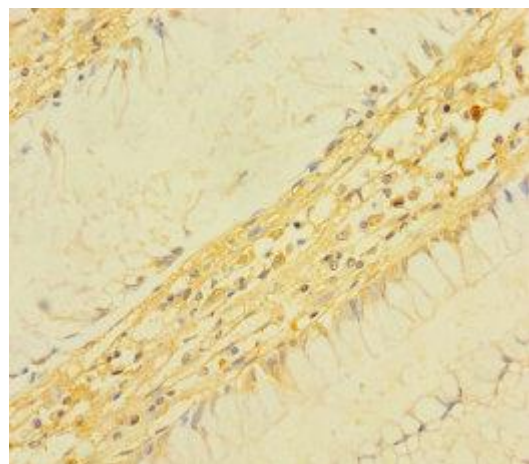
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

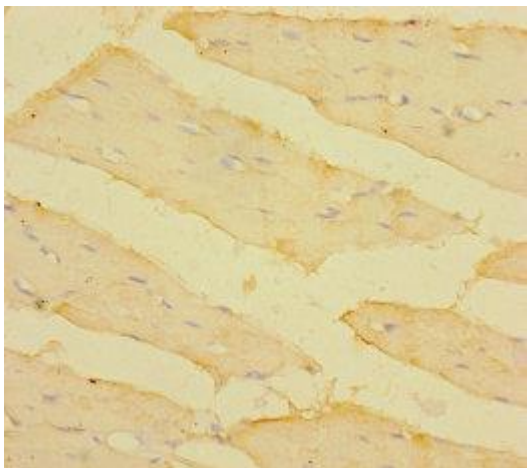
Transmembrane 4 L6 family member 5, Tetraspan transmembrane protein L6H, TM4SF5

Application

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



Immunohistochemistry of paraffin-embedded human colon cancer at dilution of 1:100



Immunohistochemistry of paraffin-embedded human skeletal muscle tissue at dilution of 1:100