

TMEM143 Polyclonal Antibody

(Catalog # A63588)

Background

mitochondrion

Description

TMEM143 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

Q96AN5

Purification

Affinity purified

Immunogen

Recombinant Human Transmembrane protein 143 protein (8-278AA)

Storage

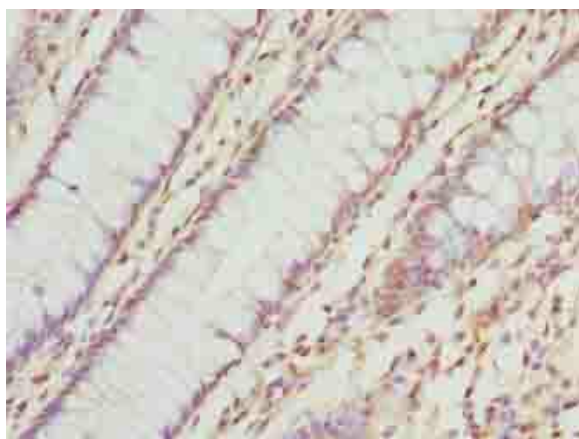
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

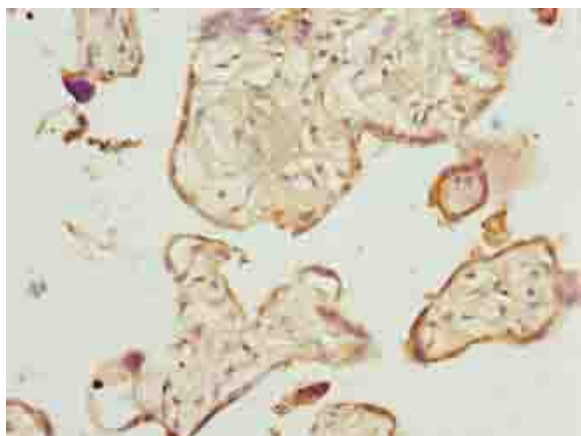
Transmembrane protein 143, TMEM143, UNQ5922/PRO19813

Application

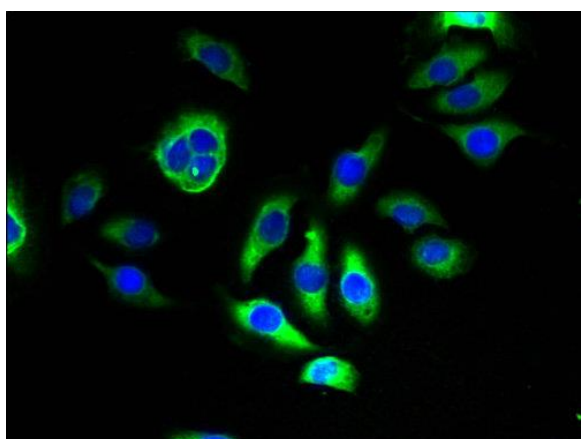
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



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



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



Immunofluorescent analysis of A549 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)