
EMC6 Polyclonal Antibody

(Catalog #A58926)

Background

ER membrane protein complex, integral component of endoplasmic reticulum membrane, integral component of membrane, integral component of omegasome membrane, autophagosome assembly, protein folding in endoplasmic reticulum

Description

EMC6 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

Q9BV81

Purification

>95%, Protein G purified

Immunogen

Recombinant Human ER membrane protein complex subunit 6 protein (2-47AA)

Storage

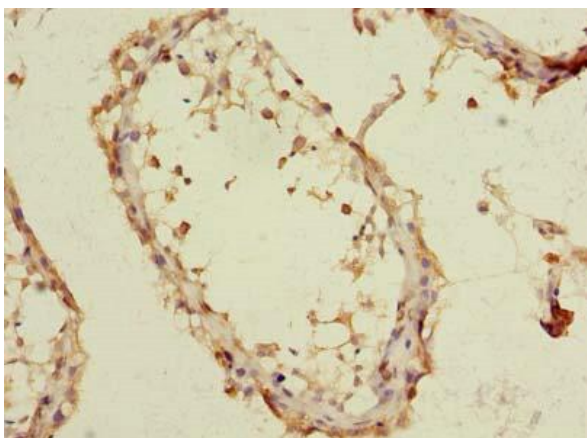
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

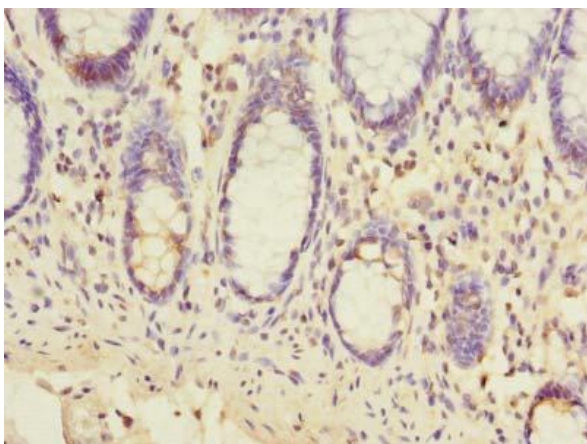
ER membrane protein complex subunit 6, EMC6, TMEM93, Transmembrane protein 93

Application

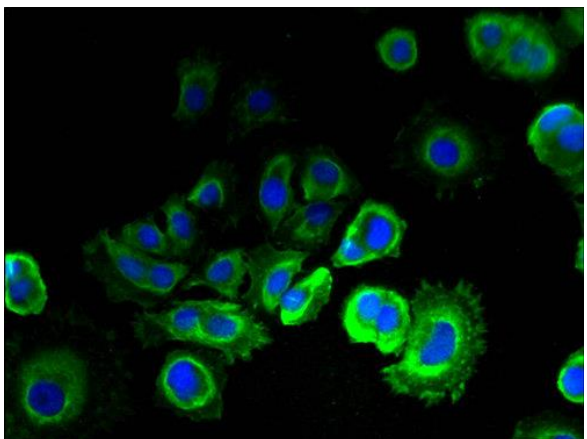
ELISA, IHC, IF; Recommended dilution: IHC:1:20-1:200, IF:1:50-1:200



Immunohistochemistry of paraffin-embedded human testis at dilution of 1:100



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



Immunofluorescent analysis of MCF-7 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)