

COL8A1 Polyclonal Antibody

(Catalog # A62310)

Background

Macromolecular component of the subendothelium. Major component of the Descemet's membrane (basement membrane) of corneal endothelial cells. Also component of the endothelia of blood vessels. Necessary for migration and proliferation of vascular smooth muscle cells and thus, has a potential role in the maintenance of vessel wall integrity and structure, in particular in atherogenesis.

Description

COL8A1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

P27658

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Collagen alpha-1(VIII) chain protein (572-744AA)

Storage

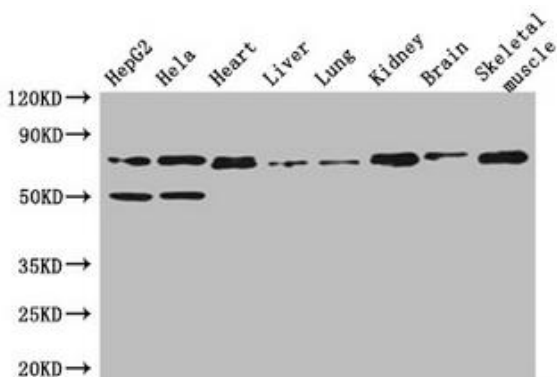
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Collagen alpha-1(VIII) chain, Endothelial collagen, COL8A1, C3orf7

Application

ELISA, WB, IF; Recommended dilution: WB:1:500-1:5000, IF:1:50-1:200



Western Blot

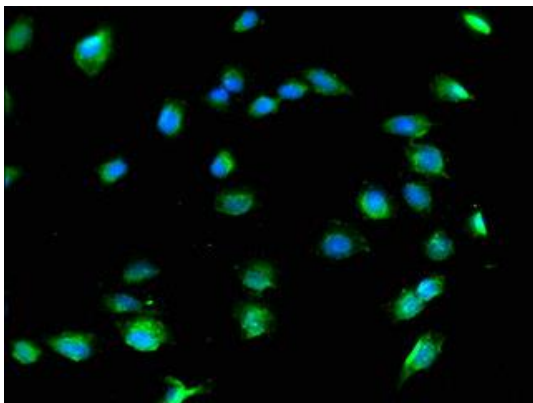
Positive WB detected in: HepG2 whole cell lysate, HeLa whole cell lysate, mouse heart tissue, mouse liver tissue, mouse lung tissue, mouse kidney tissue, mouse brain tissue, mouse skeletal muscle tissue

All lanes: COL8A1 Polyclonal Antibody at 3ug/ml
Secondary

Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 74 kDa

Observed band size: 74, 50 kDa



Immunofluorescent analysis of U251 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)