

CTGF Polyclonal Antibody

(Catalog # A52384)

Background

Major connective tissue mitogen secreted by vascular endothelial cells. Promotes proliferation and differentiation of chondrocytes. Mediates heparin- and divalent cation-dependent cell adhesion in many cell types including fibroblasts, myofibroblasts, endothelial and epithelial cells. Enhances fibroblast growth factor-induced DNA synthesis.

Description

CTGF 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

P29279

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Connective tissue growth factor protein (253-349AA)

Storage

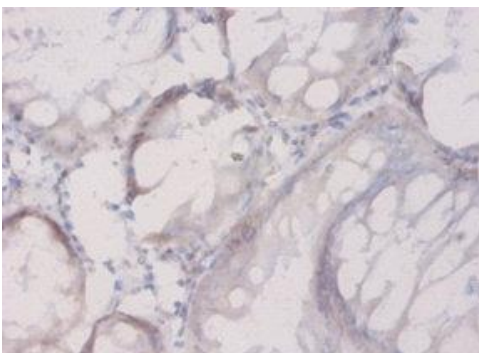
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

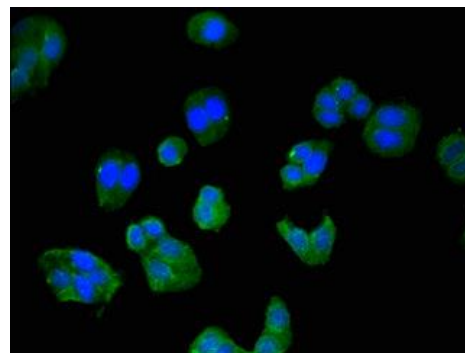
CCN family member 2, Hypertrophic chondrocyte-specific protein 24, Insulin-like growth factor-binding protein 8, CTGF, CCN2, HCS24, IGFBP8

Application

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



Immunohistochemistry of paraffin-embedded human colorectal carcinoma at dilution of 1:50



Immunofluorescent analysis of HepG2 cells at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L)