

CEACAM1 Polyclonal Antibody

(Catalog # A58526)

Background

extracellular exosome, integral component of membrane, integral component of plasma membrane, membrane, plasma membrane, protein homodimerization activity, angiogenesis, blood coagulation, cell migration, homophilic cell adhesion via plasma membrane adhesion molecules

Description

CEACAM1 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

P13688

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Carcinoembryonic antigen-related cell adhesion molecule 1 protein (35-428AA)

Storage

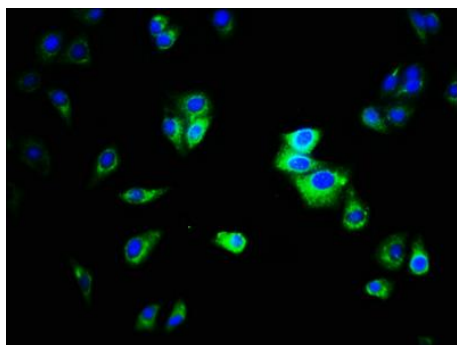
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

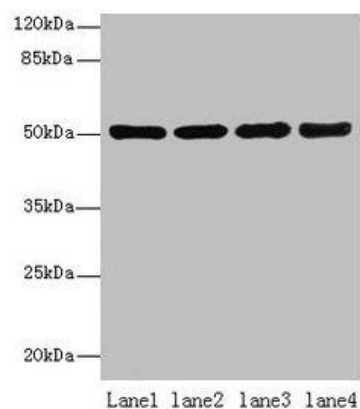
Carcinoembryonic antigen-related cell adhesion molecule 1, CEACAM1, BGP, BGP1, Biliary glycoprotein 1, BGP-1, CD_antigen: CD66a

Application

ELISA, WB, IF; Recommended dilution: WB:1:1000-1:5000, IF:1:200-1:500



Immunofluorescence staining of HepG2 cells with CEACAM1 Antibody at 1:333, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG(H+L).



Western blot

All lanes: CEACAM1 antibody at 10ug/ml

Lane 1: HT29 whole cell lysate

Lane 2: HCT116 whole cell lysate

Lane 3: COLO205 whole cell lysate

Lane 4: HepG2 whole cell lysate

Secondary

Goat polyclonal to rabbit IgG at 1/10000 dilution

Predicted band size: 58, 46, 36, 39, 51, 47, 28, 44, 52, 40 kDa

Observed band size: 51 kDa