

CSMD3 Polyclonal Antibody

(Catalog # A69991)

Background

Involved in dendrite development.

Description

CSMD3 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human

Isotype

IgG

Uniprot ID

Q7Z407

Purification

Protein G purified

Immunogen

Recombinant Human CUB and sushi domain-containing protein 3 protein (233-380AA)

Storage

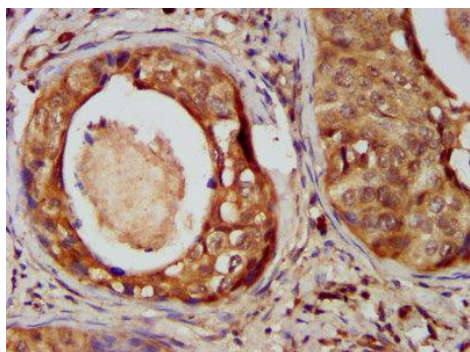
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

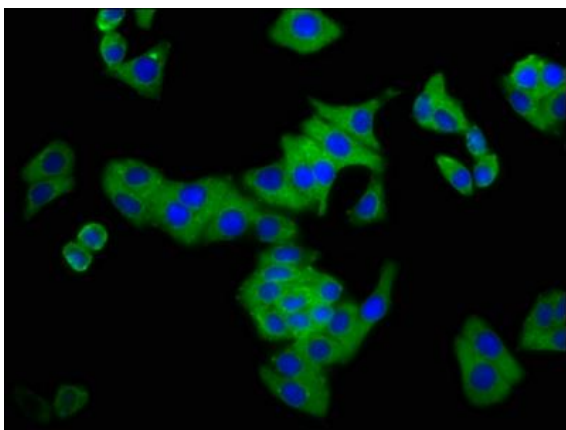
CUB and sushi domain-containing protein 3, CUB and sushi multiple domains protein 3, CSMD3, KIAA1894

Application

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



IHC image of CSMD3 Polyclonal Antibody diluted at 1:400 and staining in paraffin-embedded human cervical cancer performed on a Leica Bond™ system. After dewaxing and hydration, antigen retrieval was mediated by high pressure in a citrate buffer (pH 6.0). Section was blocked with 10% normal goat serum 30min at RT. Then primary antibody (1% BSA) was incubated at 4°C overnight. The primary is detected by a biotinylated secondary antibody and visualized using an HRP conjugated SP system.



Immunofluorescence staining of HepG2 cells with CSMD3 Polyclonal Antibody at 1:133, 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).