

Claudin-18.2 Recombinant Monoclonal Antibody [9C2]

(Catalog # A74527)

Description

Claudin-18.2 Recombinant Monoclonal Antibody [9C2]. Unconjugated. Raised in: HEK293F Cell.

Formulation

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

Specificity

Human

Isotype

hlgG1

Uniprot ID

P56856-2

Purification

Affinity Chromatography

Immunogen

Recombinant Human Claudin-18.2 protein

Storage

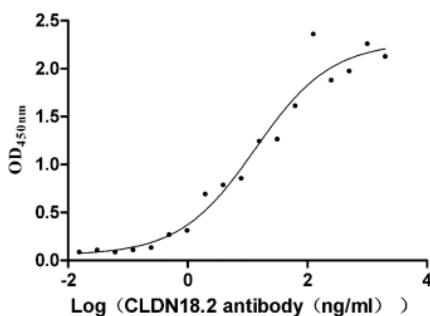
Upon receipt, store at -20°C or -80°C. Avoid repeated freeze.

Alternative Names

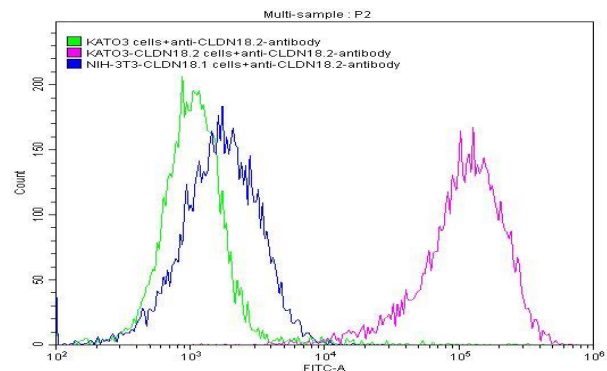
Claudin-18, CLDN18

Application

ELISA, FC; Recommended dilution: FC:1:50-1:200



The Binding Activity of Human CLDN18. 2 with Anti-CLDN18. 2 recombinant antibody. Activity: Measured by its binding ability in a functional ELISA. Immobilized Human CLDN18. 2) at 5 ug/mL can bind Anti-CLDN18. 2 recombinant antibody, the EC₅₀ is 6. 554-27. 87 ng/mL



Untransfected KATO3 cells (green line), transfected Human CLDN18. 2 KATO3 stable cells (red line) and transfected Human CLDN18. 1 NIH-3T3 stable cells (blue line) were stained with anti-CLDN18. 2 antibody (2µg/1*10⁶ cells), washed and then followed by FITC-conjugated anti-Human IgG Fc antibody and analyzed with flow cytometry.