

## CD9 Monoclonal Antibody [RMC969A]

(Catalog # A68409)

### Background

Involved in platelet activation and aggregation. Regulates paranodal junction formation. Involved in cell adhesion, cell motility and tumor metastasis. Required for sperm-egg fusion.

### Description

CD9 Monoclonal Antibody [RMC969A]. Unconjugated. Raised in: Rabbit.

### Formulation

Rabbit IgG in phosphate buffered saline, pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol. .

### Specificity

Human

### Isotype

IgG

### Uniprot ID

P21926

### Purification

Affinity Purified

### Immunogen

A synthesized peptide

### Storage

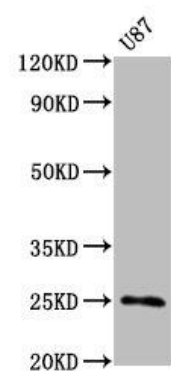
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

CD9 antigen, 5H9 antigen, Cell growth-inhibiting gene 2 protein, Leukocyte antigen MIC3, Motility-related protein, MRP-1, Tetraspanin-29, Tspan-29, p24, CD9, CD9, MIC3, TSPAN29, GIG2

### Application

ELISA, WB, IHC, IF, FC; Recommended dilution: WB:1:500-1:5000, IHC:1:50-1:500, IF:1:30-1:200



#### Western Blot

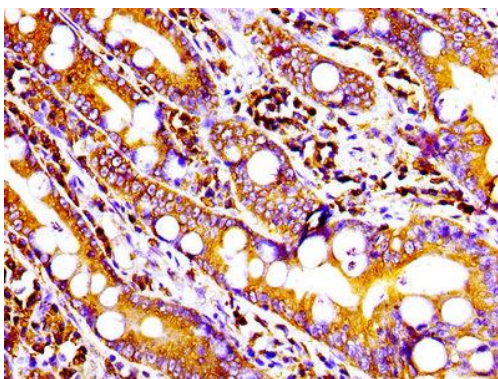
Positive WB detected in: U87 whole cell lysate

All lanes: CD9 Monoclonal Antibody [RMC969A] at 0.55 ug/ml  
Secondary

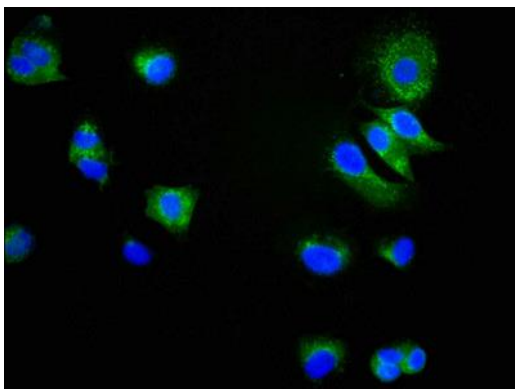
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 25 KDa

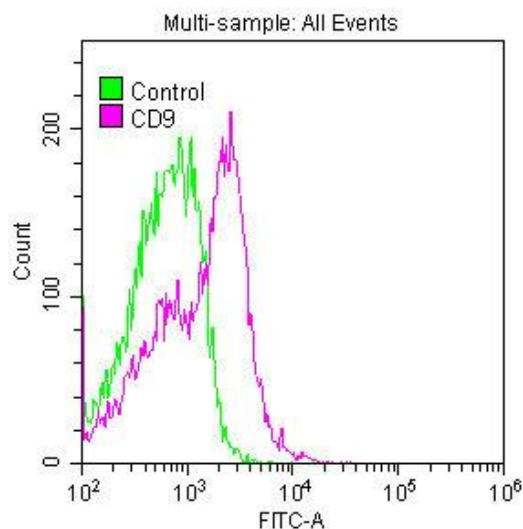
Observed band size: 25 KDa



IHC image of CD9 Monoclonal Antibody [RMC969A] diluted at 1:100 and staining in paraffin-embedded human small intestine tissue 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 MCF-7 cells with CD9 Monoclonal Antibody [RMC969A] at 1:34, 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) conditions. Acquisition of >10,000 events was performed.



Overlay histogram showing Jurkat cells stained with CD9 Monoclonal Antibody [RMC969A] (red line) at 1:50. The cells were fixed with 70% ethylalcohol (18h) and then permeabilized with 0.3% Triton X-100 for 2 min. The cells were then incubated in 1x PBS /10% normal goat serum to block non-specific protein-protein interactions followed by primary antibody for 1 h at 4°C. The secondary antibody used was FITC goat anti-rabbit IgG (H+L) at 1/200 dilution for 1 h at 4°C. Control antibody (green line) was used under the same