

CD81 Monoclonal Antibody, clone RMC960A

(Catalog A68436)

Background

May play an important role in the regulation of lymphoma cell growth. Interacts with a 16-kDa Leu-13 protein to form a complex possibly involved in signal transduction. May act as the viral receptor for HCV.

Description

CD81 Monoclonal Antibody, clone RMC960A. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse, Rat

Isotype

IgG

Uniprot ID

P60033

Purification

Affinity Purified

Immunogen

A synthesized peptide

Storage

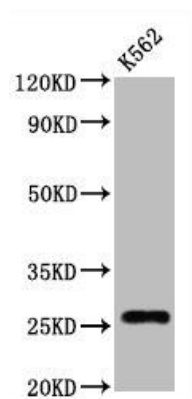
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

CD81 antigen, 26 kDa cell surface protein TAPA-1, Target of the antiproliferative antibody 1, Tetraspanin-28, Tspan-28, CD81

Application

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



Western Blot

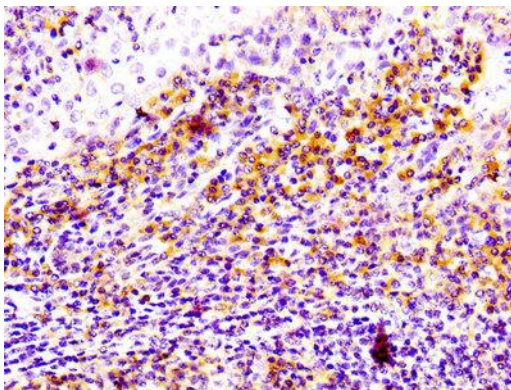
Positive WB detected in: K562 whole cell lysate

All lanes: CD81 Monoclonal Antibody [RMC960A] at 1.25 ug/ml
Secondary

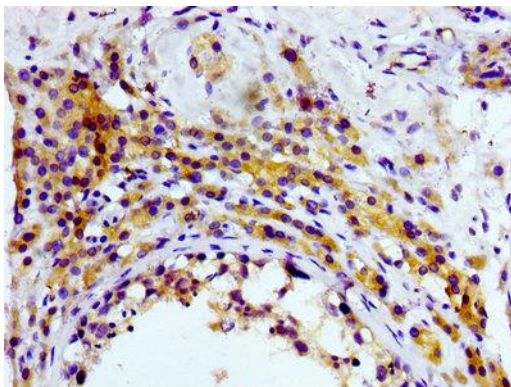
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 26 kDa

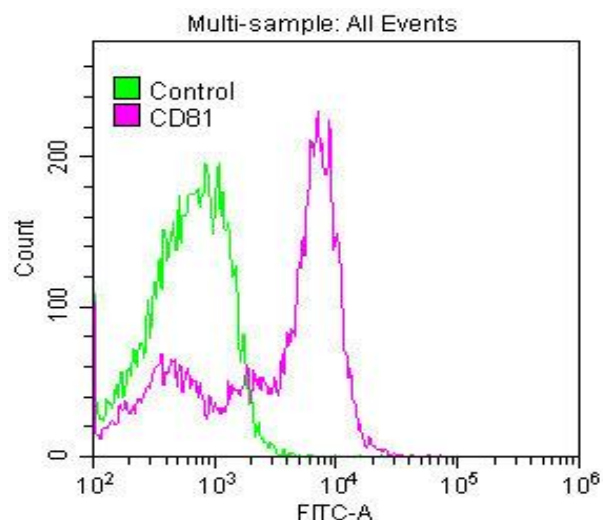
Observed band size: 26 kDa



IHC image of CD81 Monoclonal Antibody [RMC960A] diluted at 1:100 and staining in paraffin-embedded human tonsil 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.



IHC image of CD81 Monoclonal Antibody [RMC960A] diluted at 1:100 and staining in paraffin-embedded human testis 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.



Overlay histogram showing Jurkat cells stained with CD81 Monoclonal Antibody [RMC960A] (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 conditions. Acquisition of >10,000 events was performed.