CD46 monoclonal antibody (AT2G9)

Membrane Cofactor Protein (CD46) is a type 1 membrane protein that plays an important inhibitory role in the complement system. CD46 is implicated in the development and/or progression of selected cancer types.

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

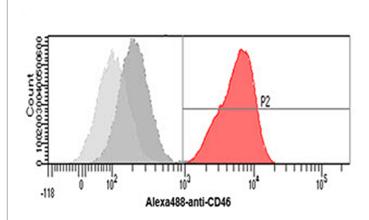
Order Online »

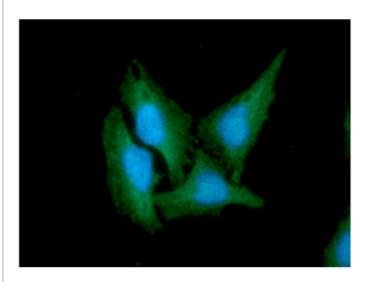
ENZ-ABS722-0100

100µl

Manuals, SDS & CofA

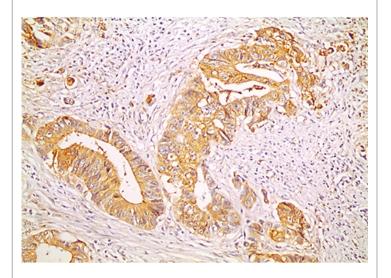
View Online »

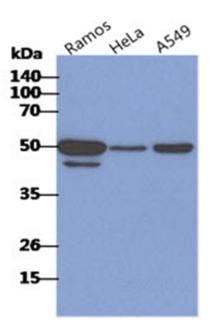




Flow cytometry analysis of CD46 in PBMC cells. The cell was stained with ENZ-ABS722 at 2-5 μ g for 1×10⁶ cells (red). A Goat anti mouse IgG (Alexa fluor 488) was used as the secondary antibody. Mouse monoclonal IgG was used as the isotype control (dark gray), cells without incubation with primary and secondary antibody was used as the negative control (light gray).

ICC/IF analysis of CD46 in HeLa cells. The cell was stained with ENZ-ABS722 (1:100). The secondary antibody (green) was used Alexa Fluor 488. DAPI was stained the cell nucleus (blue).





Paraffin embedded sections of human colorectal cancer tissue were incubated with anti-human CD46 (1:100) for 2 hours at room temperature. Antigen retrieval was performed in 0.1M sodium citrate buffer and detected using Diaminobenzidine (DAB).

The cell lysates of Ramos, HeLa and A549 (40µg) were resolved by SDS-PAGE, transferred to PVDF membrane and probed with anti-human CD46 antibody (1:1000). Proteins were visualized using a goat anti-mouse secondary antibody conjugated to HRP and an ECL detection system.

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Membrane cofactor protein isoform 14, TLX, Trophoblast

leukocyte common antigen

Application ELISA, Flow Cytometry, ICC, IF, IHC, WB

Clone AT2G9

Formulation Liquid. In PBS, pH 7.4, containing 0.02% sodium azide

and 10% glycerol.

Host Mouse

Immunogen Recombinant human CD46 (aa 35-313) purified from E.

coli.

Isotype IgG2ak

Purity Detail Protein A affinity purified

Recommendation Dilutions/ConditionsWestern Blot (1:1,000).Suggested dilutions/conditions may

not be available for all applications. Optimal conditions must be determined individually for each application.

Source Purified from ascites.

Species Reactivity Human

UniProt ID P15529

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com