CD2 recombinant monoclonal antibody (118)

The CD2 antibody is useful for immunological studies, especially in T cell activation, adhesion mechanisms, and the analysis of immune cell interactions.

CD2 is a surface glycoprotein found primarily on T cells and natural killer (NK) cells, playing a critical role in cell adhesion and signal transduction during immune responses.

This antibody is covered by our Worry-Free Guarantee.

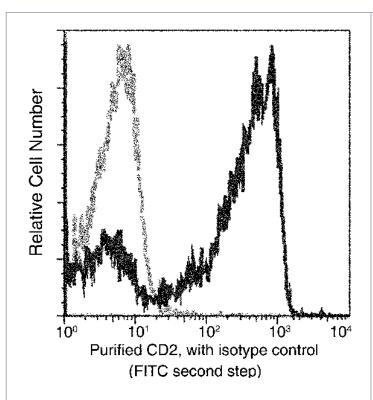
Ordering Information

Order Online »

ENZ-ABS1029-0100 100μl

Manuals, SDS & CofA

View Online »



Flow cytometric analysis of Human CD2 expression on human whole blood lymphocytes. Cells were stained with purified anti-Human CD2, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of viable lymphocytes.

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 LFA-2

Application Flow Cytometry

Clone 118

Formulation Liquid. 0.2 µm filtered solution in PBS. Preservative free.

Carrier-free.

Host Rabbit

Immunogen Recombinant human CD2.

lgG lgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:25-1:100)Suggested

dilutions/conditions may not be available for all applications. Optimal conditions must be determined

individually for each application.

Species Reactivity Human

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

