

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](#).

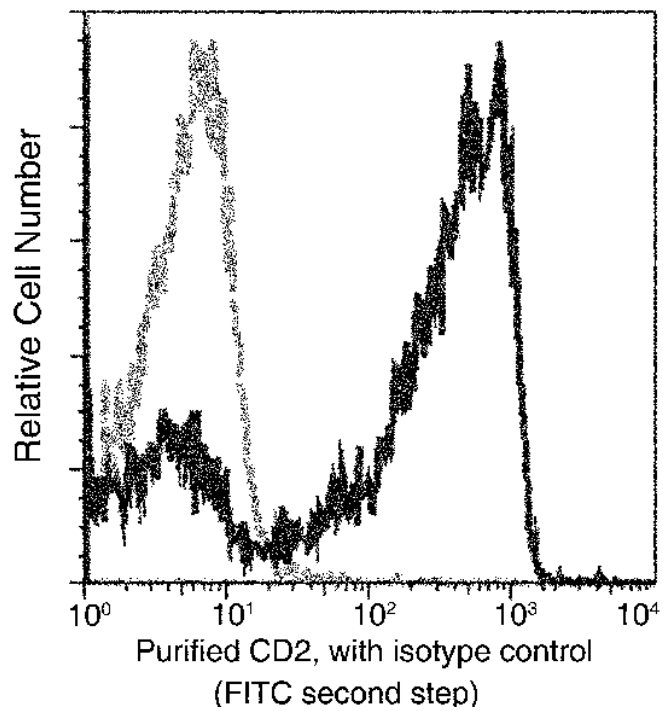
Ordering Information

[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.

Isotype IgG

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](#).



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)