

CD155 recombinant monoclonal antibody (305)

CD155, also known as poliovirus receptor (PVR), is a cell adhesion molecule involved in immune cell interactions and tumor immunity. It is a ligand for immune checkpoint receptors like DNAM-1 and TIGIT, making it a significant focus in cancer immunotherapy research.

This antibody is covered by our [Worry-Free Guarantee](#).

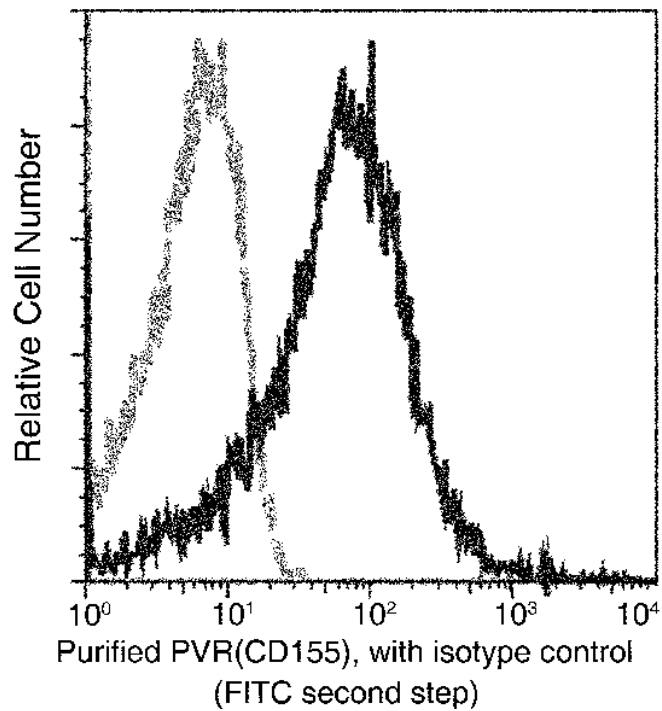
Ordering Information

[Order Online »](#)

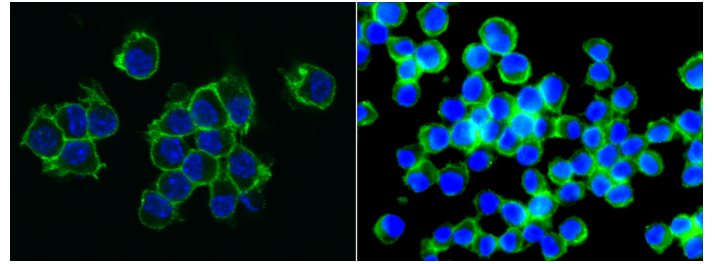
ENZ-ABS1039-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse CD155 expression on BABL/c splenocytes. Cells were stained with purified anti-Mouse CD155, then a FITC-conjugated second step antibody. The fluorescence histograms were derived from gated events with the forward and side light-scatter characteristics of intact cells.



Immunofluorescence staining of Mouse CD155 in raw264.7 cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with Rabbit anti-Mouse CD155 monoclonal antibody (1:100) at 4°C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated (left panel, captured by laser confocal scanning microscope; right panel, captured by fluorescence microscope) Goat Anti-rabbit IgG secondary antibody, counterstained with DAPI (blue). Positive staining was localized to plasma membrane.

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 Poliovirus Receptor, PVR, Nectin-like protein 5, NECL-5

Application Flow Cytometry, ICC, IF

Clone 305

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

Host Rabbit

Immunogen Recombinant mouse CD155 / PVR.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions Immunocytochemistry / Immunofluorescence (1:50-1:1'000) 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 Mouse

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800 942 9436
info-usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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