

# CD63 monoclonal antibody (12)

This monoclonal antibody detects CD63, a tetraspanin protein commonly used as a marker for exosomes and lysosomal trafficking.

CD63 is a member of the tetraspanin family and is involved in intracellular trafficking, cell adhesion, and immune signaling. It is widely used as an exosome marker and has been linked to tumor progression and immune responses.

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

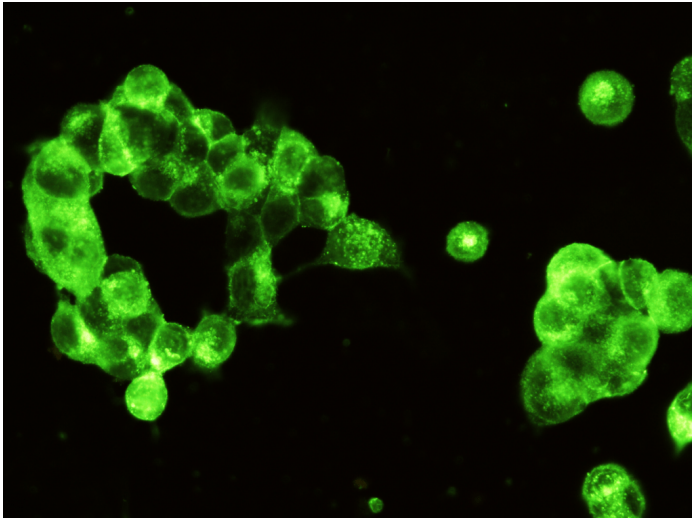
## Ordering Information

[Order Online »](#)

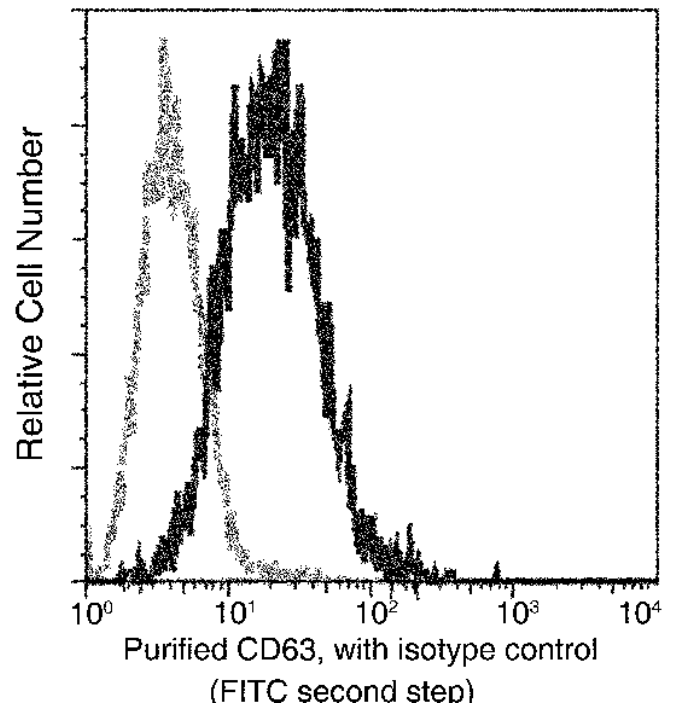
ENZ-ABS1067-0100	100µl
------------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of CD63 in MCF7 cells. Cells were fixed with 4% PFA, permeabilized with 0.3% Triton X-100 in PBS, blocked with 10% serum, and incubated with CD63 monoclonal antibody (1:60) at 4°C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-mouse IgG secondary antibody(green).



Flow cytometric analysis of Human CD63 expression on MCF-7 cells. Cells were stained with purified anti-Human CD63, then a FITC-conjugated second step antibody. The histogram were derived from gated events with the forward and side light-scatter characteristics of intact cells.

# 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	Tetraspanin-30, Tspan-30
Application	Flow Cytometry, ICC, IF
Clone	12
Formulation	Liquid. 0.2 µm filtered solution in PBS. Carrier-free. Preservative-free.
Host	Mouse
Immunogen	Recombinant human CD63 / Tspan-30 / Tetraspanin-30 protein.
Isotype	IgG
Purity Detail	Protein A purified.
Recommendation Dilutions/Conditions	Immunocytochemistry/Immuofluorescence (1:20-1:100) 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 <a href="#">Worry-Free Guarantee</a> .



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)