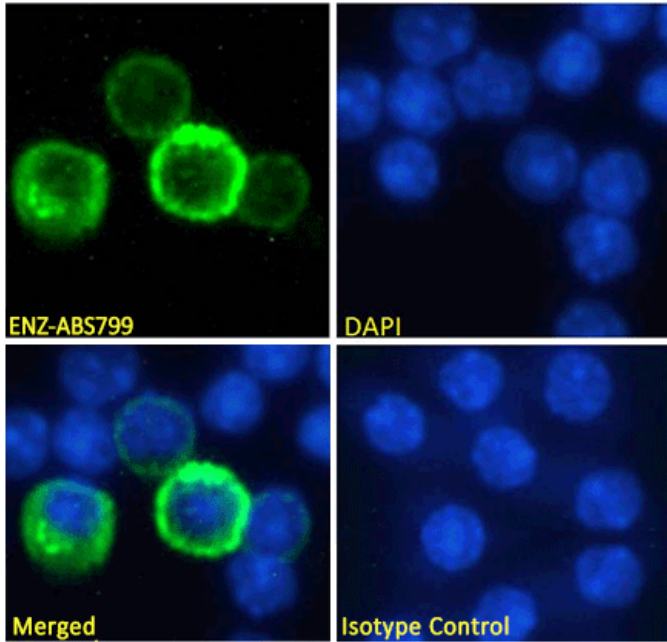


CD27 Recombinant monoclonal antibody (LG.3A10)

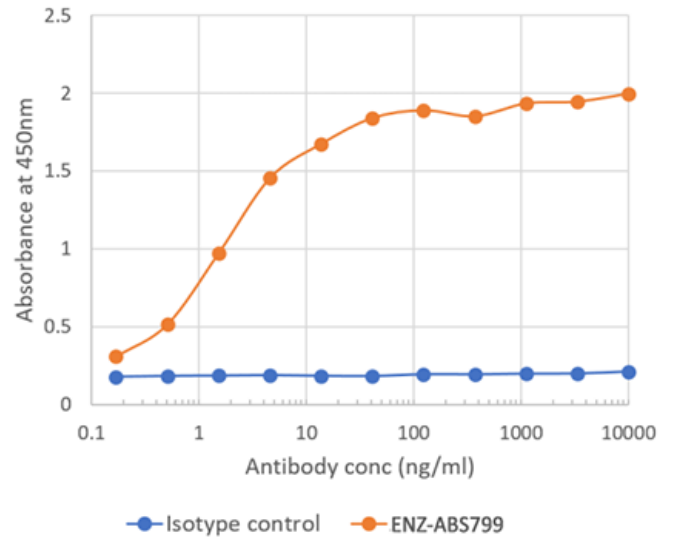
This antibody is covered by our [Worry-Free Guarantee](#).

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of fixed mouse splenocytes with ENZ-ABS799. Immunofluorescence analysis of paraformaldehyde fixed mouse (*Mus musculus*) splenocytes. Panels show from left-right, top-bottom ENZ-ABS799, DAPI, merged channels and an isotype control. The isotype control was stained with an anti-Fluorescein antibody (Ab102-23.0) followed by Alexa Fluor® 488 secondary antibody.



Binding curve of CD27 antibody to mouse CD27-Fc fusion protein.

Handling & Storage

Use/Stability	Stable at 4°C for up to 3 months. For longer storage, aliquot and store at -20°C.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TNFRSF7, CD27L receptor, T-cell activation antigen CD27
Application	ELISA, IF
Clone	LG.3A10
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	This antibody was produced by injecting Armenian hamsters intraperitoneally with a mixture of two murine CD27 tag-expressing Armenian hamster fibroblast clones (AR0524 and AR0530) in Freund's adjuvant.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Source	This chimeric rabbit antibody was made using the variable domain sequences of the original Hamster IgG format, for improved compatibility with existing reagents, assays and techniques.
Species Reactivity	Mouse
UniProt ID	P41272
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .

Last modified: May 29, 2024



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)