

CD27 Recombinant monoclonal antibody (RM27-3E5)

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

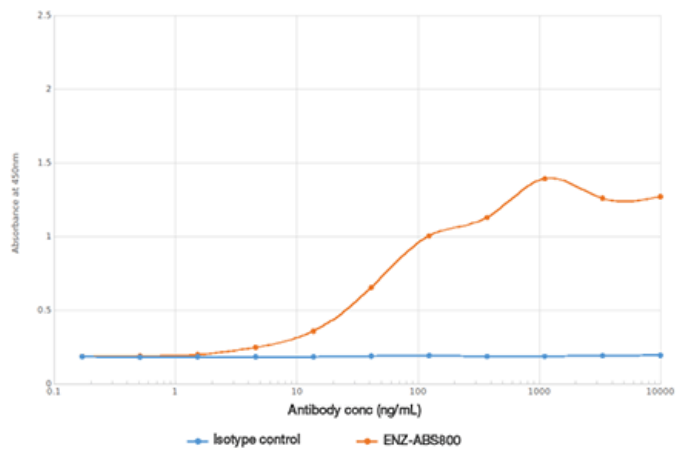
Ordering Information

[Order Online »](#)

ENZ-ABS800-0200	200µg
-----------------	-------

Manuals, SDS & CofA

[View Online »](#)



ELISA of CD27 antibody on 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
Clone	RM27-3E5
Formulation	Liquid. In PBS containing 0.02% Proclin 300.
Host	Rabbit
Immunogen	This antibody was raised by immunising mouse with mouse CD27-human IgG1 Fc fusion protein. Isolated popliteal lymph node cells were then fused with P3U1 myeloma cells to produce stable hybridomas.
Isotype	IgGk
Purity Detail	Protein A affinity purified.
Source	This chimeric rabbit antibody was made using the variable domain sequences of the original Rat IgG2a 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)