

CD137L recombinant monoclonal antibody (014)

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

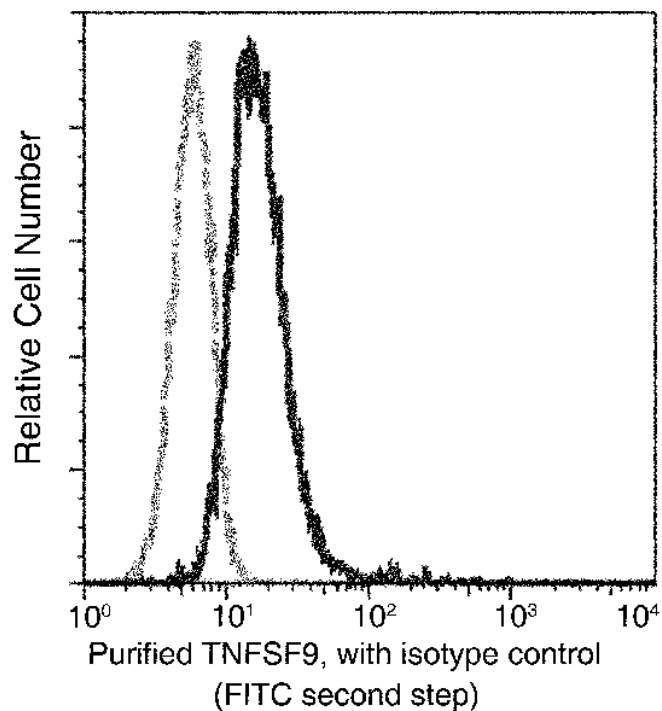
Ordering Information

[Order Online »](#)

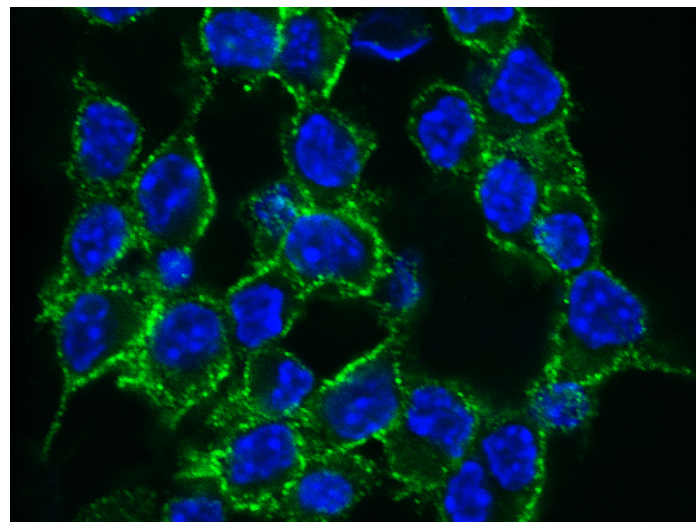
ENZ-ABS1005-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis of Mouse TNFSF9 expression on Raw264.7 cells. Cells were stained with purified anti-Mouse TNFSF9, 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.



Confocal immunofluorescence analysis of Mouse TNFSF9 in RAW264.7 cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with Rabbit anti-Mouse TNFSF9 monoclonal antibody (1:100) at 4 °C overnight. Then cells were stained with the Alexa Fluor® 488-conjugated Goat Anti-rabbit IgG secondary antibody (green) and counterstained with DAPI (blue). Positive staining was localized to plasma membrane.

Handling & Storage

Handling Protected from light. Avoid freeze-thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 4-1BBL, TNFSF 9

Application ELISA, Flow Cytometry, ICC, IF

Clone 014

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

Host Rabbit

Immunogen Recombinant mouse TNFSF9 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000)Immunocytochemistry (1:50-1:1'000)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](#)



Last modified: September 17, 2024

ENZO LIFE SCIENCES

Phone: 800.942.0430
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-
be@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