

TREM-2 recombinant monoclonal antibody (014)

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

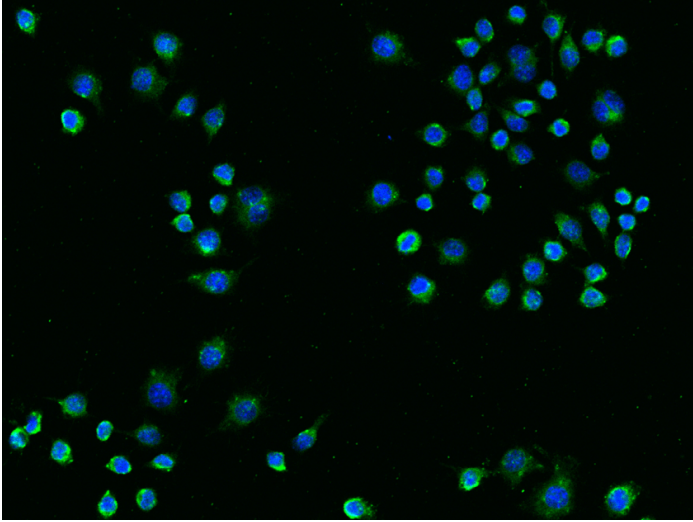
Ordering Information

[Order Online »](#)

ENZ-ABS1009-0100	100µl
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Immunofluorescence staining of mTREM2 in RAW264.7 cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-mouse mTREM2 monoclonal antibody (dilution ratio 1:60) 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 Cytoplasm and cell 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 Triggering receptor expressed on myeloid cells 2,
Triggering receptor expressed on monocytes 2

Application ELISA, ICC, IF

Clone 014

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

Host Rabbit

Immunogen Recombinant mouse TREM-2/TREM2 protein.

Isotype IgG

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Immunocytochemistry (1:20-1:100)Immunofluorescence (1:20-1:100)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: 800 642 0430
[info-](mailto:info-usa@enzolifesciences.com)
usa@enzolifesciences.com

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-](mailto:info-fr@enzolifesciences.com)
be@enzolifesciences.com

France
Phone: +33 472 440 655
[info-](mailto:info-fr@enzolifesciences.com)
fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
[info-](mailto:info-de@enzolifesciences.com)
de@enzolifesciences.com

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