

TREM-1 polyclonal antibody

Triggering Receptor Expressed on Myeloid cells-1 (TREM-1) is a receptor found on the surface of neutrophils and monocytes that amplifies inflammatory responses. It plays a role in sepsis, infectious diseases, and inflammatory conditions, making it a valuable target in immunology and inflammation research.

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

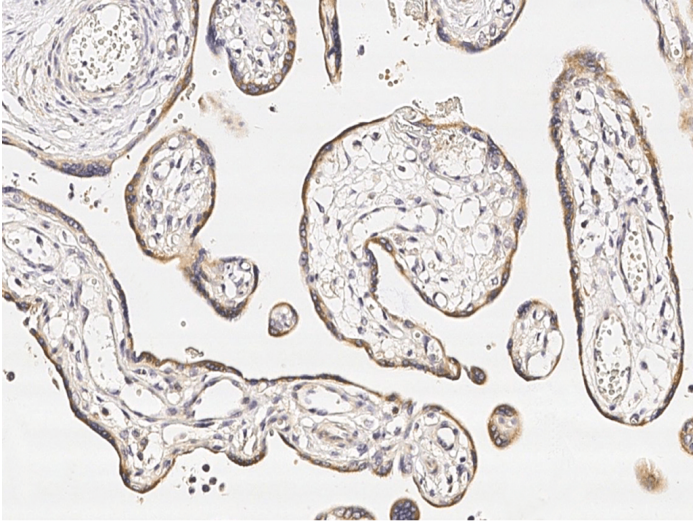
Ordering Information

[Order Online »](#)

| | |
|------------------|-------|
| ENZ-ABS1041-0100 | 100µl |
|------------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human TREM1 in human placenta with rabbit polyclonal antibody (1:1000, formalin-fixed paraffin embedded sections).

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 Triggering receptor expressed on myeloid cells 1,
Triggering receptor expressed on monocytes 1, CD354

Application ELISA, IHC (PS)

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

Host Rabbit

Immunogen Recombinant human TREM-1.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:500-1:2'000)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 [Worry-Free Guarantee](#)

Last modified: October 7, 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-be@enzolifesciences.com)

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