

CD118 monoclonal antibody (8E5E4D3)

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

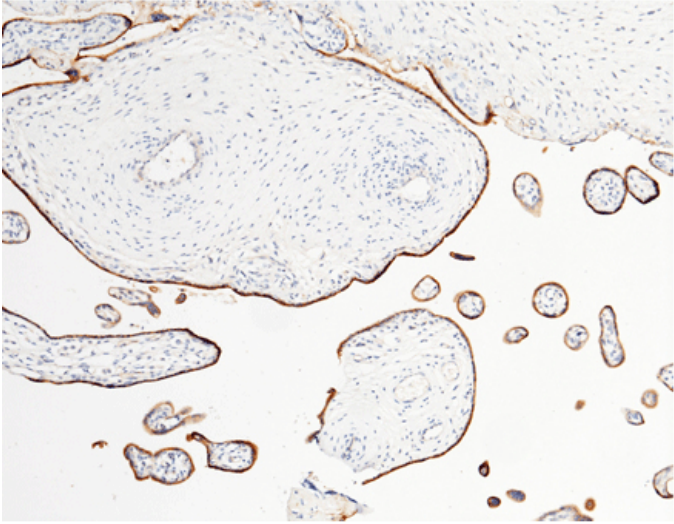
Ordering Information

[Order Online »](#)

| | |
|-----------------|-------|
| ENZ-ABS955-0100 | 100µl |
|-----------------|-------|

Manuals, SDS & CofA

[View Online »](#)



Immunochemical staining of human LIFR in human placenta with mouse monoclonal antibody (1:60, formalin-fixed paraffin embedded sections). Positive staining was localized to trophoblast.

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 LIFR, Leukemia inhibitory factor receptor

Application ELISA, IHC (PS)

Clone 8E5E4D3

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

Host Mouse

Immunogen Recombinant human LIFR protein.

Isotype IgG1

Purity Detail Protein A purified.

Recommendation Dilutions/Conditions ELISA (1:1'000-1:2'000)Immunohistochemistry (paraffin sections) (1:50-1:200)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](#).



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)