

CD105 polyclonal antibody

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

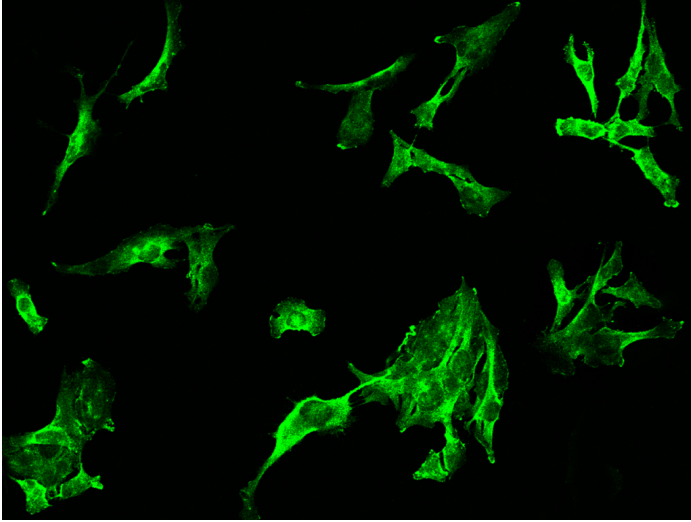
Ordering Information

[Order Online »](#)

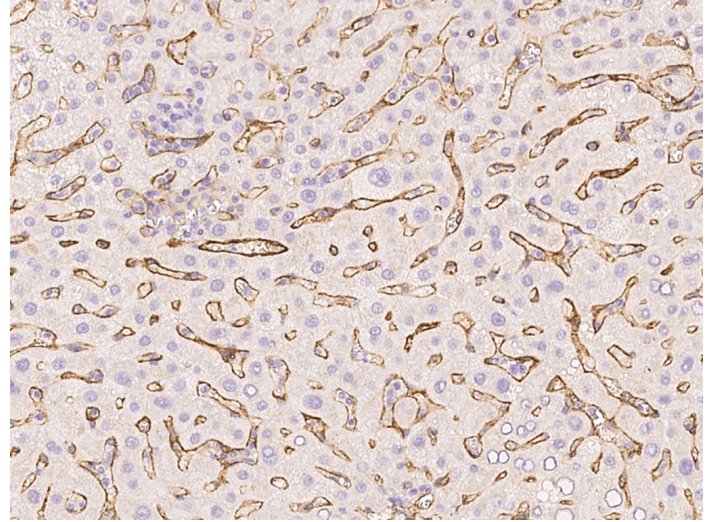
ENZ-ABS937-0100	100µl
-----------------	-------

Manuals, SDS & CofA

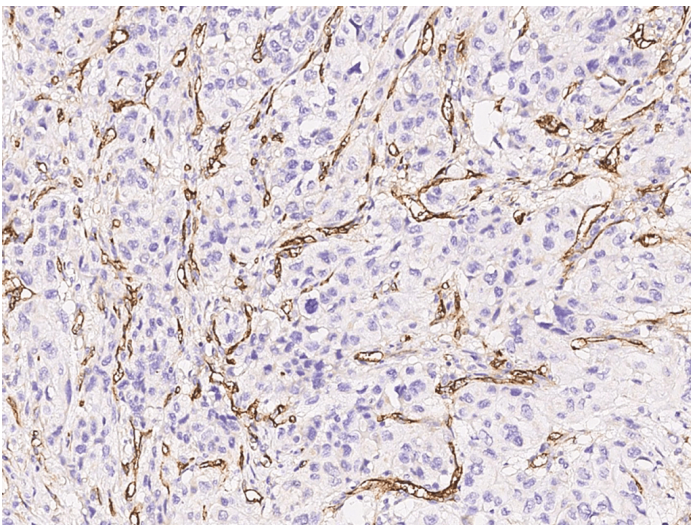
[View Online »](#)



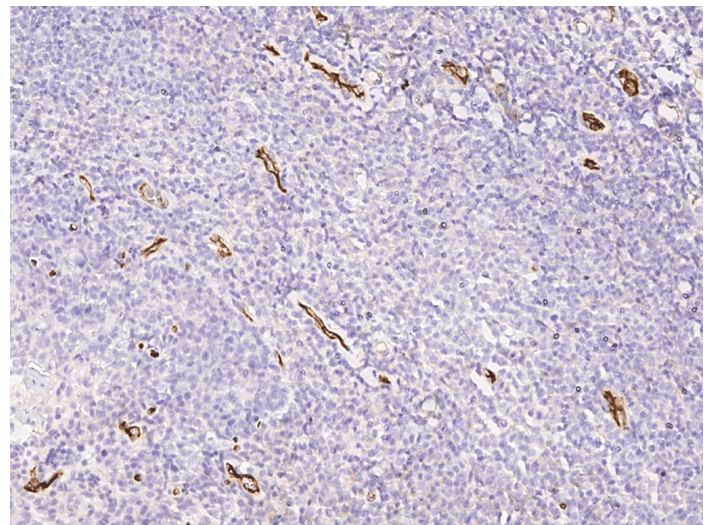
Immunofluorescence staining of CD105 in U251MG cells. Cells were fixed with 4% PFA, blocked with 10% serum, and incubated with rabbit anti-Human CD105 polyclonal antibody (dilution ratio 1:1000) at 4 °C overnight. Then cells were stained with the Alexa Fluor®488-conjugated Goat Anti-rabbit IgG secondary antibody (green). Positive staining was localized to cytoplasm and cell membrane.



Immunochemical staining of human CD105 in human liver with rabbit polyclonal antibody at 1:5000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human CD105 in human renal carcinoma with rabbit polyclonal antibody at 1:5000 dilution, formalin-fixed paraffin embedded sections.



Immunochemical staining of human CD105 in human tonsil with rabbit polyclonal antibody at 1:5000 dilution, formalin-fixed paraffin embedded sections.

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 Endoglin

Application ELISA, ICC, IF, IHC (PS)

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

Host Rabbit

Immunogen Recombinant human Endoglin/CD105 protein.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions ELISA (1:5'000-1:10'000)Immunohistochemistry (paraffin sections) (1:2'000-1:10'000)Immunocytochemistry (1:300-1:10'000)Immunofluorescence (1:300-1:10'000)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: September 17, 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-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

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

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

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