

ANGPTL4 polyclonal antibody

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

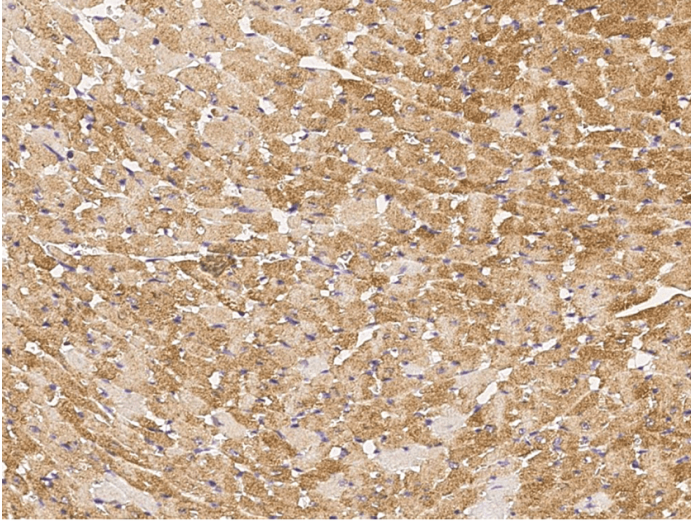
Ordering Information

[Order Online »](#)

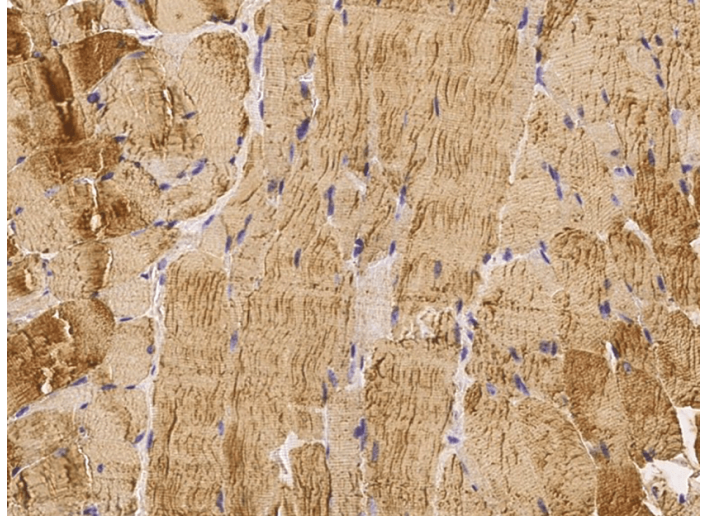
ENZ-ABS1015-0100	100µl
------------------	-------

Manuals, SDS & CofA

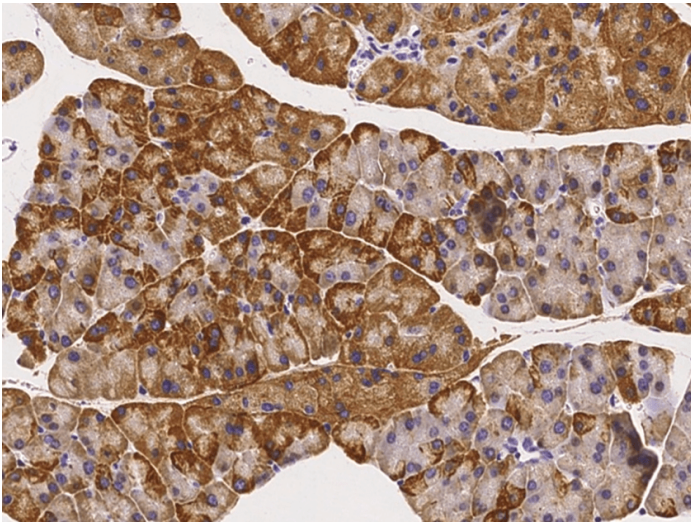
[View Online »](#)



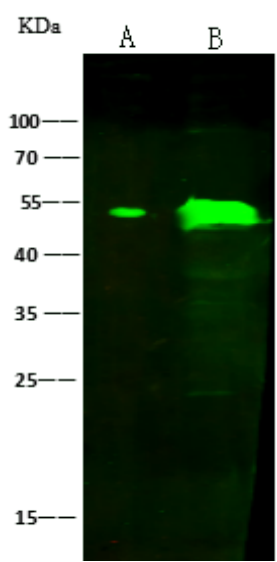
Immunohistochemical staining of mouse ANGPTL4 in mouse heart with rabbit polyclonal antibody (1:2000, formalin-fixed paraffin embedded sections).



Immunohistochemical staining of mouse ANGPTL4 in mouse skeletal muscle with rabbit polyclonal antibody (1:2000, formalin-fixed paraffin embedded sections).



Immunohistochemical staining of mouse ANGPTL4 in mouse pancreas with rabbit polyclonal antibody (1:2000, formalin-fixed paraffin embedded sections).



Anti-ANGPTL4 rabbit polyclonal antibody at 1:500 dilution. Lane A: Mouse liver tissue lysate Lane B: Mouse skeletal muscle tissue lysate. Lysates/proteins at 30 μ g per lane. Secondary Goat Anti-Rabbit IgG (H&L) /Dylight 800 at 1/10000 dilution. Developed using the Odyssey technique. Performed under reducing conditions. Observed band size:54 kDa.

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 Angiotensin-like protein 4

Application ELISA, IHC (PS), WB

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

Host Rabbit

Immunogen Recombinant mouse ANGPTL4 / angiotensin-like protein. 4 protein.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/Conditions Western Blot (1:500-1:1'000)Immunohistochemistry (paraffin sections) (1:1'000-1:5'000)ELISA (1:5'000-1:10'000)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,
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-
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)