ANGPTL4 polyclonal antibody

This antibody is covered by our Worry-Free Guarantee.

Ordering Information

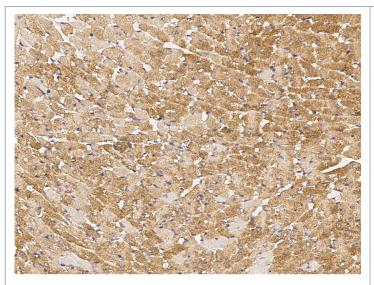
Order Online »

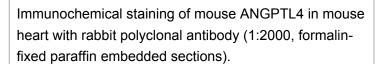
ENZ-ABS1015-0100

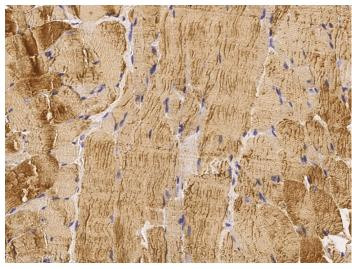
100µl

Manuals, SDS & CofA

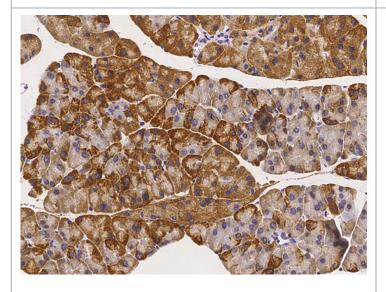
View Online »



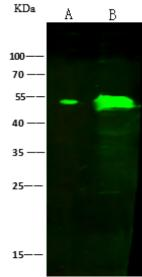




Immunochemical staining of mouse ANGPTL4 in mouse skeletal muscle with rabbit polyclonal antibody (1:2000, formalin-fixed paraffin embedded sections).



Immunochemical 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 µg 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 Angiopoietin-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 / angiopoietin-like protein.

4 protein.

Isotype IgG

Purity Detail Protein A and antigen affinity purified.

Recommendation Dilutions/ConditionsWestern 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 infousa@enzolifesciences.com European Sales Office ENZO LIFE SCIENCES (ELS) AG Phone: +41 61 926 8989 infoeu@enzolifesciences.com Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Germany Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com