

Adiponectin receptor 2 (mouse) polyclonal antibody

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

Manuals, SDS & CofA

[View Online »](#)

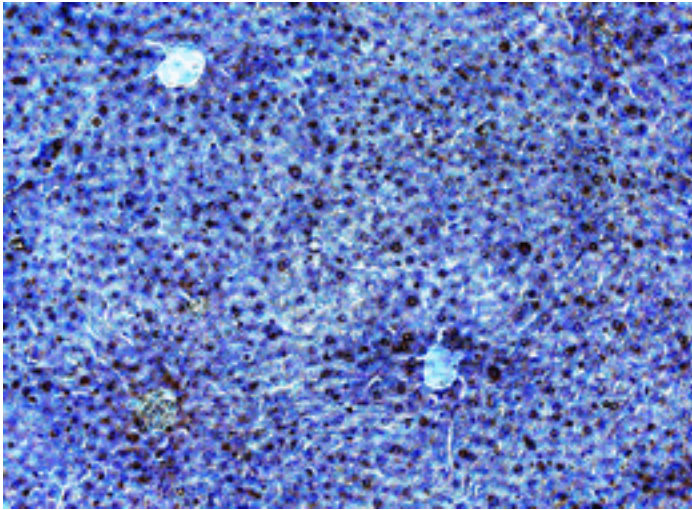


Figure 1: Immunohistochemistry on paraffin sections of mouse liver using PAb to Adiponectin Receptor 2 (mouse) (Prod. No. [ALX-210-647](#)).¹,Adiponectin receptor 2 (mouse) polyclonal antibody Immunohistochemistry,Adiponectin receptor 2 (mouse) polyclonal antibody Immunohistochemistry,yes ALX-210-647,210-

647_2.gif,http://static.enzolifesciences.com/fileadmin/files/image/210-647_2.gif,**Figure 2:** Immunohistochemistry on paraffin sections of mouse muscle using PAb to Adiponectin Receptor 2 (mouse) (Prod. No. [ALX-210-647](#)).¹

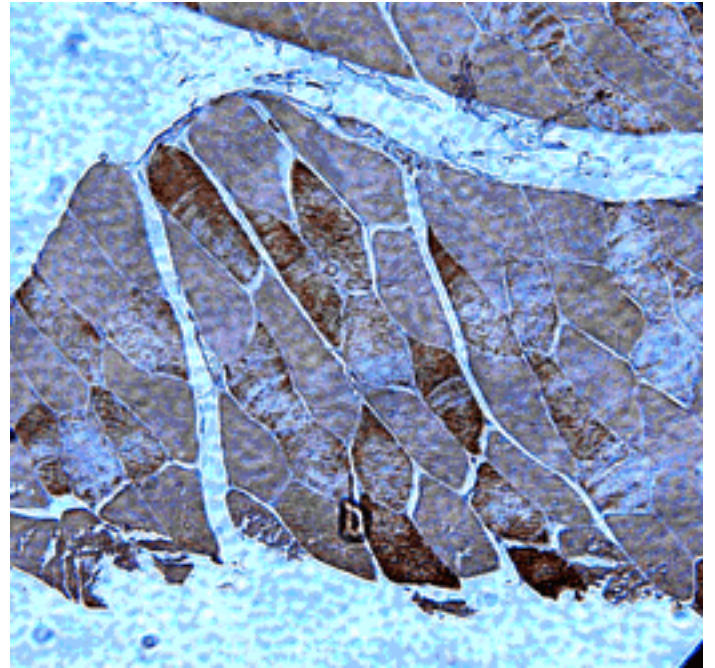


Figure 2: Immunohistochemistry on paraffin sections of mouse muscle using PAb to Adiponectin Receptor 2 (mouse) (Prod. No. ALX-210-647).

Handling & Storage

Handling	For maximum product recovery after thawing, centrifuge the vial before opening the cap. Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C.
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	AdipoR2
Application	ICC, IHC (PS), WB
Formulation	Liquid. In PBS containing 1mg/ml BSA and 0.1% sodium azide.
Host	Goat
Immunogen	Synthetic peptide corresponding to aa 11-40 (C11TRTPEPDIRLRKGHQLDDTRGSNNDNYQG40) of mouse adiponectin receptor 2 (AdipoR2).
Purity Detail	Epitope affinity-purified IgG.
Recommendation Dilutions/Conditions	Immunocytochemistry (acetone fixed cells, 1:400)Immunohistochemistry (paraffin sections, 1:250)Western Blot (1:1,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Mouse
UniProt ID	Q8BQS5

Worry-free Guarantee

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



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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

Last modified: May 29, 2024