

Adiponectin receptor 1 polyclonal antibody

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

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-210-6450	200µg
--------------	-------

Manuals, SDS & CofA

[View Online »](#)

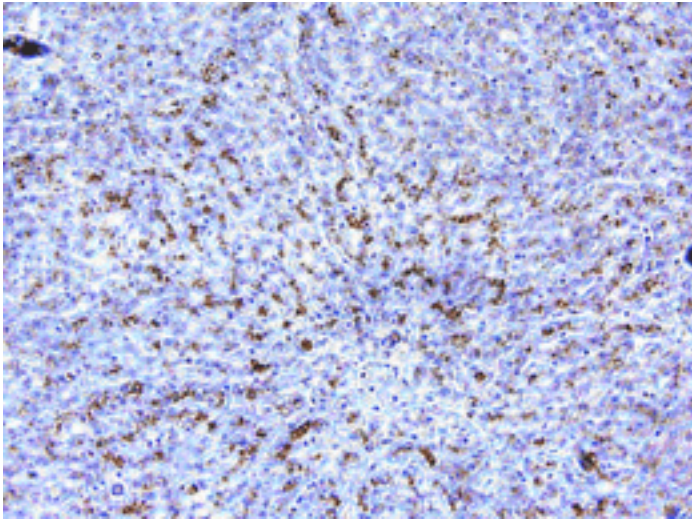


Figure 1: Immunohistochemistry on paraffin sections of human liver using PAb to Adiponectin Receptor 1 (Prod. No. [ALX-210-645](#)).¹, Adiponectin receptor 1 polyclonal antibody Immunohistochemistry, Adiponectin receptor 1 polyclonal antibody Immunohistochemistry, yes ALX-210-645, 210-

645_2.gif, http://static.enzolifesciences.com/fileadmin/files/image/210-645_2.gif, **Figure 2:** Immunohistochemistry on paraffin sections of human muscle using PAb to Adiponectin Receptor 1 (Prod. No. [ALX-210-645](#)).¹

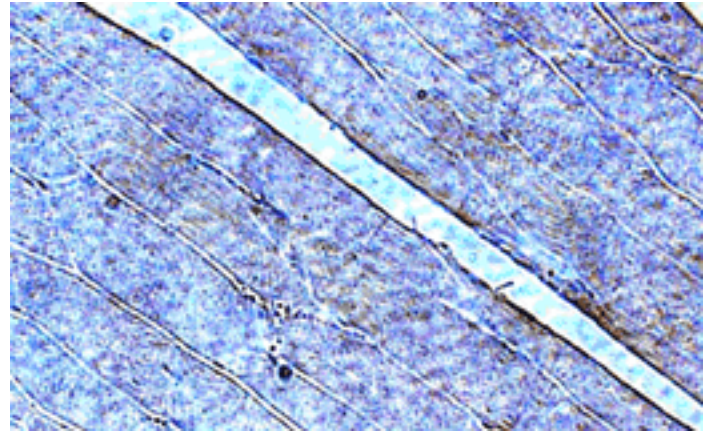


Figure 2: Immunohistochemistry on paraffin sections of human muscle using PAb to Adiponectin Receptor 1 (Prod. No. ALX-210-645).

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	AdipoR1, CGI-45
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 345-369 (V345AAAFVHFYGVSNLQEFRYGLEGGC369) of human, mouse or rat adiponectin receptor 1 (AdipoR1).
Purity Detail	Epitope affinity-purified.
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	Human, Mouse, Rat
UniProt ID	Q96A54

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