

# ACRP30headless (human) monoclonal antibody (Ne.Na)

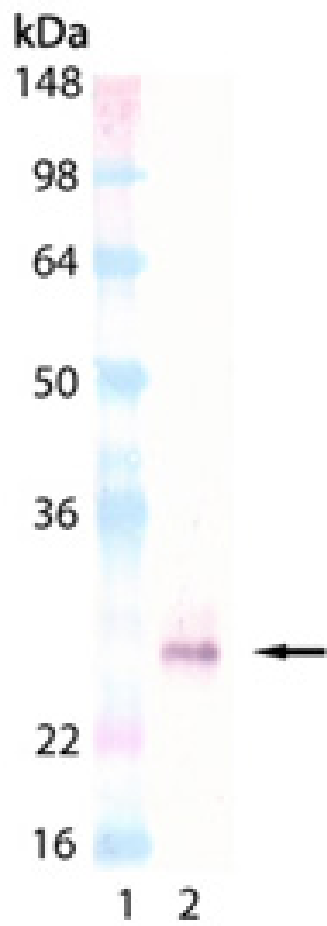
This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

Manuals, SDS & CofA

[View Online »](#)



Western blot analysis of ACRP30*headless* (human), mAb (Ne.Na) (Prod. No. ALX-804-144): Lane 1: MW Marker; Lane 2: ACRP30*headless* (human) (rec) (Prod. No. ALX-203-002).

## Handling & Storage

<b>Use/Stability</b>	Stable for at least 1 year after receipt when stored as recommended.
<b>Handling</b>	Avoid freeze/thaw cycles.
<b>Short Term Storage</b>	+4°C
<b>Long Term Storage</b>	-20°C
<b>Shipping</b>	Blue Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

<b>Alternative Name</b>	Adipocyte complement related protein of 30kDa (tail), apM1 (tail), Adiponectin (tail), AdipoQ (tail), GBP28 (tail)
<b>Application</b>	ELISA, IP, WB
<b>Clone</b>	Ne.Na
<b>Crossreactivity</b>	Does not cross-react with mouse ACRP30headless. Does not cross-react with full length adiponectin.
<b>Formulation</b>	Liquid. In PBS containing 0.02% sodium azide.
<b>Host</b>	Mouse
<b>Immunogen</b>	Recombinant human ACRP30headless (tail adiponectin).
<b>Isotype</b>	IgG1
<b>Recommendation Dilutions/Conditions</b>	ELISA (capture)Western Blot (1:1,000, colorimetric)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
<b>Source</b>	Purified from concentrated hybridoma tissue culture supernatant.
<b>Species Reactivity</b>	Human

UniProt ID

Q15848

Worry-free Guarantee

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

Last modified: May 29, 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-be@enzolifesciences.com](mailto:info-be@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)