

ACRP30headless (human) monoclonal antibody (Ne.Na)

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

Citations: 1

[View Online »](#)

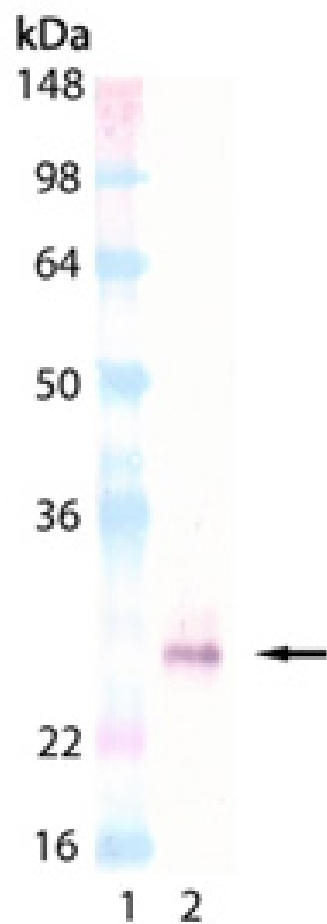
Ordering Information

[Order Online »](#)

ALX-804-144-C100	100µg
------------------	-------

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

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

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Adipocyte complement related protein of 30kDa (tail), apM1 (tail), Adiponectin (tail), AdipoQ (tail), GBP28 (tail)
Application	ELISA, IP, WB
Clone	Ne.Na
Crossreactivity	Does not cross-react with mouse ACRP30headless. Does not cross-react with full length adiponectin.
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Recombinant human ACRP30headless (tail adiponectin).
Isotype	IgG1
Recommendation Dilutions/Conditions	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.
Source	Purified from concentrated hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	Q15848

Worry-free Guarantee

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



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)