

ACRP30headless (mouse), (recombinant)

Control compound for rmsMEGACD40L[®] (Prod. No. ALX-522-120).

Ordering Information

[Order Online »](#)

ALX-203-003-C010	10µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

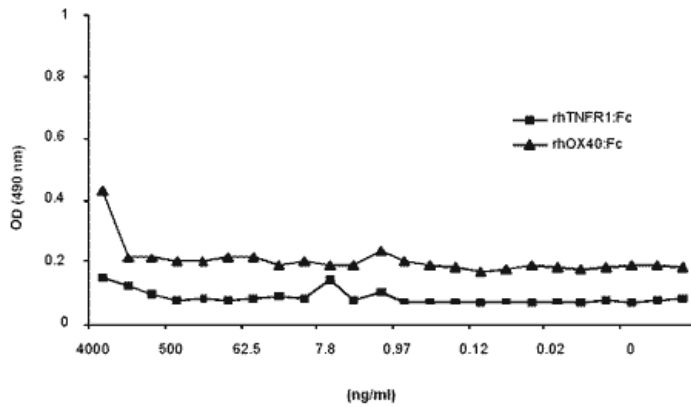


Figure: ACRP30*headless* (mouse) does not bind to rhTNF-R1 and rhOX40:Fc in an ELISA assay.

Method: A 96 well ELISA plate was coated O/N at RT with 100 μ l of 1 μ g/ml rhTNF-R1:Fc (Prod. No. [ALX-522-013](#)) and 1 μ g/ml rhOX40:Fc (Prod. No. ALX-522-022) in PBS. 100 μ l of rmsACRP30*headless* at 8 μ g/ml were mixed with 100 μ l 0.5% FCS in PBS in the first well and a series of 1:1 dilutions was prepared for the following wells. Binding of rmsACRP30*headless* to the receptor:Fc was detected after 1 hr incubation at 37°C using a goat anti-human HRP conjugated antibody (Jackson lab; 1:1⁰⁰⁰⁰)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored as recommended.
Handling	Avoid freeze/thaw cycles. After reconstitution, 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	Tail-Adp, Tail Adiponectin, Adipocyte complement related protein of 30kDa headless
Application Notes	Negative control for fusion proteins containing mouse ACRP <i>headless</i> .
Endotoxin Content	<0.1EU/μg protein (LAL-test).
Formulation	Lyophilized. Contains PBS.
MW	~25kDa
Purity	≥90% (SDS-PAGE)
Purity Detail	Purified by multi-step chromatography.
Reconstitution	Reconstitute with 100μl sterile water. Further dilutions should be made with medium containing 5% fetal calf serum or a carrier protein.
Source	Produced in HEK 293 cells. The collagen domain of mouse ACRP30 (aa 18-111) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Technical Info / Product Notes	FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	Q60994



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