

Adiponectin (mouse), (recombinant)

Homeostatic factor for regulating glucose levels, lipid metabolism, and insulin sensitivity

Citations: 12

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-059-C050	50µg
------------------	------

Manuals, SDS & CofA

[View Online »](#)

- Mimics serum adiponectin by forming high molecular weight (HMW) and hexameric species
- Expressed in mammalian cell line to guarantee proper post-translational modifications
- High purity and low endotoxin content to ensure experimental success

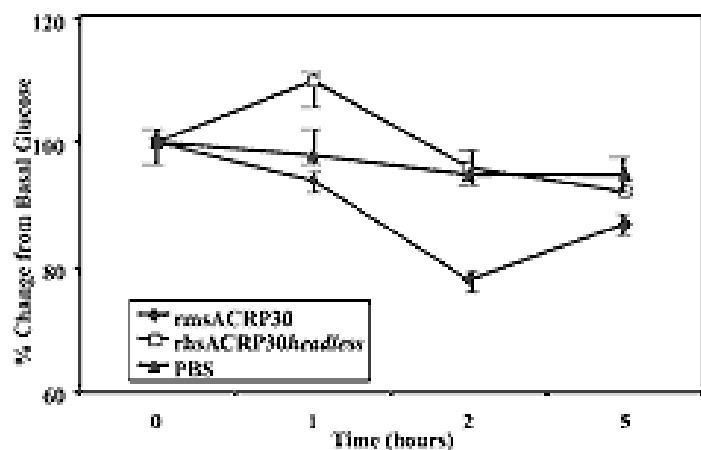


Figure 1: Bioactivity of rmsAdiponectin. **Method:** 2 groups of 4 mice each were injected with PBS (vehicle), rmAdiponectin (Prod. No. [ALX-522-059](#)) and rhACRP30 *headless* (tail adiponectin) (Prod. No. ALX-203-002) at 1µg/g body weight as a tail vein injection. Subsequently

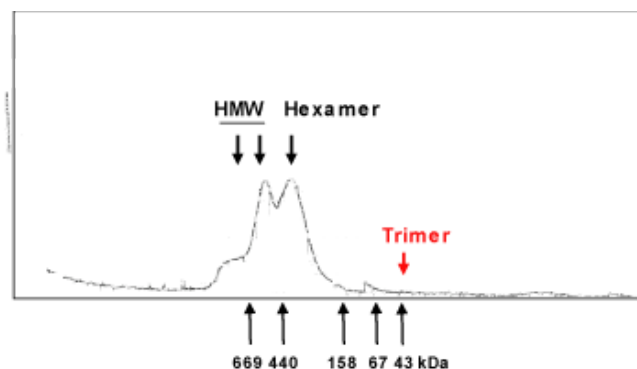


Figure 2: Superdex 200 analysis of Adiponectin (mouse) (recombinant) (Prod. No. [ALX-522-059](#)) produced in mammalian cells (HEK 293). The different forms of adiponectin are indicated by arrows: High molecular weight (HMW; representing 12-18-mer species) and hexamer. The red arrow indicates where should elute the trimeric state of adiponectin. *For further information on adiponectin forms, see Tsao et al., 2002 and Pajvani et al., 2003, in General Literature References.* “””, Adiponectin (mouse)

Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at -20°C.
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	Adipocyte complement related protein of 30kDa, ACRP30, AdipoQ, apM1, GBP28
Concentration	0.5mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~35kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 100µl sterile water.
Source	Produced in HEK 293 cells. Mouse adiponectin (aa 18-247) is fused at the N-terminus to a linker peptide (8 aa) and a FLAG [®] -tag.
Technical Info / Product Notes	<p>Historical data has shown that adiponectin inhibits serum glucose production in mice and protects mouse hepatocytes from TNF-α-induced cell death <i>in vitro</i>.</p> <p>FLAG is a registered trademark of Sigma-Aldrich Co.</p>
UniProt ID	Q60994

Last modified: May 29, 2024



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-fr@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