

Adiponectin (human), (recombinant)

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

Citations: 16

[View Online »](#)

Ordering Information

[Order Online »](#)

ALX-522-063-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

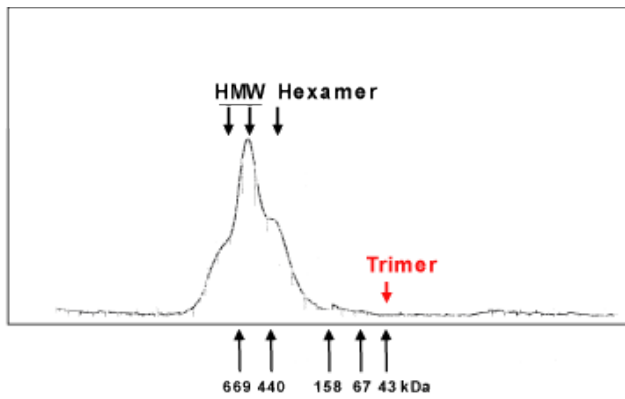


Figure: Superdex 200 analysis of Adiponectin (human) (recombinant) (Prod. No. [ALX-522-063](#)) produced in mammalian cells (HEK 293). The different forms of adiponectin are indicated by arrows: High molecular weight (HMW; representing 12-18-mer species) and hexamer. The red arrow indicates where should 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 (human)

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	1mg/ml after reconstitution.
Formulation	Lyophilized. Contains PBS.
MW	~33kDa (SDS-PAGE).
Purity	≥90% (SDS-PAGE)
Reconstitution	Reconstitute with 50µl sterile distilled water. Further dilutions should be made with medium containing 5% fetal calf serum or other carrier protein.
Source	Produced in HEK 293 cells. Human adiponectin (aa 16-244) is fused at the N-terminus to a linker peptide (14 aa) and a FLAG [®] -tag.
Technical Info / Product Notes	Historical results have shown that adiponectin activates AMPK. FLAG is a registered trademark of Sigma-Aldrich Co.
UniProt ID	Q15848



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