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 posttranslational modifications
- High purity and low endotoxin content to ensure experimental success

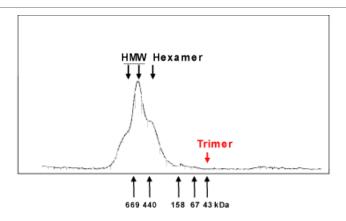


Figure: Superdex 200 analysis of Adiponectin (human) (recombinant) (Prod. No. ALX-522-063) produced in mammalian cells (HEK 293). The different forms of adiponectin are indicated by arrows: High molecular weight (HMW; representing 12-18-mer species) and hexamer. The red arrow indicates where should elute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

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 NotesHistorical results have shown that adiponectin activates

AMPK.

FLAG is a registered trademark of Sigma-Aldrich Co.

UniProt ID Q15848

