ACRP30headless (mouse), (recombinant)

Control compound for rmsMEGACD40 L^{\circledR} (Prod. No. ALX-522-120).

Ordering Information

Order Online »

ALX-203-003-C010

10µg

Manuals, SDS & CofA

View Online »

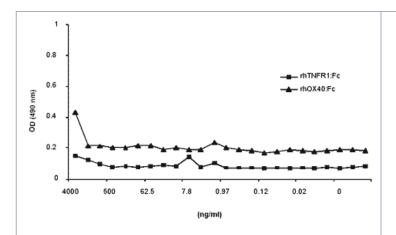


Figure: ACRP30*headless* (mouse) does not bind to rhTNF-R1 and rhOX40:Fc in an ELISA assay. **Method:** A 96 well ELISA plate was coated O/N at RT with 100 μl of 1 μg/ml rhTNF-R1:Fc (Prod. No. ALX-522-013) and 1 μg/ml rhOX40:Fc (Prod. No. ALX-522-022) in PBS. 100 μl of rmsACRP30*headless* at 8 μg/ml were mixed with 100 μl 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; 1:1"""

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 NotesNegative control for fusion proteins containing mouse

ACRPheadless.

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

Last modified: May 29, 2024



info-

eu@enzolifesciences.com