

ACRP30headless Monoclonal Antibody [NE.NA]

(Catalog No. A-0538)

Background

ACRP30 (adipocyte complement-related protein of 30kD) is a fat-derived secreted protein that is expressed only in adipose tissue. It is believed to be involved in lipid metabolism and glucose homeostasis. ACRP30 encodes a protein comparable to collagens X, VIII and the collagen-like complement factor C1q. Modifications in this gene have been linked to adiponectin deficiency. Multiple alternatively spliced variants encoding the same protein exists for this gene.

Description

Mouse monoclonal antibody to ACRP30headless, Clone NE.NA

Formulation

Liquid. In PBS containing 0.02% sodium azide.

Immunogen

Recombinant human ACRP30headless (tail adiponectin)

Specificity

Human

Isotype

IgG1

Purification

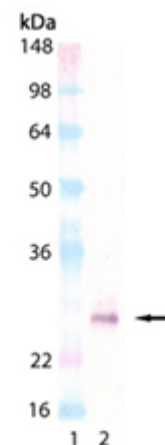
Serum

Storage

Store at 4°C for short term. Aliquot and store at -20°C for long term. Avoid multiple freeze/thaw cycles.

Application

ELISA, IP, WB (1:1000, colorimetric)



▲ Western blot analysis of ACRP30headless Monoclonal Antibody [NE.NA]; Lane 1: MW Marker; Lane 2: ACRP30headless (human) (rec).

Ordering Information

Products

ACRP30headless Monoclonal Antibody [NE.NA]

Size

100 µg

Cat. No.

A-0538-100

This product is for research purposes only. Not intended for use in diagnostic procedures.