ApoA-1 (human), (recombinant)

ApoA-1 participates in the reverse transport of cholesterol from tissues to the liver for excretion by promoting cholesterol efflux from tissues and by acting as a cofactor for the lecithin cholesterol acyltransferase (LCAT). ApoA-1 is also involved in activating spermatozoa motility.

Ordering Information

Order Online »

ENZ-PRT163-0010

10µg

Manuals, SDS & CofA

View Online »

Handling & Storage

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Apolipoprotein A-I, MGC117399

Formulation Lyophilized from 0.22µm filtered solution in PBS, pH 7.4.

Generally 5-8% mannitol or trehalose is added as a

protectant before lyophilization.

MW ~29 kDa

Purity ≥95% (SDS-PAGE)

Source Produced in HEK 293 cells.

UniProt ID P02647

Last modified: May 29, 2024



info-

eu@enzolifesciences.com