
LPA Polyclonal Antibody, HRP Conjugated

(Catalog # A50834)

Background

Apo(a) is the main constituent of lipoprotein(a) (Lp(a)). It has serine proteinase activity and is able of autoproteolysis. Inhibits tissue-type plasminogen activator 1. Lp(a) may be a ligand for megalin/Gp 330.

Description

LPA Polyclonal Antibody, HRP Conjugated. HRP. Raised in: Rabbit.

Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, PH 7.4.

Specificity

Human

Isotype

IgG

Uniprot ID

P08519

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Apolipoprotein(a) protein (3999-4112AA)

Storage

Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Apolipoprotein(a), Apo(a), Lp(a), LPA

Application

ELISA; Recommended dilution: ELISA 1:500-1:1000
