

---

## Apoe Polyclonal Antibody, FITC Conjugated

(Catalog # A53278)

---

### Background

Mediates the binding, internalization, and catabolism of lipoprotein particles. It can serve as a ligand for the LDL (apo B/E) receptor and for the specific apo-E receptor (chylomicron remnant) of hepatic tissues.

### Description

Apoe Polyclonal Antibody, FITC Conjugated. FITC. Raised in: Rabbit.

### Formulation

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

### Specificity

Mouse

### Isotype

IgG

### Uniprot ID

P08226

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Mouse Apolipoprotein E protein (19-311AA)

### 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

Apoe

### Application

ELISA; Recommended dilution: ELISA 1:100-1:500

---