

**Purified Anti-Human Apolipoprotein E
(ApoE) Antibody *E276***Catalog number: V101080
Unit size: 250 ug**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	apoE (APOE, AD2, APO-E, LDLCQ5, LPG, apolipoprotein E, ApoE4)
Clone	E276

Biological Properties

Preparation	Antibody purified by affinity chromatography
Application	Flow Cytometry (FACS)

Applications
