

**FITC Anti-Human Apolipoprotein B (ApoB)
Antibody *LDL 11***

Catalog number: V101125

Unit size: 250 µg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human, nhp
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	apoB (APOB, FLDB, LDLQC4, apoB-100, apoB-48, apolipoprotein B, FCHL2)
Clone	LDL 11
Conjugate	FITC

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with FITC under optimal conditions
Application	Flow Cytometry (FACS)

Spectral Properties

Conjugate	FITC
-----------	------

Applications
