

**Purified Anti-Human Apolipoprotein B  
(ApoB) Antibody \*LDL 20/17\***

Catalog number: V101115

Unit size: 250 µg

**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, nhp
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a
Immunogen	apoB (APOB, FLDB, LDLQC4, apoB-100, apoB-48, apolipoprotein B, FCHL2)
Clone	LDL 20/17

**Biological Properties**

---

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

**Applications**

---