

**Purified Anti-Human Apolipoprotein M
(ApoM) Antibody *M5***Catalog number: V101130
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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	apoM (APOM, G3a, HSPC336, NG20, apo-M, apolipoprotein M)
Clone	M5

Biological Properties

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

Applications
