

**Biotin Anti-Human Apolipoprotein M
(ApoM) Antibody *M7***

Catalog number: V101135

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
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG2a
Immunogen	apoM (APOM, G3a, HSPC336, NG20, apo-M, apolipoprotein M)
Clone	M7
Conjugate	Biotin

Biological Properties

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

Spectral Properties

Conjugate	Biotin
-----------	--------

Applications
