

**Biotin Anti-Human Apolipoprotein D (ApoD)
Antibody *D263***Catalog number: V101110
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 IgG1
Immunogen	apoD (APOD, Apod, apolipoprotein D)
Clone	D263
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