

**Purified Anti-Human Apolipoprotein D
(ApoD) Antibody *D544***Catalog number: V101105
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	apoD (APOD, Apod, apolipoprotein D)
Clone	D544

Biological Properties

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

Applications
