

**Purified Anti-Human Apolipoprotein J
(Clusterin) Antibody *J29***

Catalog number: V101095

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 IgG2b
Immunogen	apoJ (CLU, AAG4, APO-J, APOJ, CLI, CLU1, CLU2, CLU, AAG4, APO-J, APOJ, CLI, CLU1, CLU2, KUB1, NA1/NA2, SGP-2, SGP2, SP-40, TRPM-2, TRPM2, clusterin)
Clone	J29

Biological Properties

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

Applications
