

**Biotin Anti-Human Apolipoprotein J  
(Clusterin) Antibody \*J84\***

Catalog number: V101100

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 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	J84
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**

---