

PRODUCT INFORMATION

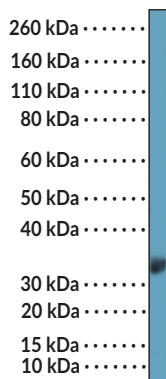
Clusterin (C-Term) Rabbit Monoclonal Antibody (Clone RM437)

Item No. 35616

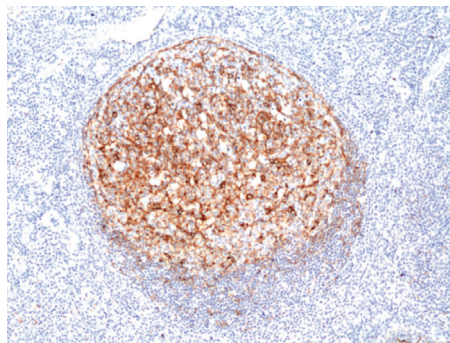
Overview and Properties

Contents:	This vial contains 100 µl of protein A-affinity purified monoclonal antibody.
Synonyms:	Apolipoprotein J, ApoJ, Ku70-binding Protein 1, SGP-2, Sulfated Glycoprotein 2, Testosterone-repressed Prostate Message 2, TRPM-2
Immunogen:	Peptide from the C-terminal region of clusterin
Cross Reactivity:	(+) Clusterin
Species Reactivity:	(+) Human
Form:	Liquid
Storage:	-20°C (as supplied)
Stability:	≥1 year
Storage Buffer:	PBS with 50% glycerol, 1% BSA, and 0.09% sodium azide
Clone:	RM437
Host:	Rabbit
Isotype:	IgG
Applications:	Immunohistochemistry (IHC) and Western blot (WB); the recommended starting dilution is 1:100-1:400 for IHC and 1:1,000-1:2,000 for WB. Other applications were not tested, therefore optimal working concentration/dilution should be determined empirically.

Images



WB of human brain tissue lysate using Clusterin (C-Term) Rabbit Monoclonal Antibody (Clone RM437) at a dilution of 1:1,000.



Immunohistochemical staining of formalin-fixed and paraffin-embedded human tonsil tissue using Clusterin (C-Term) Rabbit Monoclonal Antibody (Clone RM437) at a dilution of 1:100.

WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 01/17/2024

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

Clusterin is a secreted heterodimeric glycoprotein.¹ It is proteolytically processed to remove a signal peptide and cleave the α and β chains.^{1,2} This cleavage creates small- and medium-sized isoforms found in the nucleus or cytosol and mitochondria, respectively, and a large-sized isoform composed of α - and β chains linked *via* disulfide bonds near the center that is secreted outside the cell.² Clusterin is ubiquitously expressed in tissues and found in physiological fluids. It is a multifunctional protein involved in a wide range of processes, including immune modulation, autophagosome formation, apoptosis, and extracellular matrix remodeling, among others.³ Clusterin is upregulated by cellular stress and acts as an extracellular chaperone and scavenger to prevent protein aggregation. SNPs in *CLU*, the gene encoding clusterin, are associated with a higher rate of cognitive decline in patients with mild cognitive decline or Alzheimer's disease.⁴ Clusterin is also associated with inflammatory and cardiovascular diseases, where it has either protective or detrimental effects.² Cayman's Clusterin (C-Term) Rabbit Monoclonal Antibody (Clone RM437) can be used for immunohistochemistry (IHC) and Western blot (WB) applications.

References

1. Jones, S.E. and Jomary, C. Molecules in focus: Clusterin. *Int. J. Biochem. Cell Biol.* **34**(5), 427-31 (2002).
2. Rodríguez-Rivera, C., Garcia, M.M., Molina-Álvarez, M., *et al.* Clusterin: Always protecting. Synthesis, function and potential issues. *Biomed. Pharmacother.* **134**, 111174 (2021).
3. Rohne, P., Prochnow, H., and Kock-Brandt, C. The CLU-files: Disentanglement of a mystery. *Biomol. Concepts* **7**(1), 1-15 (2016).
4. Thambisetty, M., Beason-Held, L.L., An, Y., *et al.* Alzheimer risk variant CLU and brain function during aging. *Biol. Psychiatry* **73**(5), 399-405 (2013).