

Clusterin (human) monoclonal antibody (CLI-9)

Clusterin (SGP-2/sulfated glycoprotein-2; CLI; Apolipoprotein J (Apo J); TRPM-2) is the first identified extracellular mammalian chaperone and appears to be implicated in such diverse processes as sperm maturation, regulation of complement activation, programmed cell death, tissue remodelling and lipid transport. Increases in clusterin mRNA and protein have been consistently detected in apoptotic cell death paradigms, establishing clusterin gene expression as a prominent marker of apoptotic cell loss. A dual role has been shown for clusterin in that cells can be protected from damage and dead cells can be more easily removed after some types of cellular damage.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 3

[View Online »](#)

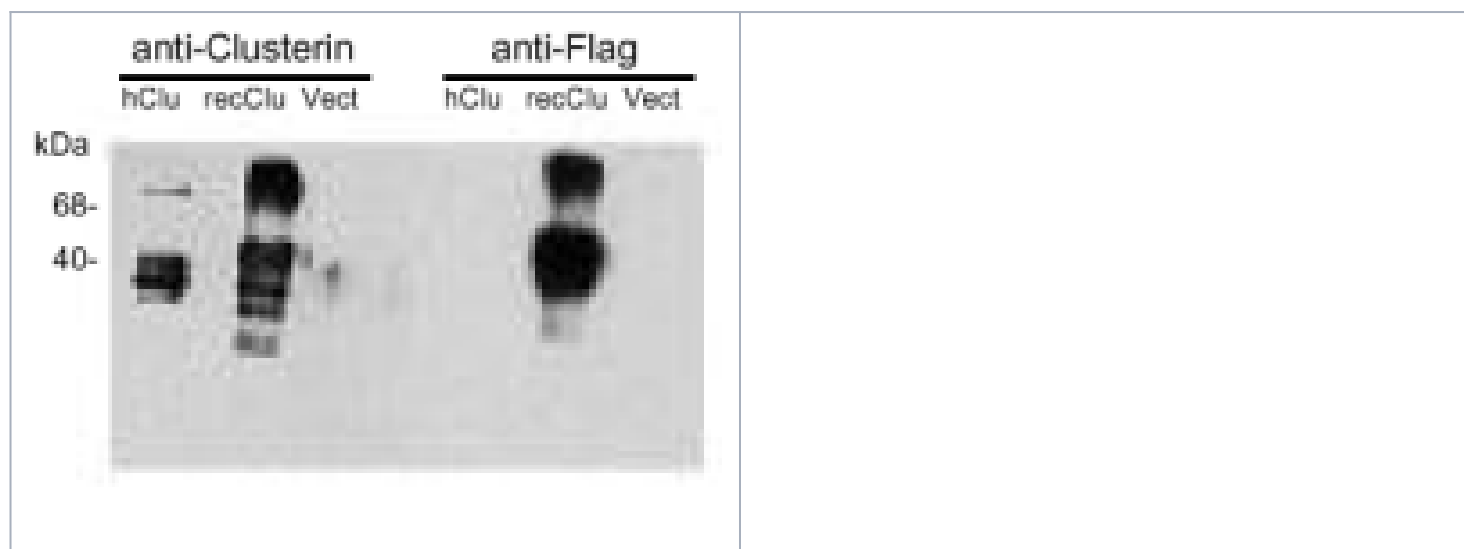
Ordering Information

[Order Online »](#)

ALX-804-126-C100	100µg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	Stable for at least 6 months after receipt when stored at +4°C undiluted.
Handling	Avoid freeze/thaw cycles.
Long Term Storage	+4°C
Shipping	Blue Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	TRPM-2, SGP-2, CLI, Apolipoprotein J, CLU, APO-J
Application	ICC, IHC (FS), IHC (PS), IP, WB
Clone	CLI-9
Formulation	Liquid. In PBS containing 0.02% sodium azide.
Host	Mouse
Immunogen	Purified human clusterin (TRPM-2).
Isotype	IgG1
Recommendation Dilutions/Conditions	Western Blot (1:2,000)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Source	Purified from hybridoma tissue culture supernatant.
Species Reactivity	Human
UniProt ID	P10909
Worry-free Guarantee	This antibody is covered by our Worry-Free Guarantee .



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)