# 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** 

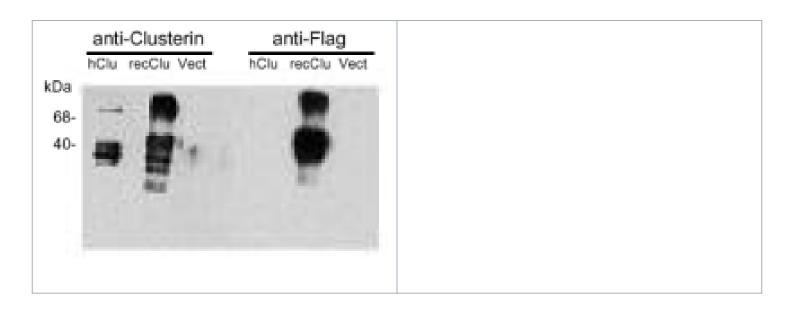
**Order Online** »

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).

lsotype lgG1

**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**

Last modified: May 29, 2024

This antibody is covered by our Worry-Free Guarantee



ENZO LIFE SCIENCES, INC. Phone: 800.942.0430 infousa@enzolifesciences.com

ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
infoeu@enzolifesciences.com

European Sales Office

Belgium, The Netherlands & Luxembourg Phone: +32 3 466 0420 infobe@enzolifesciences.com

France
Phone: +33 472 440 655
infofr@enzolifesciences.com

Phone: +49 7621 5500 526 infode@enzolifesciences.com UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
infouk@enzolifesciences.com