

Clusterin (secretory form) (human), (recombinant) (His-tag)

Clusterin is a secretory heterodimeric disulfide-linked glycoprotein containing 449 amino acids that is expressed in virtually all tissues, and found in all human fluids. It is involved in numerous physiological processes important for carcinogenesis and tumor growth, including apoptotic cell death, cell cycle regulation, DNA repair, cell adhesion, tissue remodeling, lipid transportation, membrane recycling, and immune system regulation. In addition, clusterin is up/ down regulated on the mRNA or protein level in many pathological and clinically relevant situations including cancer, organ regeneration, infection, Alzheimer disease, retinitis pigmentosa, myocardial infarction, renal tubular damage, autoimmunity and others.

Ordering Information

[Order Online »](#)

ALX-201-761-0010	10µg
ALX-201-761-0050	50µg

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	Lyophilized protein should be stored at $< -20^{\circ}\text{C}$, though stable at room temperature for 3 weeks. Reconstituted protein solution can be stored at $4-7^{\circ}\text{C}$ for 2-7 days. Working aliquots are stable for up to 3 months when stored at -20°C .
Handling	Avoid freeze/thaw cycles. After opening, prepare aliquots and store at -20°C .
Long Term Storage	-20°C
Shipping	Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	TRPM-2, Apolipoprotein J, APO-J, CLI, CLU, SGP-2
Endotoxin Content	$< 0.1\text{ng}/\mu\text{g}$ (1 IEU/ μg).
Formulation	Lyophilized. From a $0.2\ \mu\text{M}$ filtered solution of PBS, pH 7.4.
Purity	$\geq 90\%$ (reducing SDS-PAGE)
Reconstitution	Dissolve the lyophilized protein in 1X PBS. It is not recommended to reconstitute to a concentration less than $100\ \mu\text{g}/\text{ml}$.
Source	Human cells
UniProt ID	P10909

Last modified: May 29, 2024



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-
be@enzolifesciences.com](mailto:info-be@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)