

Rabbit Anti-Human Clusterin (Apolipoprotein J)

Instruction Manual

Catalog Number	PK-AB718-3856
Synonyms	Clusterin Antibody: Clusterin, Complement lysis inhibitor, CLI, Apolipoprotein J, ApoJ
Description	Clusterin, also known as Apolipoprotein J (ApoJ), is a ubiquitous multifunctional glycoprotein that can interact with a broad spectrum of molecules such as complement components, various receptors, and the Alzheimer's β -amyloid peptide. Clusterin expression is increased in Alzheimer's disease brain tissue and clusterin-immunoreactive amyloid plaques are found associated with phospho-tau-positive dystrophic neurites and it has been suggested that clusterin facilitates the conversion of diffuse β -amyloid deposits into amyloid and enhances tau phosphorylation in neurites around these plaques. Other reports show that clusterin expression is decreased in proliferating cells and is upregulated in quiescent and senescent cells, suggesting that it may also play a role in aging and tumorigenesis suppression. Clusterin exists in at least two distinct isoforms.
Quantity	100 μ g
Source / Host	Rabbit
Immunogen	Rabbit polyclonal Clusterin antibody was raised recombinant human Clusterin isoform 1 (Genbank accession No. NP_001822).
Purification Method	Affinity chromatography purified via peptide column.
Clone / IgG Subtype	Polyclonal antibody
Species Reactivity	Human
Specificity	
Formulation	Antibody is supplied in PBS containing 0.02% sodium azide.
Reconstitution	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 μ l or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
Storage & Stability	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
Applications	E, WB, IHC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually.
	Clusterin antibody can be used for the detection of Clusterin by Western blot at 0.5 - 1 μ g/mL. Antibody can also be used for immunohistochemistry starting at 10 μ g/mL.
Images	Available upon request.
References	Calero M, Rostagno A, Frangione B, et al. Clusterin and Alzheimer's disease. Subcell. Biochem. 2005; 38:273-98. Martin-Rehrmann MD, Hoe HS, Capuani EM, et al. Association of apolipoprotein J-positive β -amyloid plaques with dystrophic neurites in Alzheimer's disease brain. Neurotox. Res. 2005; 7:231-42. Trouwakos IP and Gonos ES. Clusterin/Apolipoprotein J in human aging and cancer. Int. J. Biochem. Cell Biol. 2002; 34:1430-48. Martinon F and Tschopp J. NLRs join TLRs as innate sensors of pathogens. TRENDS Imm. 2005; 26:447-54.
Images	Available upon request.
Related Products	Clusterin Peptide, Cat.No. PK-AB718-3856P Human brain Tissue Lysate, Cat. No. PK-AB718-1303 APP Antibody (AbNT), Cat. No. PK-AB718-2133 APP Antibody (CT), Cat. No. PK-AB718-2136

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.