

**Purified Mouse Anti-human Clusterin  
Antibody \*Hs-3, monoclonal, Cross  
Adsorbed\***Catalog number: V103340  
Unit size: 0.1 mg**Product Details**

|                    |                                                                             |
|--------------------|-----------------------------------------------------------------------------|
| Storage Conditions | 2-8°C with minimized light exposure. Do not freeze.                         |
| Expiration Date    | 12 months upon receiving                                                    |
| Concentration      | Lot specific (please consult certificate of analysis for given lot)         |
| Formulation        | Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA |

**Antibody Properties**

|                    |            |
|--------------------|------------|
| Species Reactivity | Human      |
| Class              | Primary    |
| Clonality          | Monoclonal |
| Host               | Mouse      |
| Immunogen          | Clusterin  |
| Clone              | Hs-3       |

**Biological Properties**

|             |                                                                                                                                                 |
|-------------|-------------------------------------------------------------------------------------------------------------------------------------------------|
| Preparation | Antibody purified by affinity chromatography, cross-adsorbed against dog, cat, pig, cow serum and then conjugated with under optimal conditions |
| Application | WB, IHC(P), ICC, ELISA                                                                                                                          |

**Applications**

Clusterin is a 61 kDa protein that can be found in the nucleus, protein-containing complex and intracellular membrane-bounded organelle of cells. It is sometimes referred to as Apolipoprotein J, complement cytolysis inhibitor, sulfated glycoprotein 2 and complement-associated protein SP-40,40. In humans, clusterin is the subject of intensive research in part because of the fact that it plays a role in the complement activation, classical pathway, positive regulation of intrinsic apoptotic signaling pathway and intrinsic apoptotic signaling pathway. Clusterin plays an important role in organismal processes, namely, microglial cell activation, antimicrobial humoral response and response to misfolded protein. It positively regulates apoptotic process, NF-kappaB transcription factor activity and protein-containing complex assembly. On the other hand, it also suppresses protein-containing complex assembly, release of cytochrome c from mitochondria and amyloid fibril formation. Clusterin binds to unfolded protein, protein-containing complex and ubiquitin protein ligase. It has been associated with vital functions like protein heterodimerization activity. Additionally, clusterin has been found to be involved in apoptotic process, neuronal signal transduction and neuron death.