

Purified Mouse Anti-pig Acrosin Antibody
ACR-2, monoclonal, Cross AdsorbedCatalog number: V103025
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	Pig
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Acrosin
Clone	ACR-2

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against human, dog, cow serum and then conjugated with under optimal conditions
Application	FC (QC TESTED), ICC, WB

Applications

Acrosin is a protein with a molecular weight of 53 kDa, expressed in the protein-containing complex, extracellular region and acrosomal matrix of cells. In humans, acrosin has been found to be involved in organismal processes, for instance, activation of adenylate cyclase activity, acrosome reaction and single fertilization. Acrosin has been closely linked to vital functions like amidase and serine-type endopeptidase activity. Acrosin recognizes zinc ion, fucose and protease. Sequencing of Acrosin has supported it contains a primary structural unit, the peptidase S1 domain.