

Purified Mouse Anti-human IgA Secretory Component Antibody *SC-05, monoclonal*

Catalog number: V1031410

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	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	IgA Secretory Component
Clone	SC-05

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IP, WB, IHC(P), IHC(F)

Applications

Polymeric immunoglobulin receptor is a transmembrane protein with a molecular weight of 97 kDa, located in the receptor complex, integral component of plasma membrane and extracellular space of cells. Sequencing of polymeric immunoglobulin receptor has supported it contains 7 conserved structural units: extracellular, cytoplasmic, Ig-like V-type 1, Ig-like V-type 2, Ig-like V-type 3, Ig-like V-type 4 and Ig-like V-type 5 domain. It takes part in processes such as organismal processes, in particular, receptor clustering, retina homeostasis and neutrophil degranulation. Polymeric immunoglobulin receptor has been thought to be involved with important functions like polymeric immunoglobulin receptor activity, and in addition, is the subject of intensive study due to the fact that it is a member of the Fc receptor signaling pathway and epidermal growth factor receptor signaling pathway.