

Purified Rabbit Anti-Streptavidin Antibody
Polyclonal

Catalog number: V102170

Unit size: 1.0 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	, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	Streptavidin

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	Flow Cytometry (FACS)

Applications

Streptavidin is a protein with a molecular weight of 22 kDa, located in the extracellular region of cells. Sequencing of Streptavidin has shown it contains a primary structural unit, the avidin-like domain. Streptavidin binds to biotin.