

Purified Rabbit Anti-V5 Antibody
Polyclonal

Catalog number: V102185

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	, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	V5

Biological Properties

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

Applications

Identifies the 14 amino acid long epitope tag, V5, derived from paramyxovirus P/V proteins (amino acid sequence: GKPIPPLLGLDST).