

HRP Rabbit Anti-VSV-g Antibody *Polyclonal*

Catalog Number: V102210

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
Formulation	HEPES – saline with 0.002% Thimerosal and stabilizer

Antibody Properties

Class	Primary
Clonality	Polyclonal
Host	Rabbit
Immunogen	VSV-g

Biological Properties

Appearance	Liquid
Preparation	Antibody purified by affinity chromatography and then conjugated with HRP under optimal conditions
Application	FACS

Spectral Properties

Conjugate	HRP
-----------	-----

Applications

Identifies the VSV-G epitope tag, derived from vesicular stomatitis virus glycoproteins (amino acid sequence: YTDIEMNRLGK).