

HRP Rabbit Anti-T7 Antibody *Polyclonal*

Catalog Number: V102180

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	T7

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 11 amino acid long peptide derived from T7 bacteriophage gene10.