

HRP Rabbit Anti-GST Antibody *Polyclonal*Catalog number: V102135
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	GST
Conjugate	HRP

Biological Properties

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

Applications

Identifies the 26 kDa polypeptide GST tag.