

HRP-labeled goat anti-rabbit IgG (H+L)

Catalog Number: 16793

Unit Size: 1 mg

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.01% thimerosal, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat

Biological Properties

Stabilizer	0.09% sodium azide, 0.2% (w/v) BSA
Appearance	Liquid
Preparation	Goat anti-rabbit IgG (H+L) is produced in goat with pooled total rabbit IgG. The antibody is conjugated with HRP under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC
Recommended Dilutions	Suggested dilutions are only guidelines; users should titrate the product for their specific assay using appropriate controls

Application	Recommended dilution
Flow Cytometry (FACS)	1-5 µg/mL
IF	2 µg/mL
IHC	1-10 µg/mL

Spectral Properties

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

Applications

The enzyme labels can be visualized by means of enzyme histochemical methods via chromogenic (producing color) reactions, or via fluorogenic (producing fluorescence) reactions. The most used enzyme is horse radish peroxidase (HRP). HRP can be visualized by chromogenic reactions such as diaminobenzidine (e.g., DAB), fluorescence (e.g., ADHP) or chemiluminescence (e.g., luminol). The highly purified and cross-absorbed purified IgGs are conjugated to HRP in high yield under a neutral condition to maximally retain the activities of both IgG and HRP using our proprietary Buccutite™ protein crosslinking technology.