

PE/Cy5 Goat Anti-Rabbit IgG (H+L) Antibody *Cross Adsorbed*

Catalog Number: 50355

Unit Size: 200 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	~0.2 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 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 and affinity purified with rabbit IgG coupled beads. The antibody is conjugated with PE/Cy5 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	PE (phycoerythrin)-Cy5
-----------	------------------------

Excitation Wavelength 565 nm

Emission Wavelength 666 nm

Applications

PE/Cy5 Goat Anti-Rabbit IgG (H+L) Antibody *Cross Adsorbed* is a high-performance secondary antibody designed to bind both heavy and light chains of rabbit immunoglobulin G (IgG). Conjugated to PE/Cy5, this antibody delivers exceptionally bright red fluorescence with excitation/emission maxima of ~565/666 nm. The PE/Cy5 fluorophore is prepared using AAT Bioquest's proprietary labeling technology to ensure superior signal intensity compared to similar antibodies from other sources. This antibody is ideal for flow cytometry, fluorescence microscopy, ELISA, and other immunoassays requiring sensitive and specific detection of rabbit IgG. Its affinity-purified formulation provides consistent performance, and its compatibility with 561 nm yellow-green laser lines and common bandpass filters (e.g., 670/30 nm) ensures seamless integration into standard fluorescence detection workflows.