

PE/Cy5 Goat Anti-human IgG (H+L) Antibody

Catalog Number: 50224

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	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

Biological Properties

Stabilizer	0.09% sodium azide, 0.2% (w/v) BSA
Appearance	Liquid
Preparation	Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG. 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/Cy5
-----------	--------

Excitation Wavelength 565 nm

Emission Wavelength 666 nm

Applications

AAT Bioquest's anti-human secondary antibodies have well-characterized specificity for human immunoglobulins and are useful in the detection, sorting or purification of its specified target. This PE/Cy5-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar PE/Cy5 goat anti-human IgG antibodies from other commercial sources, and thus can significantly increase assay sensitivities. Secondary antibodies offer increased versatility enabling users to use many detection systems (e.g. HRP, AP, fluorescence). They can also provide greater sensitivity through signal amplification as multiple secondary antibodies can bind to a single primary antibody. This antibody was purified through affinity chromatography and conjugated to PE/Cy5 (ex/em = 565/666 nm). It is compatible with the 561 nm laser and 670/30 nm bandpass filter (for example, as in the BD FACSJazz™).