

Biotinylated goat anti-human IgG (H+L) *Cross Adsorbed*

Catalog Number: 50266, 50267

Unit Size: 200 ug, 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), 2 mg/mL BSA

Unit Details

Reconstitution Volume	50266 (200 ug)	50267 (1 mg)
	200 uL ddH ₂ O	1 mL ddH ₂ O

Antibody Properties

Species Reactivity	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

Biological Properties

Stabilizer	2 mg/mL BSA
Appearance	Solid
Preparation	Goat anti-human IgG (H+L) is produced in goat with pooled total human IgG and affinity purified with human IgG coupled beads. The antibody is conjugated with Biotinylated under optimal conditions.
Application	Flow Cytometry (FACS), IF, IHC, ELISA, WB
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
ELISA	100 ng/mL
WB	1-10 µg/mL

Spectral Properties

Conjugate Biotin

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 Biotin-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. 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 biotin.