

Biotin Goat Anti-human IgG (H+L) Antibody
Cross AdsorbedCatalog number: 50266, 50267
Unit size: 200 ug, 1 mg**Product Details**

Storage Conditions	2-6°C and kept from light. To extend the shelf-life of this product, add an equal volume of glycerol to make a final concentration of approximately 50% glycerol and store at -20°C.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 2 mg/mL BSA

Unit Details

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

Antibody Properties

Species Reactivity	Human
Class	Secondary
Clonality	Polyclonal
Host	Goat

Biological Properties

Stabilizer	None
Appearance	White 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 purified IgG has a minimal cross-reaction to human, horse, mouse, rabbit and bovine IgG. The antibody is conjugated with Biotin under optimal condition.
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot
Soluble In	Water

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.