

Biotinylated goat anti-rabbit IgG (H+L)

Catalog number: 16794 Unit size: 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	16794 (1 mg)
Reconstitution Volume	1 mL ddH ₂ O
Antibody Properties	
Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat
Chemical Properties	
Molecular Weight	~150000
Biological Properties	
Stabilizer	None
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 Biotinylated under optimal condition.
Application	Immunofluorescence (IF), Flow Cytometry (FACS)
Soluble In	Water
Spectral Properties	
Conjugate	Biotinylated
Applications	

Tel: 408-733-1055 | Fax: 408-733-1304 | Email: support@aatbio.com | For Research Use Only (RUO) ©2020 AAT Bioquest, Inc. Last revised March 2020. For more information and tools, please visit https://www.aatbio.com Many biotin molecules can be conjugated to a secondary antibody with the additional advantage of streptavidin having a high affinity for biotin. Through this amplification step and having the streptavidin bound to labels such as HRP or fluorescent probes, proteins which are expressed at low levels are more likely to be detected. For most assays, the biotin-labeled secondary antibody is added first, sequentially followed by the streptavidin conjugated to a label e.g. HRP.