

Cy7® goat anti-rabbit IgG (H+L) *Cross Adsorbed*

Catalog Number: 16880

Unit Size: 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	1 mL ddH ₂ O
-----------------------	-------------------------

Antibody Properties

Species Reactivity	Rabbit
Class	Secondary
Clonality	Polyclonal
Host	Goat

Chemical Properties

Molecular Weight	~150000
------------------	---------

Biological Properties

Stabilizer	2 mg/mL BSA
Appearance	Solid
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 Cy7® 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	Cy7®
Excitation Wavelength	756 nm
Emission Wavelength	779 nm

Applications

AAT Bioquest's anti-rabbit secondary antibodies are affinity-purified antibodies with well-characterized specificity for rabbit immunoglobulins and are useful in the detection, sorting or purification of its specified target. This Cy7-labeled secondary antibody was prepared using AAT Bioquest's proprietary labeling technology. It demonstrated much brighter signal compared to the similar Cy7 goat anti-rabbit IgG antibodies from other commercial sources, 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.