

iFluor™ 633 goat anti-rabbit IgG (H+L)

Catalog number: 16638, 16808

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 16638 (200 ug) 16808 (1 mg)

Reconstitution Volume 200 uL ddH₂O 1 mL ddH₂O

Antibody Properties

Species Reactivity Rabbit

Class Secondary

Clonality Polyclonal

Host Goat

Chemical Properties

Molecular Weight ~150000

Biological Properties

Stabilizer None

Appearance Black 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 iFluor™ 633 under optimal

condition.

Application Immunofluorescence (IF), Flow Cytometry (FACS)

Soluble In Water

Spectral Properties

Conjugate iFluor™ 633

Excitation Wavelength 640 nm

Applications

AAT Bioquest's iFluor™ dyes are optimized for labeling proteins, in particular, antibodies. These dyes are bright, photostable and have minimal quenching on proteins. They can be well excited by the major laser lines of fluorescence instruments (e.g., 350, 405, 488, 555 and 633 nm). iFluor™ 633 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of ~638 nm and ~655 nm respectively. These spectral characteristics make them an excellent alternative to Alexa Fluor® 633 goat anti-rabbit IgG (H+L) conjugate (Alexa Fluor® is the trademark of Invitrogen).