

**iFluor® 840 goat anti-rabbit IgG (H+L)**

Catalog Number: 48062, 48063

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	<b>48062 (200 ug)</b>	<b>48063 (1 mg)</b>
	200 uL ddH <sub>2</sub> O	1 mL ddH <sub>2</sub> 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. The antibody is conjugated with iFluor® 840 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	iFluor® 840
Excitation Wavelength	836 nm
Emission Wavelength	879 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, 532-561, 633-647, and 808 nm). iFluor® 840 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of 836 nm and 879 nm, respectively. These unique spectral characteristics makes iFluor® 840 goat anti-rabbit IgG (H+L) conjugates ideal for various NIR imaging applications, including Western blotting, ELISA, protein arrays, tissue section imaging, and *in vivo* imaging.