

## iFluor™ 820 goat anti-rabbit IgG (H+L)

Catalog number: 48058, 48059 Unit size: 200 µg, 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	PBS, 2 mg/mL BSA	
Unit Details			
Unit	48058 (200 μg) 48059 (1 mg)		
Reconstitution Volume	200 μL ddH <sub>2</sub> O 1 mL ddH <sub>2</sub> O		
Antibody Properties			
Species Reactivity	Rabbit	Rabbit	
Class	Secondary		
Clonality	Polyclonal	Polyclonal	
Host	Goat	Goat	
Chemical Properties			
Molecular Weight	~150000		
<b>Biological Properties</b>			
Stabilizer	None	None	
Appearance	Green solid	Green 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™ 820 under optimal condition.		
Application	Flow Cytometry (FACS), ELISA, HC, Western Blot		
Soluble In	Water		
Spectral Properties			
Conjugate	iFluor™ 820	iFluor™ 820	
Excitation Wavelength	822 nm		

Tel: 408-733-1055 | Fax: 408-733-1304 | Email: support@aatbio.com | For Research Use Only (RUO) ©2021 AAT Bioquest, Inc. Last revised October 2021. For more information and tools, please visit https://www.aatbio.com

## Applications

AAT Bioquest's iFluor<sup>™</sup> 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<sup>™</sup> 820 goat anti-rabbit IgG (H+L) conjugate has fluorescence excitation and emission maxima of <sup>®</sup>822 nm and <sup>®</sup>850 nm, respectively. These unique spectral characteristics makes iFluor<sup>™</sup> 820 goat anti-rabbit IgG (H+L) conjugate har structure spectral characteristics makes iFluor<sup>™</sup> 820 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.