

# Rabbit Anti-ICG [Indocyanine Green] Antibody, Biotinylated

Catalog Number: 46840

Unit Size: 100 ug

## 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	PBS, BSA

## Unit Details

---

Reconstitution Volume	0.1 mL ddH <sub>2</sub> O
-----------------------	---------------------------

## Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	ICG coupled to carrier protein

## Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy5 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 647 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

## Spectral Properties

---

Conjugate	Biotinylated
-----------	--------------

## Applications

---

Rabbit Anti-ICG [Indocyanine Green] Antibody, Biotinylated is a polyclonal antibody that specifically binds to ICG-labeled molecules and is directly labeled with biotin. This biotinylated format allows sensitive signal detection using streptavidin- or avidin-linked systems in diverse assay types. The antibody may cross-react with other heptamethine cyanine dyes used in near-infrared imaging. It is ideal for signal amplification in low-abundance target detection, validating ICG conjugation, and enabling multiplexed assays in various applications.