

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 350-Labeled

Catalog Number: 46822

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 ddH2O
-----------------------	--------------

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	ICG coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy3 coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated molecules and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This direct labeling enables single-step fluorescence detection without requiring a secondary antibody. The antibody may cross-react with other related cyanine dyes commonly used in near-infrared imaging. Its blue emission avoids spectral interference with NIR fluorophores, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.