

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 555-Labeled

Catalog Number: 46828

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. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 555 nm. This direct labeling format enables one-step fluorescence detection of ICG-labeled molecules without the need for secondary antibodies. 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 performing multiplexed fluorescence detection with minimal spectral overlap.