

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 488-Labeled

Catalog Number: 46825

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 iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated molecules and is directly labeled with iFluor® 488, a green-fluorescent dye excited at around 490 nm. This configuration enables rapid, single-step fluorescence detection without requiring a secondary antibody. The antibody may cross-react with other heptamethine cyanine dyes used in near-infrared imaging. Its green emission is spectrally distinct from NIR dyes, making it ideal for signal amplification in low-abundance target detection, conjugation verification, and multiplexed fluorescence workflows.