

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 750-Labeled

Catalog Number: 46830

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 Cy5 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® 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation around 750 nm. This direct labeling format supports sensitive, single-step detection of ICG-labeled targets without the need for secondary antibodies. The antibody may cross-react with other heptamethine cyanine dyes used in NIR fluorescence imaging. It is ideal for signal amplification in low-abundance target detection, validating ICG conjugation, and performing multiplexed fluorescence assays in the near-infrared spectrum.