

Rabbit Anti-ICG [Indocyanine Green] Antibody

Catalog Number: 46820

Unit Size: 1 mg

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, 0.05% Sodium Azide

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	ICG coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with DNP 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)

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody is a polyclonal antibody that specifically recognizes ICG-conjugated molecules. It enables sensitive detection of ICG-labeled antibodies, proteins, and probes in immunoassays, imaging, and other fluorescence-based applications. The antibody may also exhibit cross-reactivity with related heptamethine cyanine dyes commonly used in near-infrared (NIR) platforms. It is ideal for signal amplification in low-abundance target detection, validating ICG dye conjugation, and supporting multiplexed NIR detection workflows.