

Rabbit Anti-ICG [Indocyanine Green] Antibody

Catalog Number: 46820

Unit Size: 1 mg

Product Details

Storage Conditions	Minimize light exposure, Refrigerated (2-8 °C)
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	PBS

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

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.