

Rabbit Anti-ICG [Indocyanine Green] Antibody, Biotinylated

Catalog Number: 46840

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

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, Biotinylated is a polyclonal antibody that specifically binds to ICG-labeled molecules and is directly labeled with biotin. This biotinylated format allows sensitive signal detection using streptavidin- or avidin-linked systems in diverse assay types.

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 enabling multiplexed assays in various applications.