

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 555-Labeled

Catalog Number: 46828

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	N/A
------------------	-----

Soluble In	Water
------------	-------

Spectral Properties

Excitation Wavelength	557 nm
-----------------------	--------

Emission Wavelength	570 nm
---------------------	--------

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 555 nm. This direct labeling format enables one-step fluorescence detection of ICG-labeled molecules without the need for secondary antibodies.

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 performing multiplexed fluorescence detection with minimal spectral overlap.