

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 350-Labeled

Catalog Number: 46822

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	345 nm
-----------------------	--------

Emission Wavelength	450 nm
---------------------	--------

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated molecules and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This direct labeling enables single-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with other related cyanine dyes commonly used in near-infrared imaging. Its blue emission avoids spectral interference with NIR fluorophores, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.