

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 750-Labeled

Catalog Number: 46830

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	757 nm
-----------------------	--------

Emission Wavelength	779 nm
---------------------	--------

Applications

Rabbit Anti-ICG [Indocyanine Green] Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to ICG-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation around 750 nm. This direct labeling format supports sensitive, single-step detection of ICG-labeled targets without the need for secondary antibodies.

The antibody may cross-react with other heptamethine cyanine dyes used in NIR fluorescence imaging. It is ideal for signal amplification in low-abundance target detection, validating ICG conjugation, and performing multiplexed fluorescence assays in the near-infrared spectrum.