

Rabbit Anti-Fluorescein [FITC] Antibody, iFluor® 488-Labeled

Catalog Number: 46661

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	491 nm
-----------------------	--------

Emission Wavelength	516 nm
---------------------	--------

Applications

Rabbit Anti-Fluorescein [FITC] Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC and is directly labeled with iFluor® 488, a green-emitting fluorophore with excitation around 490 nm. This labeling format allows for sensitive, one-step fluorescence detection of FITC-conjugated targets without requiring a secondary antibody.

The antibody may cross-react with related dyes including Alexa Fluor® 488, DyLight® 488, and other xanthene-based fluorophores. It is ideal for signal amplification in detecting low-abundance targets, verifying fluorescein conjugation, and enabling multiplex fluorescence assays in applications such as flow cytometry, microscopy, and ELISA.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.