

Rabbit Anti-Fluorescein [FITC] Antibody

Catalog Number: 46750

Unit Size: 1 mg

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-Fluorescein [FITC] Antibody is a polyclonal antibody that specifically binds to fluorescein-based dyes such as FITC (fluorescein isothiocyanate), a widely used green fluorophore in molecular biology. This antibody enables detection of FITC-labeled biomolecules including antibodies, oligonucleotides, and proteins.

It may cross-react with structurally related dyes like Alexa Fluor® 488, and DyLight® 488. The antibody is ideal for signal amplification in applications where FITC-labeled targets are present at low levels, and is suitable for use in various techniques.

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