

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 555-Labeled

Catalog Number: 46598

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	557 nm
-----------------------	--------

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

Applications

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission maxima of ~556/568 nm. This configuration allows for one-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with similar fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. The spectral distinctness of iFluor® 555 makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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