

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled

Catalog Number: 46512

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

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

Applications

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima around 749/775 nm. The direct labeling enables single-step detection of Alexa Fluor® 488-labeled molecules without the need for a secondary antibody.

The antibody may cross-react with structurally related green fluorophores such as iFluor® 488, ATTO 488, and rhodamine 110. With its long-wavelength emission, iFluor® 750 is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows..

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.