

Rabbit Anti-Alexa Fluor 488 Antibody, iFluor® 555-Labeled

Catalog Number: 46505

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-Alexa Fluor 488 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 488-conjugated targets and is directly labeled with iFluor® 555, a bright, orange-red fluorophore excited at 550–560 nm. This direct labeling allows for single-step detection of Alexa Fluor 488-tagged biomolecules without requiring a secondary antibody.

The antibody may cross-react with similar green fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. Its emission in the orange-red spectrum enables spectral separation from Alexa Fluor® 488, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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