

Rabbit Anti-Alexa Fluor 555 Antibody

Catalog Number: 46530

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-Alexa Fluor 555 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated biomolecules. It offers high-affinity detection of Alexa Fluor® 555-labeled proteins, nucleic acids, or other tagged targets in immunoassays and imaging workflows.

The antibody may cross-react with spectrally similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. It is compatible with standard secondary antibody systems and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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