

Rabbit Anti-Alexa Fluor 488 Antibody

Catalog Number: 46500

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 488 Antibody is a polyclonal antibody that binds to conjugates labeled with Alexa Fluor® 488 dye. It offers high-affinity binding, making it ideal for applications requiring signal amplification or dye conjugation verification.

The antibody may cross-react with similar fluorophores such as iFluor® 488, ATTO® 488, and rhodamine 110. It 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.