

Rabbit Anti-Alexa Fluor 750 Antibody

Catalog Number: 46590

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 750 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules. It is ideal for detecting Alexa Fluor® 750-labeled antibodies, proteins, or probes in fluorescence-based applications.

The antibody may cross-react with structurally related fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. It is compatible with techniques including Western blotting, immunofluorescence microscopy, ELISA, and flow cytometry.

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