

Rabbit Anti-Alexa Fluor 647 Antibody

Catalog Number: 46560

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 647 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules. It is ideal for detecting Alexa Fluor® 647-labeled antibodies, proteins, or probes.

The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO® 647 and DyLight® 650. It is compatible with applications including flow cytometry, immunofluorescence, Western blotting, and ELISA.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.