

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 350-Labeled

Catalog Number: 46592

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	345 nm
-----------------------	--------

Emission Wavelength	450 nm
---------------------	--------

Applications

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitable blue fluorophore with excitation/emission of ~346/442 nm. This direct labeling format enables one-step detection of Alexa Fluor® 750-labeled targets without the need for secondary antibodies.

The antibody may cross-react with structurally related dyes such as iFluor® 750, ATTO® 740, and DyLight® 755. Its blue emission is well separated from near-infrared signals, making it ideal for multiplex assays and low-background imaging workflows.

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