

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 555-Labeled

Catalog Number: 46543

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	557 nm
-----------------------	--------

Emission Wavelength	570 nm
---------------------	--------

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation around 555 nm. This direct labeling format enables single-step detection, improving signal sensitivity and reducing background interference.

The antibody may cross-react with structurally similar dyes such as iFluor® 555, ATTO 550, and Cy3. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation, and performing multiplex fluorescence assays.

Alexa Fluor® and Cy3 are registered trademarks of Thermo Fisher Scientific.