

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 488-Labeled

Catalog Number: 46535

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	491 nm
Emission Wavelength	516 nm

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 488, a bright green-emitting fluorophore with excitation/emission maxima around 491/517 nm. This direct labeling format eliminates the need for a secondary antibody, enabling fast and simplified detection of Alexa Fluor® 555-labeled biomolecules.

The antibody may cross-react with structurally related dyes such as iFluor® 555, ATTO 550, and Cy3. Its green emission makes it compatible with multicolor fluorescence assays that require distinct spectral separation from red and far-red fluorophores. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.