

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 350-Labeled

Catalog Number: 46532

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 555 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 350, a UV-excitable blue fluorophore with excitation/emission maxima around 346/442 nm. This direct format enables one-step fluorescence detection of Alexa Fluor® 555 without the need for a secondary antibody.

The antibody may cross-react with related fluorophores such as iFluor® 555, ATTO 550, and Cy3. Its spectral separation from Alexa Fluor® 555 makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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