

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled

Catalog Number: 46542

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	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima of ~752/778 nm. This direct labeling enables single-step fluorescence detection without the need for secondary antibodies.

The antibody may cross-react with similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. The near-infrared emission of iFluor® 750 allows for low-background detection and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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