

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 647-Labeled

Catalog Number: 46538

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	656 nm
Emission Wavelength	670 nm

Applications

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. This direct format eliminates the need for secondary antibodies, allowing one-step detection of Alexa Fluor® 555-labeled molecules.

The antibody may cross-react with similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. Its far-red emission makes it ideal for multicolor fluorescence workflows by avoiding overlap with orange-emitting dyes. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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