

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 647-Labeled

Catalog Number: 46622

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 350 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. This configuration allows for one-step fluorescence detection without the need for secondary antibodies.

The antibody may cross-react with structurally similar fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. The far-red emission of iFluor® 647 avoids spectral overlap with blue fluorophores, making this reagent especially effective in signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

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