

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 750-Labeled

Catalog Number: 46572

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 647 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission of ~752/778 nm. This direct labeling format allows single-step detection of Alexa Fluor® 647-labeled targets without requiring secondary antibodies.

The antibody may cross-react with structurally similar fluorophores such as iFluor® 647, Cy5, ATTO 647, and DyLight® 650. Its near-infrared emission reduces background and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection

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