

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

Catalog Number: 46602

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 750 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. This direct labeling allows for one-step detection of Alexa Fluor® 750 targets without the need for a secondary antibody.

The antibody may cross-react with similar near-infrared fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. Its spectral compatibility makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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