

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

Catalog Number: 46573

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 647 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation around 656 nm and emission near 670 nm. This direct labeling format enables one-step fluorescence detection with high sensitivity and minimal background.

The antibody may cross-react with structurally related fluorophores such as iFluor® 647, Cy5, ATTO® 647, and DyLight® 650. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.