

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

Catalog Number: 46603

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 750 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation around 752 nm and emission near 778 nm. This direct labeling format allows single-step fluorescence detection with high sensitivity and minimal background.

The antibody may cross-react with structurally similar dyes such as iFluor® 750, ATTO 740, and DyLight® 755. 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.