

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 750-Labeled

Catalog Number: 46632

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 350 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission at around 752/778 nm. This direct format allows single-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with structurally related fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its NIR emission provides high signal-to-noise ratios and is well-suited for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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