

## 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 multicolor fluorescence imaging, deep-tissue analysis, and high-sensitivity detection workflows.

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