

## 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. ATTO® is a registered trademark of ATTO-TEC GmbH.*