

## Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 555-Labeled

Catalog Number: 46598

Unit Size: 100 ug

### Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, BSA

### Unit Details

---

Reconstitution Volume	0.1 mL ddH <sub>2</sub> O
-----------------------	---------------------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 750 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 750 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 555 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

### Applications

---

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission maxima of ~556/568 nm. This configuration allows for one-step fluorescence detection without the need for a secondary antibody. The antibody may cross-react with similar fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. The spectral distinctness of iFluor® 555 makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*