

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

Catalog Number: 46542

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 ddH2O
-----------------------	--------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 555 coupled to carrier protein

### Biological Properties

---

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

### Spectral Properties

---

Conjugate	iFluor® 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

### Applications

---

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission maxima of ~752/778 nm. This direct labeling enables single-step fluorescence detection without the need for secondary antibodies. The antibody may cross-react with similar fluorophores such as iFluor® 555, ATTO 550, and Cy3. The near-infrared emission of iFluor® 750 allows for low-background detection and is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.