

## Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 555-Labeled

Catalog Number: 46628

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, 0.2% (w/v) BSA

### Unit Details

---

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

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 350 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 350 coupled carrier protein in rabbit and affinity purified with Protein G 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 350 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 555, a bright orange fluorophore with excitation/emission at around 555/565 nm. This direct labeling format enables one-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its orange emission is spectrally distinct from Alexa Fluor® 350, making it ideal for multicolor imaging, flow cytometry, and other fluorescence-based assays.

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