

## Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 555-Labeled

Catalog Number: 46568

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 647 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 647 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 647 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with iFluor® 555, a bright orange-red fluorophore with excitation/emission of ~555/565 nm. This direct format enables one-step detection of Alexa Fluor® 647-labeled targets without the need for secondary antibodies.

The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO 647, and DyLight® 650. The orange-red fluorescence of iFluor® 555 is spectrally distinct from far-red dyes, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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