

## Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 647-Labeled

Catalog Number: 46622

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 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 A beads. The antibody is conjugated with iFluor® 647 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 647
Excitation Wavelength	656 nm
Emission Wavelength	670 nm

### Applications

---

Rabbit Anti-Alexa Fluor 350 Antibody, iFluor® 647-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 647, a far-red fluorophore with excitation/emission maxima of ~656/670 nm. This configuration allows for one-step fluorescence detection without the need for secondary antibodies. The antibody may cross-react with structurally similar fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. The far-red emission of iFluor® 647 avoids spectral overlap with blue fluorophores, making this reagent especially effective in signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows. *Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.*