

## Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 350-Labeled

Catalog Number: 46652

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 594 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 594 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 350 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

### Applications

---

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This configuration allows single-step fluorescence detection without requiring a secondary antibody. The antibody may cross-react with fluorophores similar to Alexa Fluor® 594, such as iFluor® 594, ATTO 594, and Texas Red®. Its blue emission is well-separated from orange-red dyes, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific.*