

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

Catalog Number: 46663

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 Volume 0.1 mL ddH2O

### 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 G beads. The antibody is conjugated with iFluor® 594 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 594
Excitation Wavelength	587 nm
Emission Wavelength	603 nm

### Applications

---

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 594-Labeled is a polyclonal antibody that selectively targets Alexa Fluor® 594-conjugated biomolecules. It is directly conjugated to iFluor® 594, a bright red-emitting fluorophore with excitation/emission maxima of ~590/617 nm, enabling robust one-step detection of Alexa Fluor® 594-labeled targets in immunofluorescence, microscopy, and flow cytometry.

This antibody may cross-react with spectrally similar dyes such as DyLight® 594, and ATTO 594. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

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