

## Rabbit Anti-Alexa Fluor 594 Antibody

Catalog Number: 46650

Unit Size: 1 mg

### 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, 0.05% Sodium Azide

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 594 coupled to carrier protein

### Biological Properties

---

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Alexa Fluor 594 coupled carrier protein in rabbit and affinity purified with Protein A beads.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Applications

---

Rabbit Anti-Alexa Fluor 594 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules. It is designed for detecting Alexa Fluor® 594-labeled proteins, antibodies, or probes. The antibody may cross-react with similar fluorophores such as iFluor® 594, ATTO 594, and Texas Red®. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection. *Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific.*