

## Rabbit Anti-Alexa Fluor 594 Antibody

Catalog Number: 46650

Unit Size: 1 mg

### Product Details

---

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

### Chemical Properties

---

Appearance	Liquid
Molecular Weight	N/A
Soluble In	PBS

### Spectral Properties

---

Excitation Wavelength	N/A
Emission Wavelength	N/A

### 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. ATTO® is a registered trademark of ATTO-TEC GmbH.*