

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

Catalog Number: 46652

Unit Size: 100 ug

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