

## Rabbit Anti-Alexa Fluor 555 Antibody, Biotinylated

Catalog Number: 46550

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	N/A
Emission Wavelength	N/A

### Applications

---

Rabbit Anti-Alexa Fluor 555 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-based systems in various assay formats.

The antibody may cross-react with related fluorophores such as iFluor® 555, ATTO 550, and Cy3. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

*Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.*