

## Rabbit Anti-Alexa Fluor 750 Antibody, Biotinylated

Catalog Number: 46610

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 750 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with biotin. This format enables flexible detection using streptavidin- or avidin-based systems across a wide range of assay types.

The antibody may cross-react with related near-infrared fluorophores such as iFluor® 750, ATTO® 740, and DyLight® 755. Its broad compatibility makes it well-suited for immunofluorescence, ELISA, blotting, and multiplexed detection workflows.

*Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.*