

## 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. Its broad assay compatibility makes it suitable for immunofluorescence, ELISA, blotting, and multiplexed applications.

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