

## Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 350-Labeled

Catalog Number: 46532

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 555 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 350, a UV-excitable blue fluorophore with excitation/emission maxima around 346/442 nm. This direct format enables one-step fluorescence detection of Alexa Fluor® 555 without the need for a secondary antibody.

The antibody may cross-react with related fluorophores such as iFluor® 555, ATTO® 550, and Cy3. Its spectral separation from Alexa Fluor® 555 makes it well suited for multicolor fluorescence applications including microscopy, flow cytometry, and ELISA.

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