

## Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 555-Labeled

Catalog Number: 46568

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	557 nm
-----------------------	--------

Emission Wavelength	570 nm
---------------------	--------

### Applications

---

Rabbit Anti-Alexa Fluor 647 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with iFluor® 555, a bright orange-red fluorophore with excitation/emission of ~555/565 nm. This direct format enables one-step detection of Alexa Fluor® 647-labeled targets without the need for secondary antibodies.

The antibody may cross-react with related fluorophores such as iFluor® 647, Cy5, ATTO® 647, and DyLight® 650. The orange-red fluorescence of iFluor® 555 is spectrally distinct from far-red dyes, supporting multiplex workflows in ELISA, flow cytometry, and fluorescence imaging.

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