

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

Catalog Number: 46633

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 350 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with iFluor® 350, a UV-excitable blue fluorophore with excitation around 350 nm and emission near 442 nm. This direct labeling enables one-step fluorescence detection, reducing assay complexity and minimizing background signal.

The antibody may cross-react with structurally similar dyes such as iFluor® 350, AMCA, and Cascade Blue®. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in microscopy, ELISA, and flow cytometry.

*Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.*