

## Rabbit Anti-Alexa Fluor 350 Antibody

Catalog Number: 46620

Unit Size: 1 mg

### 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 350 Antibody is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules. It is designed for detecting Alexa Fluor® 350-labeled antibodies, proteins, or other probes in fluorescence-based assays.

The antibody may cross-react with related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. It is suitable for use in fluorescence microscopy, ELISA, and other applications requiring blue-channel detection.

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