

# Rabbit Anti-Alexa Fluor 350 Antibody, Biotinylated

Catalog Number: 46640

Unit Size: 100 ug

## Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, BSA

## Unit Details

---

Reconstitution Volume	0.1 mL ddH <sub>2</sub> O
-----------------------	---------------------------

## Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 350 coupled to carrier protein

## Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 350 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

## Spectral Properties

---

Conjugate	Biotin
-----------	--------

## Applications

---

Rabbit Anti-Alexa Fluor 350 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 350-conjugated molecules and is directly labeled with biotin. This configuration enables flexible detection using streptavidin- or avidin-conjugated reporters in a variety of assay formats. The antibody may cross-react with structurally related blue fluorophores such as iFluor® 350, AMCA, and Cascade Blue®. Its versatility makes it ideal for signal amplification, conjugation

verification, and multiplexed fluorescence detection. *Alexa Fluor® and Cascade Blue® are registered trademarks of Thermo Fisher Scientific.*