

## Rabbit Anti-Coumarin Antibody, iFluor® 350-Labeled

Catalog Number: 46791

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, 0.2% (w/v) BSA

### Unit Details

---

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

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Coumarin coupled to carrier protein

### Biological Properties

---

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

### Spectral Properties

---

Conjugate	iFluor® 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

### Applications

---

Rabbit Anti-Coumarin Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to coumarin and coumarin-derived fluorophores and is directly labeled with iFluor® 350, a bright blue fluorophore excited at around 350 nm. This direct labeling format enables single-step fluorescence detection without secondary antibodies.

The antibody may cross-react with related blue-emitting dyes such as AMC, AMCA, and other coumarin analogs. The spectral match between the coumarin target and iFluor® 350 label enhances signal detection sensitivity, making it ideal for validating coumarin conjugation and for use in multiplexed fluorescence assays.