

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

Catalog Number: 46791

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-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.