

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, BSA

Unit Details

Reconstitution Volume	0.1 mL ddH2O
-----------------------	--------------

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 A 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.