

Rabbit Anti-Coumarin Antibody

Catalog Number: 46780

Unit Size: 1 mg

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, 0.05% Sodium Azide

Antibody Properties

Clonality	Polyclonal
Host	Rabbit
Immunogen	Coumarin coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Coumarin coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 350 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Rabbit Anti-Coumarin Antibody is a polyclonal antibody that specifically recognizes coumarin and coumarin-derived fluorophores commonly used as blue-emitting labels in biological assays. It provides high-affinity binding for detecting coumarin-conjugated biomolecules. The antibody may cross-react with related blue-emitting dyes such as AMC (7-amino-4-methylcoumarin), AMCA, and other coumarin derivatives. It is ideal for signal amplification in detecting low-abundance coumarin conjugates, validating dye incorporation, and enabling multiplexed fluorescence assays.