

Rabbit Anti-Coumarin Antibody, Biotinylated

Catalog Number: 46810

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 ICG 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	Biotinylated
-----------	--------------

Applications

Rabbit Anti-Coumarin Antibody, Biotinylated is a polyclonal antibody that specifically binds to coumarin and coumarin-derived fluorophores and is directly labeled with biotin. The biotinylation enables flexible detection using streptavidin- or avidin-based signal amplification systems across multiple assay platforms. The antibody may cross-react with structurally similar blue fluorophores such as AMC, AMCA, and other coumarin analogs. It is ideal for signal amplification in low-abundance analyte

detection, validating coumarin conjugation efficiency, and supporting multiplexed fluorescence workflows in ELISA, blotting, and microscopy.