

Rabbit Anti-Coumarin Antibody, Biotinylated

Catalog Number: 46810

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	N/A
Emission Wavelength	N/A

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.