

Rabbit Anti-Cy3 Antibody

Catalog Number: 46880

Unit Size: 1 mg

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-Cy3 Antibody is a polyclonal antibody that specifically binds to Cy3-conjugated biomolecules. It is designed to recognize the cyanine dye Cy3 commonly used for fluorescent labeling in microscopy, flow cytometry, and molecular biology assays.

The antibody may also cross-react with structurally similar fluorophores such as Alexa Fluor® 555, ATTO 550, and iFluor® 555. It is ideal for signal amplification in low-abundance target detection, verifying Cy3 conjugation efficiency, and enabling multiplex fluorescence analysis in various platforms.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.