

Rabbit Anti-Cy5 Antibody

Catalog Number: 46910

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-Cy5 Antibody is a polyclonal antibody that specifically binds to Cy5-labeled molecules. It is designed to detect Cy5-conjugated proteins, oligonucleotides, or other biomolecules in fluorescence-based assays with high sensitivity.

The antibody may cross-react with structurally similar dyes including iFluor® 647, Alexa Fluor® 647, DyLight® 650, and ATTO 647. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and supporting multiplex fluorescence workflows in microscopy, flow cytometry, and blotting platforms.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.