

Rabbit Anti-Cy7 Antibody

Catalog Number: 46940

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-Cy7 Antibody is a polyclonal antibody that specifically binds to molecules labeled with the near-infrared dye Cy7. It is designed for detecting Cy7-conjugated antibodies, proteins, or nucleic acids in fluorescence-based applications.

The antibody may cross-react with spectrally similar fluorophores such as Alexa Fluor® 750, DyLight® 755, and iFluor® 750. It is ideal for signal amplification in low-abundance target detection, verifying conjugation efficiency, and enabling multiplexed fluorescence assays.

Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.