

Rabbit Anti-Cy7 Antibody, Biotinylated

Catalog Number: 46960

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-Cy7 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy7-conjugated targets and is directly labeled with biotin. This biotinylated format enables flexible signal detection using streptavidin- or avidin-based systems across a range of assay platforms.

The antibody may cross-react with other near-infrared dyes 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 performing multiplex fluorescence assays.

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