

Rabbit Anti-Cy5 Antibody, Biotinylated

Catalog Number: 46930

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-Cy5 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy5-conjugated molecules and is directly labeled with biotin. This format enables flexible signal detection using streptavidin- or avidin-based reporters in both colorimetric and fluorescence-based assays.

The antibody may cross-react with structurally related dyes such as 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 multiplexed analysis across ELISA, blotting, and imaging workflows.

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