

Rabbit Anti-Cy3 Antibody, Biotinylated

Catalog Number: 46900

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

Chemical Properties

Appearance	Solid
------------	-------

Molecular Weight	~150000
------------------	---------

Soluble In	water
------------	-------

Spectral Properties

Excitation Wavelength	N/A
-----------------------	-----

Emission Wavelength	N/A
---------------------	-----

Applications

Rabbit Anti-Cy3 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Cy3-labeled targets and is directly conjugated to biotin. This biotinylated format enables flexible signal detection using streptavidin- or avidin-linked reporters across a wide range of fluorescence and enzymatic assays.

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

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