

Rabbit Anti-Cy3 Antibody, iFluor® 555-Labeled

Catalog Number: 46891

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	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-Cy3 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Cy3-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 550 nm. The direct labeling format enables one-step fluorescence detection, reducing background and improving workflow efficiency.

The antibody may cross-react with related dyes such as Alexa Fluor® 555, ATTO 550, and iFluor® 555. Its spectral alignment with Cy3 allows enhanced signal amplification without additional detection reagents, making it suitable for multiplex assays, microscopy, and flow cytometry.

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