

Rabbit Anti-Cy7 Antibody, iFluor® 750-Labeled

Catalog Number: 46951

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	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-Cy7 Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to Cy7-conjugated targets and is directly labeled with iFluor® 750, a near-infrared fluorophore with excitation/emission around 752/778 nm. This direct labeling enables sensitive, single-step detection of Cy7-labeled molecules without secondary antibodies.

The antibody may cross-react with similar dyes such as Alexa Fluor® 750, DyLight® 755, and other far-red fluorophores. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in deep-tissue imaging, flow cytometry, and high-content analysis.

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