

Rabbit Anti-Alexa Fluor 555 Antibody, iFluor® 555-Labeled

Catalog Number: 46543

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-Alexa Fluor 555 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 555-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation around 555 nm. This direct labeling format enables single-step detection, improving signal sensitivity and reducing background interference.

The antibody may cross-react with structurally similar dyes such as iFluor® 555, ATTO® 550, and Cy3. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation, and performing multiplex fluorescence assays.

Alexa Fluor® and Cy3 are registered trademarks of Thermo Fisher Scientific. ATTO® is a registered trademark of ATTO-TEC GmbH.