

Rabbit Anti-Alexa Fluor 750 Antibody, iFluor® 555-Labeled

Catalog Number: 46598

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 750 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission maxima of ~556/568 nm. This configuration allows for one-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with similar fluorophores such as iFluor® 750, ATTO® 740, and DyLight® 755. The spectral distinctness of iFluor® 555 makes it ideal for use in multiplex immunofluorescence, ELISA, and flow cytometry panels alongside far-red dyes.

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