

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 555-Labeled

Catalog Number: 46658

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 594 Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated targets and is directly labeled with iFluor® 555, an orange-emitting fluorophore with excitation/emission at around 556/570 nm. This direct format allows single-step fluorescence detection without the need for a secondary antibody.

The antibody may cross-react with structurally related dyes such as iFluor® 594, ATTO 594, and Texas Red®. It is ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® and Texas Red® are registered trademarks of Thermo Fisher Scientific.