

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 350-Labeled

Catalog Number: 46652

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	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated molecules and is directly labeled with iFluor® 350, a bright, UV-excitable blue fluorophore. This configuration allows single-step fluorescence detection without requiring a secondary antibody.

The antibody may cross-react with fluorophores similar to Alexa Fluor® 594, such as iFluor® 594, ATTO 594, and Texas Red®. Its blue emission is well-separated from orange-red dyes, making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

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