

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

Catalog Number: 46663

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	587 nm
Emission Wavelength	603 nm

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, iFluor® 594-Labeled is a polyclonal antibody that selectively targets Alexa Fluor® 594-conjugated biomolecules. It is directly conjugated to iFluor® 594, a bright red-emitting fluorophore with excitation/emission maxima of ~590/617 nm, enabling robust one-step detection of Alexa Fluor® 594-labeled targets in immunofluorescence, microscopy, and flow cytometry.

This antibody may cross-react with spectrally similar dyes such as DyLight® 594, and ATTO® 594. It is ideal for signal amplification in low-abundance target detection, verifying dye conjugation efficiency, and performing multiplex fluorescence assays in various applications.

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