

Rabbit Anti-Alexa Fluor 594 Antibody, Biotinylated

Catalog Number: 46670

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	N/A
Emission Wavelength	N/A

Applications

Rabbit Anti-Alexa Fluor 594 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 594-conjugated targets and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-linked systems in a wide range of assay formats.

The antibody may cross-react with structurally similar fluorophores 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.