

Rabbit Anti-Alexa Fluor 488 Antibody, Biotinylated

Catalog Number: 46520

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 488 Antibody, Biotinylated is a polyclonal antibody that specifically recognizes Alexa Fluor® 488-conjugated molecules and is directly conjugated to biotin. The biotin tag enables flexible signal amplification using streptavidin, or other avidin-based systems, making it suitable for diverse assay formats.

The antibody may cross-react with structurally related dyes such as iFluor® 488, ATTO 488, and rhodamine 110. Its broad binding and versatile detection makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.

Alexa Fluor® is a registered trademark of Thermo Fisher Scientific.