

Rabbit Anti-Alexa Fluor 750 Antibody, Biotinylated

Catalog Number: 46610

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 750 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with biotin. This format enables flexible detection using streptavidin- or avidin-based systems across a wide range of assay types.

The antibody may cross-react with related near-infrared fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. Its broad compatibility makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

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