

Rabbit Anti-Alexa Fluor 647 Antibody, Biotinylated

Catalog Number: 46580

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 647 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 647-conjugated molecules and is directly labeled with biotin. This configuration enables flexible signal detection using streptavidin- or avidin-based systems in a wide range of assay formats.

The antibody may cross-react with structurally similar fluorophores such as iFluor® 647, Cy5, ATTO® 647, and DyLight® 650. It is suitable for ELISA, Western blotting, immunofluorescence microscopy, and multiplexed detection applications.

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