

Rabbit Anti-DNP Antibody, Biotinylated

Catalog Number: 46870

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-DNP Antibody, Biotinylated is a polyclonal antibody that specifically recognizes DNP (dinitrophenyl) conjugates and is directly labeled with biotin. DNP is a common hapten used in immunoassay development. The biotin tag enables flexible detection using streptavidin- or avidin-based systems for a wide range of assay formats.

This antibody is suitable for detecting DNP-labeled haptens, peptides, or proteins making it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection.