

Rabbit Anti-DNP Antibody

Catalog Number: 46850

Unit Size: 1 mg

Product Details

Storage Conditions	Refrigerated (2-8 °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 is a polyclonal antibody that specifically binds to dinitrophenyl (DNP)-labeled molecules. It is ideal for detecting DNPTagged proteins, peptides, or other biomolecules in a range of biochemical assays including immunoassays and redox-related studies.

The antibody offers strong affinity for DNP epitopes and may cross-react with structurally related compounds. It is well-suited for applications such as protein labeling detection and multiplexed fluorescence assays.