

Mouse Monoclonal Anti-DNP Antibody, Biotinylated

Catalog Number: 46876

Unit Size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.2% (w/v) BSA

Unit Details

Reconstitution Volume	0.1 mL ddH ₂ O
-----------------------	---------------------------

Antibody Properties

Clonality	Monoclonal (IgG2A)
Host	Mouse
Immunogen	DNP

Biological Properties

Stabilizer	BSA
Appearance	Solid
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	Biotin
-----------	--------

Applications

Mouse Monoclonal Anti-DNP Antibody, Biotinylated is a monoclonal 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.

The monoclonal format ensures high specificity, low background, and lot-to-lot consistency, making it ideal for reproducible results. This antibody is suitable for detecting DNP-labeled haptens, peptides, or proteins, and is well suited for signal amplification, conjugation verification, and multiplexed fluorescence detection.