

Mouse Monoclonal Anti-DNP Antibody

Catalog Number: 46871

Unit Size: 1 mg

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, 0.05% Sodium Azide

Antibody Properties

Clonality	Monoclonal (IgG2A)
Host	Mouse
Immunogen	DNP

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Applications

Mouse Monoclonal Anti-DNP Antibody is a specific monoclonal antibody that recognizes dinitrophenyl (DNP) moieties. DNP is a widely used hapten for labeling proteins, nucleic acids, and other biomolecules in various immunoassay formats. This antibody is ideal for detecting DNP-conjugated molecules in ELISA, Western blot, immunocytochemistry, and other applications requiring hapten recognition.

The monoclonal format provides batch-to-batch consistency, superior specificity, and reduced background in complex biological samples. This antibody is purified and suitable for optimal performance in both direct and indirect immunodetection assays.