

Rabbit Anti-DNP Antibody

Catalog Number: 46850

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	Polyclonal
Host	Rabbit
Immunogen	DNP coupled to carrier protein

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Preparation	It is produced with Cy5 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

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.