

Rabbit Anti-DNP Antibody, Biotinylated

Catalog Number: 46870

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	Polyclonal
Host	Rabbit
Immunogen	DNP coupled to carrier protein

Biological Properties

Stabilizer	BSA
Appearance	Solid
Preparation	Antibody is produced with immunogen in rabbit, affinity purified and conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

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.