

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, Biotinylated

Catalog Number: 46985

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C)
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.2% (w/v) BSA, 0.05% Sodium Azide

Antibody Properties

Clonality	Monoclonal
Host	Mouse
Isotype	IgG1
Immunogen	DIG (Digoxigenin)

Biological Properties

Stabilizer	0.2% (w/v) BSA, 0.05% Sodium Azide
Appearance	Liquid
Soluble In	Water
Application	ELISA, ICC, WB, IHC

Spectral Properties

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

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, Biotinylated is engineered to recognize digoxigenin-conjugated molecules with high specificity. The biotin tag allows flexible detection using a wide range of streptavidin- or avidin-based systems, including HRP, AP, or fluorophore conjugates. This antibody is ideal for workflows such as ELISA, Western blotting, in situ hybridization, and fluorescence microscopy. Its monoclonal format ensures reproducibility across applications and compatibility with colorimetric, chemiluminescent, or fluorescent readouts.