

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody

Catalog Number: 46980

Unit Size: 1 mg

Product Details

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

Antibody Properties

Clonality	Monoclonal
Host	Mouse
Isotype	IgG1
Immunogen	DIG (Digoxigenin)

Biological Properties

Stabilizer	0.05% Sodium Azide
Appearance	Liquid
Soluble In	Water
Application	ELISA, ICC, IHC, WB

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody is designed to detect digoxigenin-conjugated molecules with high specificity and minimal cross-reactivity. Digoxigenin labeling is a widely used strategy for non-radioactive detection in both nucleic acid and protein-based assays. This antibody performs reliably in diverse applications such as ELISA, Western blotting, immunocytochemistry, and fluorescence or chromogenic in situ hybridization. Its monoclonal format supports reproducible results across batches and is compatible with various detection systems using enzyme-, fluorophore-, or biotin-labeled secondary antibodies.