

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, HRP-Conjugated

Catalog Number: 46986

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C)
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.01% thimerosal, 0.2% (w/v) BSA

Antibody Properties

Clonality	Monoclonal
Host	Mouse
Isotype	IgG1
Immunogen	DIG (Digoxigenin)

Biological Properties

Stabilizer	0.01% thimerosal, 0.2% (w/v) BSA
Appearance	Liquid
Soluble In	Water
Application	ELISA, WB, ICC, IHC

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, HRP-Conjugated is designed to detect digoxigenin-labeled 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 HRP-conjugated antibody enables direct enzymatic detection using substrates such as TMB, DAB, or ECL. It is suitable for a variety of applications, including ELISA, Western blotting, immunocytochemistry, and in situ hybridization workflows where DIG-labeled targets require sensitive and reliable signal development.