

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, Alkaline Phosphatase (AP)-Conjugated

Catalog Number: 46987

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, WB, ICC, IHC

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, Alkaline Phosphatase (AP)-Conjugated is designed for the sensitive detection of digoxigenin-labeled probes in enzyme-based assays. The antibody specifically binds DIG-conjugated targets and is coupled to alkaline phosphatase, allowing visualization via chromogenic or chemiluminescent substrates. This format is suitable for applications including ELISA, Western blotting, immunocytochemistry, and in situ hybridization. The monoclonal nature of the antibody ensures consistent performance and compatibility with DIG-based detection workflows requiring AP-conjugated reagents.