

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 350-Labeled

Catalog Number: 46981

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
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	IF, ELISA, FACS

Spectral Properties

Conjugate	iFluor 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 350-Labeled is designed for the fluorescent detection of digoxigenin-conjugated molecules using blue channel emission. The iFluor® 350 dye offers bright blue fluorescence when excited near 345 nm, supporting applications requiring minimal spectral interference with green and red channels. This antibody enables direct detection of DIG-labeled probes without the need for secondary antibodies, making it ideal for fluorescence in situ hybridization (FISH), immunocytochemistry, and multiplex imaging. Its monoclonal format ensures batch-to-batch consistency, and the conjugated fluorophore provides streamlined signal readout.