

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 750-Labeled

Catalog Number: 46984

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	ELISA, IF, FACS

Spectral Properties

Conjugate	iFluor 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 750-Labeled is designed to detect digoxigenin-conjugated molecules with high specificity and near-infrared fluorescence signal. iFluor® 750 exhibits excitation and emission maxima at ~757 nm and ~779 nm, enabling sensitive detection in the NIR channel for low-background imaging. The antibody's direct conjugation format simplifies protocols by eliminating the need for secondary antibodies. It is well-suited for fluorescence in situ hybridization (FISH), immunocytochemistry, and other NIR-based applications. The monoclonal format ensures reproducibility and consistency across experiments.