

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 488-Labeled

Catalog Number: 46982

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 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 488-Labeled is designed to detect digoxigenin-conjugated molecules with high specificity and green fluorescence readout. The iFluor® 488 dye exhibits excitation and emission maxima of 491 nm and 516 nm, making it compatible with standard green/FITC detection channels used in fluorescence microscopy and flow cytometry. This antibody enables direct fluorescent detection of DIG-labeled probes without the need for secondary reagents, making it well suited for fluorescence in situ hybridization (FISH), immunocytochemistry, and other multicolor fluorescence-based applications. Its monoclonal format supports consistent, reproducible results across experimental batches.