

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 555-Labeled

Catalog Number: 46983

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	FACS, ELISA, IF

Spectral Properties

Conjugate	iFluor 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

Applications

Mouse Monoclonal Anti-DIG (Digoxigenin) Antibody, iFluor® 555-Labeled is designed to detect digoxigenin-conjugated molecules with high specificity and orange fluorescence readout. The iFluor® 555 dye has excitation and emission maxima near 557 nm and 570 nm, providing bright signal output in the orange channel for fluorescence imaging applications. This directly labeled antibody eliminates the need for secondary detection steps and is ideal for fluorescence in situ hybridization (FISH), immunocytochemistry, and other fluorescence-based workflows. Its monoclonal format ensures consistent performance across batches and experiments.