

Mouse Monoclonal Anti-DNP Antibody, iFluor® 350-Labeled

Catalog Number: 46872

Unit Size: 100 ug

Product Details

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.2% (w/v) BSA

Unit Details

Reconstitution Volume	0.1 mL ddH ₂ O
-----------------------	---------------------------

Antibody Properties

Clonality	Monoclonal (IgG2A)
Host	Mouse
Immunogen	DNP

Biological Properties

Stabilizer	BSA
Appearance	Solid
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 350
Excitation Wavelength	345 nm
Emission Wavelength	450 nm

Applications

Mouse Monoclonal Anti-DNP Antibody, iFluor® 350-Labeled is a monoclonal antibody that specifically binds to dinitrophenyl (DNP) moieties and is directly labeled with iFluor® 350, a blue-emitting fluorophore excitable around 350 nm. This direct labeling format enables sensitive, single-step fluorescence detection of DNP-conjugated proteins, peptides, and small molecules.

The monoclonal antibody provides high specificity and lot-to-lot consistency, making it ideal for verifying DNP conjugation and detecting DNP-tagged molecules in fluorescence-based applications. Its blue fluorescence emission is well suited for multiplex immunofluorescence, flow cytometry, and other assays requiring minimal spectral overlap.