

## Mouse Monoclonal Anti-DNP Antibody, iFluor® 488-Labeled

Catalog Number: 46873

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 <sub>2</sub> 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® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

### Applications

---

Mouse Monoclonal Anti-DNP Antibody, iFluor® 488-Labeled is a monoclonal antibody that specifically binds to dinitrophenyl (DNP) moieties and is directly labeled with iFluor® 488, a green-emitting fluorophore excitable around 490 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 green fluorescence emission is well suited for multiplex immunofluorescence, flow cytometry, and other assays requiring minimal spectral overlap.