

## Rabbit Anti-DNP Antibody, iFluor® 488-Labeled

Catalog Number: 46855

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, BSA

### Unit Details

---

Reconstitution Volume	0.1 mL ddH <sub>2</sub> O
-----------------------	---------------------------

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	DNP coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Cy7 coupled carrier protein in rabbit and affinity purified with Protein A beads. The antibody is conjugated with iFluor® 750 under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 488
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

### Applications

---

Rabbit Anti-DNP Antibody, iFluor® 488-Labeled is a polyclonal antibody that specifically binds to dinitrophenyl (DNP) groups and is directly labeled with iFluor® 488, a green-emitting fluorophore excited at around 490 nm. This configuration allows direct, single-step fluorescence detection without the need for secondary reagents. The antibody is ideal for detecting DNP-labeled peptides, proteins, and other biomolecules with high sensitivity. It is suitable for validating DNP conjugation efficiency and enables multiplexed detection in fluorescence-based assays.