

## Rabbit Anti-DNP Antibody, iFluor® 555-Labeled

Catalog Number: 46858

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 ddH2O
-----------------------	--------------

### 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 Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	iFluor® 555
Excitation Wavelength	557 nm
Emission Wavelength	570 nm

### Applications

---

Rabbit Anti-DNP Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to dinitrophenyl (DNP) groups and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 550 nm. This configuration allows direct, single-step fluorescence detection of DNP-conjugated molecules, eliminating the need for secondary antibodies. The antibody is optimized for detecting DNP-labeled proteins, peptides, and probes in fluorescence-based applications. It is ideal for signal amplification in low-abundance targets, verifying conjugation efficiency, and supporting multiplexed imaging or assay formats.