

## Rabbit Anti-DNP Antibody, iFluor® 350-Labeled

Catalog Number: 46852

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.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

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

### Applications

---

Rabbit Anti-DNP Antibody, iFluor® 350-Labeled is a polyclonal 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 antibody is highly suitable for validating DNP conjugation and integrating DNP-tagged molecules into multiplex fluorescence-based assays.