

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, 0.2% (w/v) 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 G 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.