

## Mouse Monoclonal Anti-DNP Antibody, HRP-Conjugated

Catalog Number: 46877

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	Phosphate-buffered saline (PBS, pH 7.2), 0.01% thimerosal, 0.2% (w/v) BSA

### Unit Details

---

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

### Antibody Properties

---

Class	Secondary
Clonality	Monoclonal
Host	Mouse
Immunogen	DNP

### Biological Properties

---

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

### Spectral Properties

---

Conjugate	HRP
-----------	-----

### Applications

---

Mouse Monoclonal Anti-DNP Antibody, HRP-Conjugated is a monoclonal antibody that specifically recognizes DNP (dinitrophenyl) conjugates and is directly conjugated to horseradish peroxidase (HRP). DNP is a widely used hapten in immunoassay development. The HRP tag enables sensitive colorimetric or chemiluminescent detection in ELISA, Western blotting, and immunohistochemical assays. The monoclonal format ensures high specificity, low background, and reproducible performance across batches. This antibody is ideal for detecting DNP-labeled haptens, peptides, or proteins in enzymatic immunoassays requiring high sensitivity and reliable signal amplification.