

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, 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. 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.