

Rabbit Anti-DNP Antibody, iFluor® 750-Labeled

Catalog Number: 46860

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 pooled total human IgG, and affinity purified with human IgG coupled beads
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

Spectral Properties

Conjugate	iFluor® 750
Excitation Wavelength	757 nm
Emission Wavelength	779 nm

Applications

Rabbit Anti-DNP Antibody, iFluor® 750-Labeled is a polyclonal antibody that specifically binds to dinitrophenyl (DNP) groups and is directly labeled with iFluor® 750, a near-infrared fluorophore excited at around 750 nm. This direct labeling format

enables high-sensitivity detection of DNP-labeled molecules without the need for a secondary antibody.

The antibody is ideal for applications requiring minimal autofluorescence and deep-tissue imaging, multiplex immunofluorescence, and low-abundance target detection. It supports signal amplification workflows and verification of DNP conjugation in biological samples.