

Rabbit Anti-DNP Antibody, iFluor® 750-Labeled

Catalog Number: 46860

Unit Size: 100 ug

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

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.