

Rabbit Anti-DNP Antibody, iFluor® 555-Labeled

Catalog Number: 46858

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	557 nm
Emission Wavelength	570 nm

Applications

Rabbit Anti-DNP Antibody, iFluor® 555-Labeled is a polyclonal antibody that specifically binds to dinitrophenyl (DNP) groups and is directly labeled with iFluor® 555, an orange-emitting fluorophore excited at around 550 nm. This configuration allows direct, single-step fluorescence detection of DNP-conjugated molecules, eliminating the need for secondary antibodies.

The antibody is optimized for detecting DNP-labeled proteins, peptides, and probes in fluorescence-based applications. It is ideal for signal amplification in low-abundance targets, verifying conjugation efficiency, and supporting multiplexed imaging or assay formats.