

Rabbit Anti-DNP Antibody, iFluor® 488-Labeled

Catalog Number: 46855

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	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.