

Rabbit Anti-DNP Antibody, iFluor® 350-Labeled

Catalog Number: 46852

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	345 nm
Emission Wavelength	450 nm

Applications

Rabbit Anti-DNP Antibody, iFluor® 350-Labeled is a polyclonal antibody that specifically binds to dinitrophenyl (DNP) moieties and is directly labeled with iFluor® 350, a blue-emitting fluorophore excitable around 350 nm. This direct labeling format enables sensitive, single-step fluorescence detection of DNP-conjugated proteins, peptides, and small molecules.

The antibody is highly suitable for validating DNP conjugation and integrating DNP-tagged molecules into multiplex fluorescence-based assays.