



Glowing products for science

DNP Recombinant Monoclonal Rat Antibody (rLO-DNP-2) – Biotium Choice

A recombinant rat monoclonal antibody that recognizes the dinitrophenol (DNP) hapten for detection of DNP conjugated proteins and nucleic acids. The antibody belongs to the Biotium Choice list of select antibodies that have been validated in-house for optimal performance.



Product Description

DNP Recombinant Monoclonal Rat Antibody (rLO-DNP-2) is a recombinant rat monoclonal antibody that recognizes the dinitrophenol (DNP) hapten for detection of DNP conjugated proteins and nucleic acids. This antibody belongs to the Biotium Choice list of select antibodies that have been validated and optimized in-house for fluorescence microscopy. The antibody is available conjugated to a selection of CF® Dyes. They are supplied in PBS, 2 mg/mL rBSA, and 0.05% sodium azide.

- Available in 8 bright and photostable CF® Dyes, including [near-IR CF®740 and CF®750](#)
- Suitable for flow cytometry, immunohistochemistry, immunofluorescence, and western

Discover Sharper Signals and Unmatched Panel Flexibility with Biotium Choice Antibodies – Powered by CF® Dyes

Biotium Choice antibodies are carefully curated and validated in-house to offer exceptional signal-to-noise. Labeled with our advanced CF® Dyes, they are our top-recommended antibodies for immunofluorescence and other applications.

Biotium Choice Antibody Features

- Robust and validated clones against common targets
- Developed and optimized for immunofluorescence and other applications
- Conjugated to bright, photostable [CF® Dye](#) for superior signal and clarity
- New antibody clones and dye conjugates continuously in development

[View our full catalog of Biotium Choice antibodies](#)

DNP Recombinant Monoclonal Rat Antibody (rLO-DNP-2) – Biotium Choice

Conjugation	Ex/Em	Conc.	Size	Catalog No.	Dye Features
CF®488A	490/516 nm	100 ug/mL	1 mL	P038-488A-1ML	CF®488A Features
CF®568	562/584 nm	100 ug/mL	1 mL	P038-568-1ML	CF®568 Features
CF®594	593/615 nm	100 ug/mL	1 mL	P038-594-1ML	CF®594 Features
CF®640R	642/663 nm	100 ug/mL	1 mL	P038-640R-1ML	CF®640R Features
CF®647	652/668 nm	100 ug/mL	1 mL	P038-647-1ML	CF®647 Features
CF®680R	680/701 nm	100 ug/mL	1 mL	P038-680R-1ML	CF®680R Features
CF®740	742/767 nm	100 ug/mL	1 mL	P038-740-1ML	CF®740 Features
CF®750	755/779 nm	100 ug/mL	1 mL	P038-750-1ML	CF®750 Features

This datasheet was generated on December 24, 2025 at 05:26:36 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/dnp-recombinant-monoclonal-rat-antibody-rlo-dnp-2-biotium-choice/>

Call us : [800-304-5357](tel:800-304-5357)

Product attributes

Antibody Number	P038
Biotium Choice Antibody	Primary
Antibody type	Biotium Choice Primary
Host species	Rat
Antibody reactivity (target)	DNP
Isotype	IgG1, kappa
Synonyms	2,4-dinitrophenyl
Antibody application notes	Flow cytometry: 1 ug/million cells, Immunofluorescence: 1-5 ug/mL, Near-IR WB: 0.1-0.5 ug/mL. Optimal concentration to be determined by end-user.
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C, Protect fluorescent conjugates from light
Shelf life	Guaranteed for at least 24 months from date of receipt when stored as recommended
Regulatory status	For research use only (RUO)
Antibody/conjugate formulation	Conjugates: PBS/2 mg/mL rBSA/0.05% azide
Secondary/tag antibody applications	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
Antibody research areas	Protein tag
Product origin	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant rat IgG produced in Chinese Hamster Ovary (CHO) cell line