

## 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 and HRP.

- Available in bright and photostable CF® Dyes, including [near-IR CF@740](#), [CF@750](#), and [CF@790](#), and HRP
- 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
<a href="#">CF@488A</a>	490/516 nm	100 ug/mL	1 mL	<a href="#">P038-488A-1ML</a>	<a href="#">CF@488A Features</a>
<a href="#">CF@568</a>	562/584 nm	100 ug/mL	1 mL	<a href="#">P038-568-1ML</a>	<a href="#">CF@568 Features</a>
<a href="#">CF@594</a>	593/615 nm	100 ug/mL	1 mL	<a href="#">P038-594-1ML</a>	<a href="#">CF@594 Features</a>
<a href="#">CF@640R</a>	642/663 nm	100 ug/mL	1 mL	<a href="#">P038-640R-1ML</a>	<a href="#">CF@640R Features</a>
<a href="#">CF@647</a>	652/668 nm	100 ug/mL	1 mL	<a href="#">P038-647-1ML</a>	<a href="#">CF@647 Features</a>
<a href="#">CF@680R</a>	680/701 nm	100 ug/mL	1 mL	<a href="#">P038-680R-1ML</a>	<a href="#">CF@680R Features</a>
<a href="#">CF@740</a>	742/767 nm	100 ug/mL	1 mL	<a href="#">P038-740-1ML</a>	<a href="#">CF@740 Features</a>
<a href="#">CF@750</a>	755/779 nm	100 ug/mL	1 mL	<a href="#">P038-750-1ML</a>	<a href="#">CF@750 Features</a>
<a href="#">CF@790</a>	783/808 nm	100 ug/mL	1 mL	<a href="#">P038-790-1ML</a>	<a href="#">CF@790 Features</a>
<a href="#">HRP</a>	N/A	500 ug/mL	200 uL	<a href="#">P038-HRP-200UL</a>	

### Product attributes

<b>Antibody Number</b>	P038
<b>Biotium Choice Antibody</b>	Primary, Tag Antibody
<b>Antibody type</b>	Biotium Choice, Primary, Tag Antibody
<b>Host species</b>	Rat
<b>Antibody reactivity (target)</b>	DNP
<b>Isotype</b>	IgG1, kappa
<b>Synonyms</b>	2,4-dinitrophenyl
<b>Antibody application notes</b>	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.
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store CF® Dye conjugates at 2 °C to 8 °C, Store HRP conjugates at -10 °C to -35 °C, Protect from light
<b>Shelf life</b>	Guaranteed for at least 24 months from date of receipt when stored as recommended
<b>Regulatory status</b>	For research use only (RUO)
<b>Antibody/conjugate formulation</b>	CF® Dye Conjugates: PBS/2 mg/mL rBSA/ 0.05% sodium azide, HRP conjugates: PBS/50% glycerol/2 mg/mL rBSA
<b>Secondary/tag antibody applications</b>	Flow cytometry, IHC, IF (cells or tissue sections), Western blot
<b>Antibody research areas</b>	Protein tag
<b>Product origin</b>	Recombinant BSA produced in Chinese hamster ovary cells, Recombinant rat IgG produced in Chinese Hamster Ovary (CHO) cell line

This datasheet was generated on June 7, 2026 at 08:37:03 PM. 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/>