

# BrdU monoclonal antibody (2B1)

The thymidine analogue BrdU (5-Bromo-2-Deoxyuridine), a derivative of uridine, can substitute for thymidine during DNA synthesis. The detection of BrdU incorporation into DNA provides a common method to quantify newly synthesized DNA and identify cells in the S-phase of the cell cycle. The use of BrdU incorporation in proliferation assays facilitates the study of DNA repair, sister chromatid exchange, and the cytokinetics of normal and neoplastic cells.

This antibody is covered by our [Worry-Free Guarantee](#).

Citations: 1

[View Online »](#)

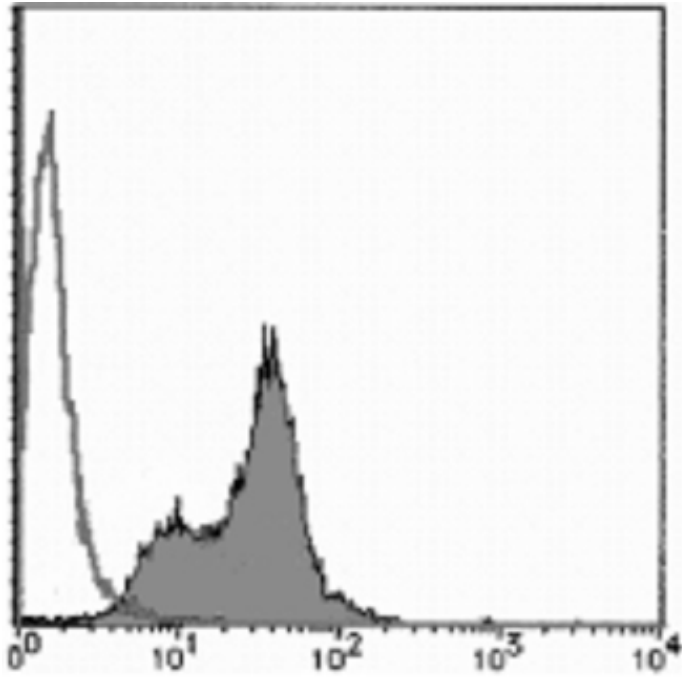
## Ordering Information

[Order Online »](#)

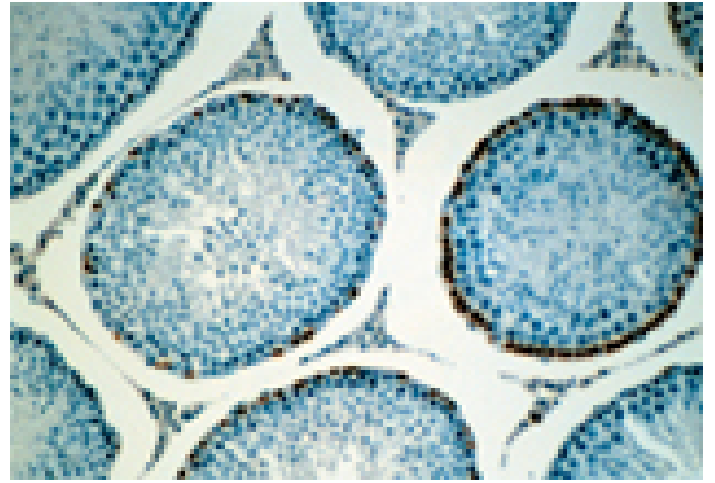
ADI-MSA-200-E	100µg
---------------	-------

## Manuals, SDS & CofA

[View Online »](#)



Flow cytometric analysis indicates the reaction of isotypic control (open histogram) and BrdU, mAb (2B1) (shaded histogram) to BrdU treated Raji cells.



Immunohistochemistry analysis of paraffin-embedded section of mouse testis stained with BrdU, mAb (2B1) to detect cells which have incorporated BrdU.

## Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

## Regulatory Status RUO - Research Use Only

## Product Details

Alternative Name	5-Bromodeoxyuridine
Application	Flow Cytometry, IHC
Clone	2B1
Formulation	Liquid. In PBS, pH 7.2, containing 50% glycerol.
Host	Mouse
Immunogen	Iododeoxyuridine-Ovalbumin.
Isotype	IgG1
Purity Detail	Protein A affinity purified.
Recommendation Dilutions/Conditions	Flow Cytometry (10µg/ml)Immunohistochemistry (10µg/ml)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined individually for each application.
Species Reactivity	Species independent
Worry-free Guarantee	This antibody is covered by our <a href="#">Worry-Free Guarantee</a>

Last modified: May 29, 2024



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)