

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 »](#)

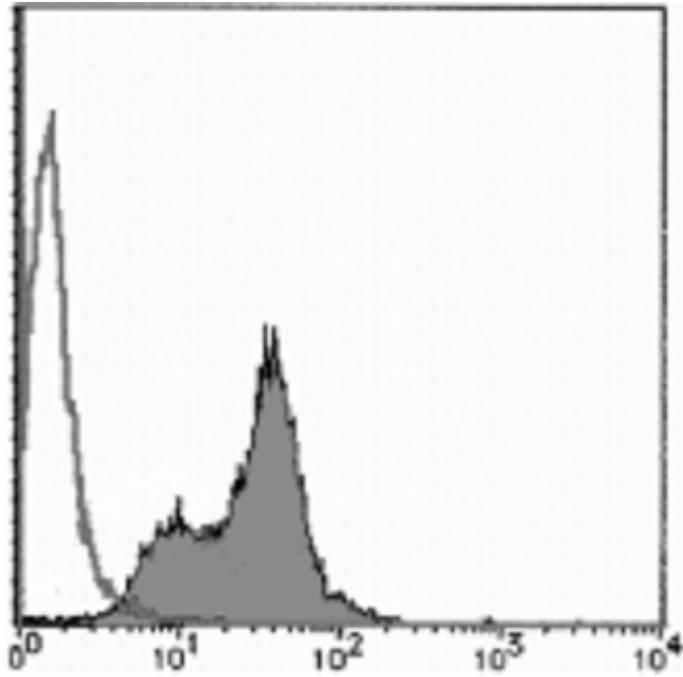
Ordering Information

[Order Online »](#)

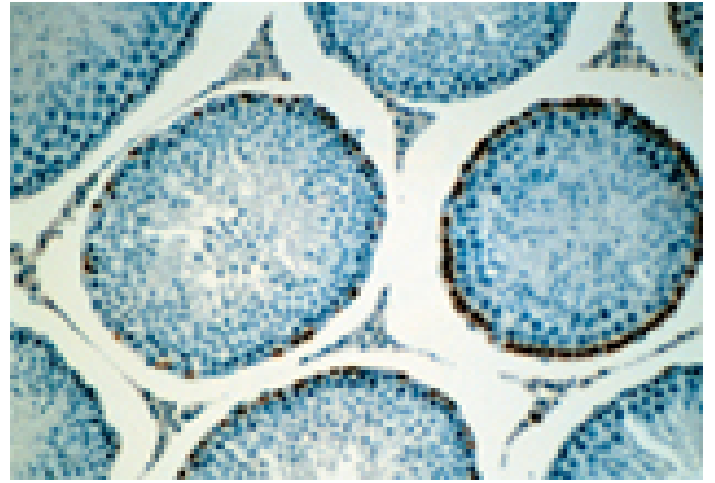
ADI-MSA-200-E	100µg
---------------	-------

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 Flow Cytometry (10µg/ml)Immunohistochemistry (10µg/ml)Suggested
Dilutions/Conditions 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 [Worry-Free Guarantee](#).



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-
fr@enzolifesciences.com](mailto:info-fr@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)