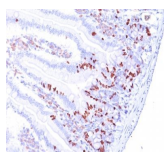


# BrdU Monoclonal Mouse Antibody (85-2C8)



## Product Description

Bromodeoxyuridine (BrdU) is a synthetic nucleoside thymidine analog. It is commonly used to detect proliferation in cells. In cells cultured in the presence of BrdU, BrdU will incorporate into newly synthesized DNA during S-phase. Antibodies specific for BrdU can then be used to detect replicated DNA.

This antibody is available purified, with BSA (0.2 mg/mL) or purified, BSA-free (1 mg/mL). **Catalog number key for antibody number 0886, Anti-BrdU (85-2C8)**

## Product attributes

<b>Antibody number</b>	#0886
<b>Antibody reactivity (target)</b>	BrdU
<b>Antibody type</b>	Primary
<b>Host species</b>	Mouse
<b>Clonality</b>	Monoclonal
<b>Clone</b>	85-2C8
<b>Isotype</b>	IgG1
<b>Molecular weight</b>	Depends on the target
<b>Synonyms</b>	Bromodeoxyuridine; 5-bromo-2-deoxyuridine
<b>Human gene symbol</b>	Not Applicable
<b>Entrez gene ID</b>	Not Applicable
<b>SwissProt</b>	Not Applicable
<b>Unigene</b>	Not Applicable
<b>Immunogen</b>	Bromodeoxyuridine (BrdU) conjugated to hemocyanine (isolated from <i>Helix pomatia</i> )
<b>Verified antibody applications</b>	IHC (FFPE) (verified)
<b>Antibody target cellular localization</b>	Nucleus
<b>Species reactivity</b>	All species
<b>Antibody application notes</b>	Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL
<b>Positive control</b>	Cells grown in presence of BrdU or tissues from experimental animals injected with BrdU
<b>Shipping condition</b>	Room temperature
<b>Storage Conditions</b>	Store at 2 to 8 °C, Protect fluorescent conjugates from light, Note: store BSA-free antibodies at -10 to -35 °C
<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>	Conjugates: 0.1 mg/mL in PBS/0.1% BSA/0.05% azide, HRP conjugates: 0.1 mg/mL in PBS/0.05% BSA, Purified: 0.2 mg/mL in PBS/0.05% BSA/0.05% azide, Purified, BSA-free: 1 mg/mL in PBS without azide
<b>Antibody research areas</b>	Cancer, Cell cycle
<b>Product origin</b>	Product may contain either bovine serum albumin (BSA) from bovine serum ( <i>Bos taurus</i> ), or recombinant BSA produced in Chinese hamster ovary cells. Inquire for the specific lot.

Antibody # prefix	Conjugation	Ex/Em (nm)	Laser line	Detection channel	Dye Features
BNC04	CF@405S	404/431	405	DAPI (microscopy), AF405	<a href="#">CF@405S Features</a>
BNC88	CF@488A	490/515	488	GFP, FITC	<a href="#">CF@488A Features</a>
BNC68	CF@568	562/583	532, 561	RFP, TRITC	<a href="#">CF@568 Features</a>
BNC94	CF@594	593/614	561	Texas Red®	<a href="#">CF@594 Features</a>
BNC40	CF@640R	642/662	633-640	Cy@5	<a href="#">CF@640R Features</a>
BNC47	CF@647	650/665	633-640	Cy@5	<a href="#">CF@647 Features</a>
BNC74	CF@740	742/767	633-685	775/50	<a href="#">CF@740 Features</a>
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified, BSA-free	N/A	N/A	N/A	

Alexa Fluor, Pacific Blue, Pacific Orange, and Texas Red are trademarks or registered trademarks of Thermo Fisher Scientific; Cy is a registered trademark of Cytiva; IRDye, LI-COR, and Odyssey are registered trademarks of LI-COR Bioscience.

This datasheet was generated on November 23, 2024 at 07:57:13 PM. Visit [product page](#) to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-brdu-bromodeoxyuridine-85-2c8/>