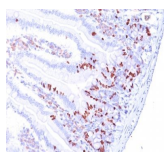


# BrdU Monoclonal Mouse Antibody (BRD469 + BRD494 + BRD.3)



## 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 1000, Anti-BrdU (BRD469 BRD494 BRD.3)**

## Product attributes

Antibody number	#1000
Antibody reactivity (target)	BrdU
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	BRD469 + BRD494 + BRD.3
Isotype	IgG1
Molecular weight	Depends on the target
Synonyms	Bromodeoxyuridine; 5-bromo-2-deoxyuridine
Human gene symbol	Not Applicable
Entrez gene ID	Not Applicable
SwissProt	Not Applicable
Unigene	Not Applicable
Immunogen	Bromodeoxyuridine (BrdU) conjugated to BSA or KLH
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	All species
Antibody application notes	Immunohistology formalin-fixed 0.5-1 ug/mL. Staining of formalin-fixed tissues, incubate sections in 4N HCl for 30 minutes at RT followed by digestion with trypsin at 1mg/mL PBS, 10 min at 37C. Immunofluorescence 0.5-1 ug/mL. Flow Cytometry 0.5-1 ug/million cells. Optimal dilution for a specific application should be determined by user
Positive control	Cells grown in presence of BrdU or tissues from experimental animals injected with BrdU
Shipping condition	Room temperature
Storage Conditions	Store at 2 to 8 °C. Protect fluorescent conjugates from light. Note: store BSA-free antibodies at -10 to -35 °C
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: 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
Antibody research areas	Cancer, Cell cycle
Product origin	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 February 13, 2025 at 04:57:53 AM. Visit product page to check for updated information before use. Product link: <https://biotium.com/product/monoclonal-anti-brdu-bromodeoxyuridine-brd469-brd494-brd-3/>