BrdU Monoclonal Mouse Antibody (BRD469)



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 0469, Anti-BrdU (BRD469)

Call us: 800-304-5357

Product attributes

Fibuuci ailiibules		
Antibody number	#0469	
Antibody reactivity (target)	BrdU	
Antibody type	Primary	
Host species	Mouse	
Clonality	Monoclonal	
Clone	BRD469	
Isotype	lgG1	
Molecular weight	307Da	
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 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, Immunofluorescence 0.5-1 ug/mL, Flow Cytometry 0.5-1 ug/million cells/0.1 mL	
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 (Bos taurus), 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	CF®405S Features
BNC88	CF®488A	490/515	488	GFP, FITC	CF®488A Features
BNC68	CF®568	562/583	532, 561	RFP, TRITC	CF®568 Features
BNC94	CF®594	593/614	561	Texas Red®	CF®594 Features
BNC40	CF®640R	642/662	633-640	Cy®5	CF®640R Features
BNC47	CF®647	650/665	633-640	Cy®5	CF®647 Features
BNC74	CF®740	742/767	633-685	775/50	CF®740 Features
BNCB	Biotin	N/A	N/A	N/A	
BNUB	Purified	N/A	N/A	N/A	
BNUM	Purified,	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 25, 2025 at 11:30:21 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/monoclonal-anti-brdu-bromodeoxyuridine-brd469/