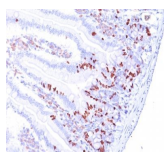


BrdU Monoclonal Mouse Antibody (MoBu-1)



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 0884, Anti-BrdU (MoBu-1)**

Product attributes

Antibody number	#0884
Antibody reactivity (target)	BrdU
Antibody type	Primary
Host species	Mouse
Clonality	Monoclonal
Clone	MoBu-1
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 hemocyanine (isolated from <i>Helix pomatia</i>)
Verified antibody applications	IHC (FFPE) (verified)
Antibody target cellular localization	Nucleus
Species reactivity	All species
Antibody application notes	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 (<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	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, 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 07:43:16 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/monoclonal-anti-brdu-bromodeoxyuridine-mobu-1/>