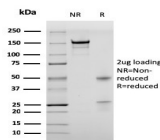


BrdU Recombinant Monoclonal Mouse Antibody (rBRD469)



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 1538, Anti-BrdU (rBRD469)**

Product attributes

| | |
|--|---|
| Antibody number | #1538 |
| Verified antibody applications | IHC (FFPE) (verified) |
| Antibody reactivity (target) | BrdU |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Monoclonal, Recombinant Monoclonal |
| Clone | rBRD469 |
| Isotype | IgG1, kappa |
| Molecular weight of antigen | Depends on the target |
| Synonyms | Bromodeoxyuridine; 5-bromo-2-deoxyuridine |
| Entrez gene ID | Not Applicable |
| SwissProt | Not Applicable |
| Unigene | Not Applicable |
| Immunogen | Bromodeoxyuridine (BrdU) conjugated to KLH |
| Antibody target cellular localization | Nucleus |
| Species reactivity | All species |
| Positive control | Cells grown in presence of BrdU or tissues from experimental animals injected with BrdU |
| Shipping condition | Room temperature |
| Storage Conditions | Note: store BSA-free antibodies at -10 to -35 °C, Store at 2 to 8 °C, Protect fluorescent conjugates from light |
| 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 |
| Shelf life | Guaranteed for at least 24 months from date of receipt when stored as recommended |
| 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 June 23, 2026 at 10:37:26 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/brdu-recombinant-monoclonal-mouse-antibody-rbrd469/>