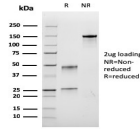


BrdU Recombinant Monoclonal Mouse Antibody (rBRD.3)



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).

Product attributes

| | |
|--|---|
| Antibody number | #3464 |
| Antibody reactivity (target) | BrdU |
| Antibody type | Primary |
| Host species | Mouse |
| Clonality | Recombinant Monoclonal |
| Clone | rBRD.3 |
| Isotype | IgG1, kappa |
| Molecular weight | 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 BSA |
| 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 |

This datasheet was generated on July 1, 2024 at 12:19:43 AM. Visit product page to check for updated information before use.
 Product link: <https://biotium.com/product/brdu-recombinant-monoclonal-mouse-antibody-rbrd-3/>