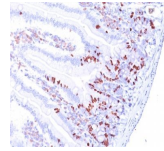


BrdU Monoclonal Mouse Antibody (BRD469 + BRD494 + BRD.3)



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).

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