

**Purified Mouse Anti-BrdU Antibody *Bu20a,
monoclonal***Catalog number: V103300
Unit size: 0.1 mg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	BrdU
Clone	Bu20a

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	FC (QC TESTED), IHC(P), IHC(F)

Applications

5-Bromo-2'-deoxyuridine (Bromodeoxyuridine, bromodeoxyuridine, BrdU, 5-bromodeoxyuridine, BUdR, BrDu, BUDR) is a pyrimidine 2'-deoxyribonucleoside compound having 5-bromouracil as the nucleobase. It has a variety of applications in the examination of key cell-based enzymes such as thymidylate kinase, sphingomyelin phosphodiesterase and thermostable purine nucleoside phosphorylase, especially in cell-lines like control fibroblasts, dividing cells and malignant cells. Harmonised basis, cell cycle and superpathway of pyrimidine deoxyribonucleotides de novo biosynthesis are, namely, just a few of the processes that it has helped researchers to better understand.