

Purified Mouse Anti-BrdU Antibody
MoBu-1, monoclonalCatalog number: V103295
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	MoBu-1

Biological Properties

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

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.