

BrdU antibody (mAb)

Catalog Nos: 61273, 61274

RRID: AB_2793579

Clone: Bu20a

Isotype: IgG1

Application(s): IHC

Reactivity: Not Species Specific

Quantities: 100 µg, 10 µg

Purification: Protein A Chromatography

Host: Mouse

Concentration: 1 µg/µl

Molecular Weight: Not Applicable

Background: BrdU (Bromodeoxyuridine) is a thymidine analogue, capable of being incorporated into DNA in place of thymidine during DNA synthesis. By pulse labelling with BrdU, the DNA synthesis in the cells can be analysed by detection with anti-BrdU antibodies. Antibodies against BrdU recognize BrdU in single-stranded DNA (produced by partial denaturation of double-stranded DNA), as well as free BrdU or BrdU coupled to a protein carrier.

Immunogen: This BrdU antibody was raised against Bromodeoxyuridine conjugated to BSA and recognizes BrdU.

Buffer: Purified IgG in PBS with 30% glycerol and 0.035% sodium azide. Sodium azide is highly toxic.

Application Notes:

This antibody is also available as an AbFlex[®] engineered recombinant antibody. For details on the corresponding AbFlex Recombinant Antibody, see Catalog No. 91269.

Storage and Guarantee: Some products may be shipped at room temperature. This will not affect their stability or performance. Avoid repeated freeze/thaw cycles by aliquoting items into single-use fractions for storage at -20°C for up to 2 years. Keep all reagents on ice when not in storage. This product is guaranteed for 12 months from date of receipt.

This product is for research use only and is not for use in diagnostic procedures.