

AbFlex® BrdU antibody (rAb)

Catalog Nos: 91269, 91270**RRID:** AB_2793827**Isotype:** IgG2a**Application(s):** IF**Reactivity:** Human, Not Species Specific**Quantities:** 100 µg, 10 µg**Purification:** Protein A Chromatography**Host:** Mouse**Concentration:** 1 µg/µl**Molecular Weight:** Not Applicable

Background: AbFlex® antibodies are recombinant antibodies (rAbs) that have been generated using defined DNA sequences to produce highly specific, reproducible antibodies. Each AbFlex antibody contains a 6xHis Tag, a Biotinylation Tag for enzymatic biotin conjugation using the biotin ligase, BirA, and a sortase recognition motif (LPXTG) to attach a variety of labels directly to the antibody including fluorophores, enzymatic substrates (HRP, AP), peptides, drugs as well as solid supports. AbFlex® BrdU antibody was expressed as full-length IgG with mouse immunoglobulin heavy and light chains (IgG2a isotype) in mammalian 293 cells.

BrdU (Bromodeoxyuridine) is a thymidine analogue, capable of being incorporated into DNA in place of thymidine during DNA synthesis. By pulse labeling with BrdU, the DNA synthesis in the cells can be analyzed 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 antibody was raised against Bromodeoxyuridine conjugated to BSA (Bovine Serum Albumin).

Buffer: Purified IgG in 140 mM Hepes, pH 7.5, 70 mM NaCl, 32 mM NaOAc, 0.035% sodium azide, 30% glycerol. Sodium azide is highly toxic.

Application Notes:

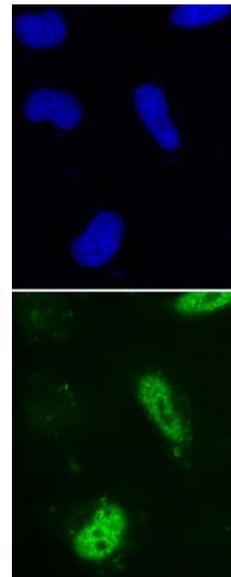
Applications Validated by Active Motif:

IF: 0.5 - 2 µg/ml

AbFlex® recombinant antibodies are genetically derived from DNA sequences of parental hybridoma clones. For details on the parental clone, see Catalog No. 61273.

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.



AbFlex® BrdU antibody (rAb) tested by Immunofluorescence
HeLa cells were stained with 0.5 µg/ml BrdU antibody, followed by anti-mouse IgG2a – Alexa Fluor 488 (lower panel) or DAPI (upper panel). Prior to staining, HeLa cells were treated for 1 hr at 37°C in cell culture medium supplemented with 10 µM BrdU. DNA was denatured with NaOH treatment for detection of BrdU incorporation in cells undergoing DNA replication at the time of labeling.