

BrdU (5-Bromo-2'-deoxyuridine)

BrdU (5-Bromo-2'-deoxyuridine) can be incorporated into DNA during cell division and subsequently detected by a BrdU antibody. The probe can be used to study cell-cycle kinetics (1).



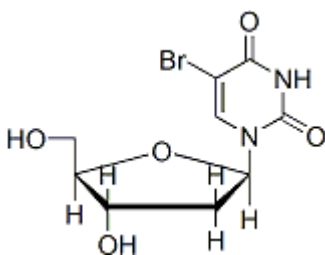
Product attributes

CAS number	59-14-3
------------	---------

Product Description

BrdU (5-Bromo-2'-deoxyuridine) can be incorporated into DNA during cell division and subsequently detected by a BrdU antibody (see our [anti-BrdU antibodies](#)). The probe can be used to study cell-cycle kinetics (1).

- White solid soluble in DMSO
- Store at -20 °C
- C₉H₁₁BrN₂O₅
- MW: 307.10
- [59-14-3]



References

1. Methods Cell Biol 41, 297 (1994).

This datasheet was generated on June 20, 2026 at 07:15:52 PM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/5-bromo-2-deoxyuridine-brdu/>