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

Call us: 800-304-5357

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 <u>anti-BrdU antibodies</u>). 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 November 25, 2025 at 11:21:34 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/5-bromo-2-deoxyuridine-brdu/