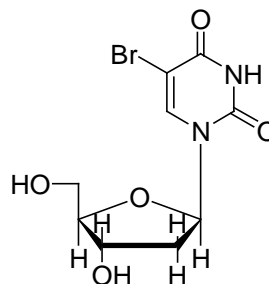


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: 5-Bromo-2'-deoxyuridine**CATALOG #** 40024**MOLECULAR INFORMATION:** C₉H₁₁BrN₂O₅
MWt: 307.10
[59-14-3]**PROPERTIES:**

Color & Form	Off-white solid
Solubility	Soluble in DMSO
Purity	≥ 95% by HPLC
Absorption	280nm (pH1.0)
Extinction Coefficient	9,000

STORAGE AND HANDLING: Store at 4°C**APPLICATION:** BrdU 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.¹Ref.: *Methods Cell Biol* **41**, 297(1994)**TOXICITY:** Not established. Not listed by NTP, IARC or OSHA.

FIRST AID:	Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing, move individual to fresh air and seek medical advice immediately.
-------------------	---

Disclaimer: <i>Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.</i>
--