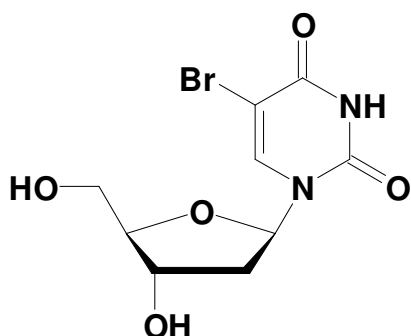


5-Br-dU

5-Bromo-2'-deoxy-uridine

Cat. No.	Amount
N-1040	100 mg



Cat. No.: N-1040

Molecular Formula: C₉H₁₁N₂O₅Br

Molecular Weight: 307.10

Purity: > 99%

Storage conditions:

Short term exposure (up to 1 week cumulative) to ambient temperature possible. Long term storage at < -20°C. If stored as recommended, Jena Bioscience guarantees optimal performance of this product for 12 months after date of delivery.

For research use only!

Used for detection of proliferating cells in living tissues.

BrdU is efficiently incorporated into DNA of replicating cells by substituting thymidine. BrdU containing DNA is subsequently detected with BrdU-specific antibodies.

Selected References:

Gratzner *et al.* (1992) Monoclonal antibody to 5-bromo and 5-iododeoxyuridine: a new reagent for detection of DNA replication. *Science* **218** :474.

Qin *et al.* (1993) Comparison of the classical autoradiographic and the immunohistochemical methods with BrdU for measuring proliferation parameters in colon cancer. *Anticancer Res* **13** :731.

Morstyn *et al.* (1986) immunohistochemical identification of proliferating cells in organ culture using bromodeoxyuridine and a monoclonal antibody. *J Histochem Cytochem* **34 (6)** :697.