

Product Information

BrdU (5-Bromo-2'-Deoxyuridine)

Catalog No.	Unit Size
A011	50 mg

Storage upon receipt:

- -20°C
- Protect from light

Product Description

Cells can naturally incorporate the thymidine analog **BrdU** into their cells during cell division, making this nucleoside analog an excellent marker for cell cycle and proliferation analysis. Analysis of incorporated BrdU can be with an anti-BrdU antibody.

Chemical structures

5-bromo-2'-deoxyuridine

Molecular formula: C₉H₁₁BrN₂O₅

Molecular weight: 307.1

CAS number: 59-14-3

