PRODUCT INFORMATION



5-Bromouridine

Item No. 30457

CAS Registry No.:	957-75-5	
Formal Name:	5-bromo-uridine	OH
Synonyms:	(-)-5-Bromouridine, BrU, BrUrd,	'
	NSC 38296	0 0 0
MF:	C ₉ H ₁₁ BrN ₂ O ₆	н Ц н
FW:	323.1	N N
Purity:	≥98%	ОН
UV/Vis.:	λ _{max} : 211, 280 nm	0
Supplied as:	A solid	
Storage:	-20°C	Br
Stability:	≥4 years	

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

5-Bromouridine is supplied as a solid. A stock solution may be made by dissolving the 5-bromouridine in the solvent of choice, which should be purged with an inert gas. 5-Bromouridine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of 5-bromouridine in these solvents is approximately 15 and 10 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of 5-bromouridine can be prepared by directly dissolving the solid in aqueous buffers. The solubility of 5-bromouridine in PBS, pH 7.2, is approximately 5 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

5-Bromouridine is a brominated analog of the nucleoside uridine (Item No. 20300).¹ It can be incorporated into RNA and subsequently detected by antibodies against bromodeoxyuridine (Item No. 15580).² 5-Bromouridine decreases the viability of HL-60 and MOLT-4 cells (LC₅₀s = 10 and 20 μ M, respectively). It induces apoptosis and halts the cell cycle at the S phase in HL-60 cells. It is photoreactive, and UV irradiation has been used to cross-link RNA containing 5-bromouridine to proteins in the study of of RNA-protein interactions.^{1,3} 5-Bromouridine can also be incorporated into RNA using 5-bromouridine 5'-triphosphate (Item No. 18140).

References

- 1. Tanner, N.K., Hanna, M.M., and Abelson, J. Binding interactions between yeast tRNA ligase and a precursor transfer ribonucleic acid containing two photoreactive uridine analogues. Biochemistry 27(24), 8852-8861 (1988).
- 2. Li, X., Patel, R., Melamed, M.R., et al. The cell cycle effects and induction of apoptosis by 5-bromouridine in cultures of human leukaemic MOLT-4 and HL-60 cell lines and mitogen-stimulated normal lymphocytes. Cell Prolif. 27(6), 307-319 (1994).
- Gott, J.M., Willis, M.C., Koch, T.H., et al. A specific, UV-induced RNA-protein cross-link using 3. 5-bromouridine-substituted RNA. Biochemistry 30(25), 6290-6295 (1991).

WARNING THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFFTY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 10/10/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD ANN ARBOR, MI 48108 · USA PHONE: [800] 364-9897 [734] 971-3335 FAX: [734] 971-3640 CUSTSERV@CAYMANCHEM.COM WWW.CAYMANCHEM.COM