

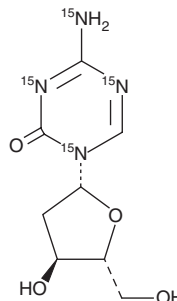
PRODUCT INFORMATION



Decitabine-¹⁵N₄

Item No. 22091

Formal Name: 4-(amino-¹⁵N)-1-((2R,4S,5R)-4-hydroxy-5-(hydroxymethyl)tetrahydrofuran-2-yl)-1,3,5-triazin-2(1H)-one-1,3,5-¹⁵N₃
Synonyms: DAC-¹⁵N₄, 5-aza-2'-Deoxycytidine-¹⁵N₄
MF: C₈H₁₂[¹⁵N]₄O₄
FW: 232.2
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Decitabine-¹⁵N₄ is supplied as a solid. A stock solution may be made by dissolving the decitabine-¹⁵N₄ in the solvent of choice. Decitabine-¹⁵N₄ is slightly soluble in the organic solvents DMSO and methanol (warmed).

Description

Decitabine-¹⁵N₄ is intended for use as an internal standard for the quantification of decitabine (Item No. 11166). Decitabine is a 2'-deoxy analog of 5-azacytidine (Item No. 11164) that induces dose-dependent inhibition of DNA methyltransferases and hypomethylation of DNA.¹⁻³ This action is useful in conditions characterized by DNA hypermethylation, as is found in myelodysplastic syndromes.⁴

References

1. Esteller, M. Epigenetics in cancer. *N. Engl. J. Med.* **358**, 1148-1159 (2008).
2. Lyko, F. and Brown, R. DNA methyltransferase inhibitors and the development of epigenetic cancer therapies. *J. Natl. Cancer Inst.* **97(20)**, 1498-1506 (2005).
3. Stresemann, C., Brueckner, B., Musch, T., *et al.* Functional diversity of DNA methyltransferase inhibitors in human cancer cell lines. *Cancer Res.* **66(5)**, 2794-2800 (2006).
4. Ornstein, M.C. and Sekeres, M.A. Combination strategies in myelodysplastic syndromes. *Int. J. Hematol.* **95**, 26-33 (2012).

WARNING

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

SAFETY 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, 12/22/2022

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