

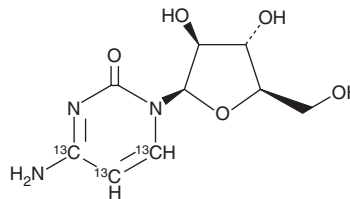
PRODUCT INFORMATION



Cytarabine-¹³C₃

Item No. 28805

Formal Name:	4-amino-1-β-D-arabinofuranosyl-2(1H)-pyrimidinone-4,5,6- ¹³ C ₃
Synonyms:	Ara-C- ¹³ C ₃ , 1-β-D-Arabinofuranosylcytosine- ¹³ C ₃
MF:	C ₆ [¹³ C ₃]H ₁₃ N ₃ O ₅
FW:	246.2
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

Cytarabine-¹³C₃ (ara-C-¹³C₃) is supplied as a solid. A stock solution may be made by dissolving the ara-C-¹³C₃ in the solvent of choice, which should be purged with an inert gas. Ara-C-¹³C₃ is soluble in the organic solvent DMSO (warmed).

Ara-C-¹³C₃ is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Ara-C-¹³C₃ is intended for use as an internal standard for the quantification of ara-C (Item No. 16069) by GC- or LC-MS. Ara-C is a nucleoside analog and prodrug form of the DNA polymerase inhibitor ara-CTP.¹ It is triphosphorylated to ara-CTP by the successive actions of deoxycytidine kinase, deoxycytidylate kinase, and nucleoside diphosphate kinase.² Ara-C inhibits proliferation of HL-60, ML-1, Raji, and Jurkat human leukemia cell lines with IC₅₀ values of 37, 17, 16, and 72 nM, respectively.³ It induces cell cycle arrest at the G₀/G₁ phase in HL-60 cells when used at concentrations of 2.5 and 15 μM.¹ Ara-C (75 mg/kg per day, i.p.) reduces tumor growth and increases tumor caspase-3 activity in an MOLM-13 mouse xenograft model.⁴ It also increases survival and reduces brain herpesvirus titers in infected rats when administered subcutaneously at doses of 80 and 320 mg/kg.⁵ Formulations containing ara-C have been used in the treatment of acute myeloid leukemia.

References

1. Li, Z., Guo, J.-R., Chen, Q.-Q., *et al.* Exploring the antitumor mechanism of high-dose cytarabine through the metabolic perturbations of ribonucleotide and deoxyribonucleotide in human promyelocytic leukemia HL-60 cells. *Molecules* **22**(3), E499 (2017).
2. Emadi, A. and Karp, J.E. The clinically relevant pharmacogenomic changes in acute myelogenous leukemia. *Pharmacogenomics* **13**(11), 1257-1269 (2012).
3. Qin, T., Youssef, E.M., Jelinek, J., *et al.* Effect of cytarabine and decitabine in combination in human leukemic cell lines. *Clin. Cancer Res.* **13**(14), 4225-4232 (2007).
4. Kelly, K.R., Espitia, C.M., Taverna, P., *et al.* Targeting PIM kinase activity significantly augments the efficacy of cytarabine. *Br. J. Haematol.* **156**(1), 129-132 (2012).
5. Renis, H.E. Antiviral activity of cytarabine in Herpesvirus-infected rats. *Antimicrob. Agents Chemother.* **4**(4), 439-444 (1973).

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, 12/09/2022

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