

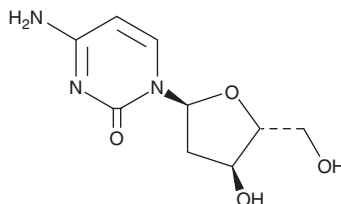
PRODUCT INFORMATION



α -2'-Deoxycytidine

Item No. 46222

CAS Registry No.: 4449-40-5
Formal Name: 4-amino-1-(2-deoxy- α -D-erythro-pentofuranosyl)-2(1H)-pyrimidinone
Synonym: 2'-Deoxy- α -Cytidine
MF: C₉H₁₃N₃O₄
FW: 227.2
Purity: $\geq 95\%$
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

α -2'-Deoxycytidine is supplied as a solid. Aqueous solutions of α -2'-deoxycytidine can be prepared by directly dissolving the solid in aqueous buffers. α -2'-Deoxycytidine is soluble (≥ 10 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

α -2'-Deoxycytidine is the unnatural α -anomer of 2'-deoxycytidine (Item Nos. 34708 | 30125). It inhibits the growth of L5178Y leukemia cells ($EC_{50} = 50$ μ g/ml).¹

Reference

1. Yamaguchi, T. and Saneyoshi, M. Synthetic nucleosides and nucleotides. XXI. On the synthesis and biological evaluations of 2'-deoxy- α -D-ribofuranosyl nucleosides and nucleotides. *Chem. Pharm. Bull. (Tokyo)* **32**(4), 1441-1450 (1984).

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, 08/24/2026

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