

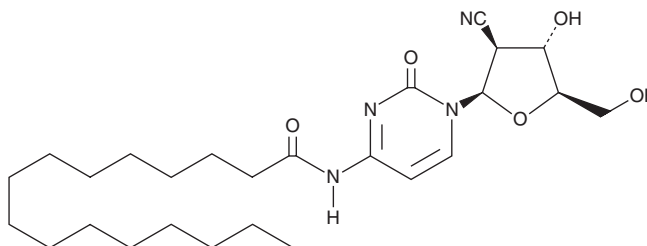
PRODUCT INFORMATION



Sapacitabine

Item No. 27321

CAS Registry No.: 151823-14-2
Formal Name: N-[1-(2-cyano-2-deoxy-b-D-arabinofuranosyl)-1,2-dihydro-2-oxo-4-pyrimidinyl]-hexadecanamide
Synonyms: CS 682, CYC-682
MF: C₂₆H₄₂N₄O₅
FW: 490.6
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

Sapacitabine is supplied as a solid. A stock solution may be made by dissolving the sapacitabine in the solvent of choice, which should be purged with an inert gas. Sapacitabine is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Sapacitabine is a prodrug form of the nucleoside derivative CNDAC (Item No. 40672).¹ *In vivo*, sapacitabine (5 and 15 mg/kg) reduces tumor volume in an MV4-11 mouse xenograft model when administered alone or in combination with the histone deacetylase (HDAC) inhibitor SAHA (vorinostat; Item No. 10009929).

Reference

- Green, S.R., Choudhary, A.K., and Fleming, I.N. Combination of sapacitabine and HDAC inhibitors stimulates cell death in AML and other tumour types. *Br. J. Cancer* **103(9)**, 1391-1399 (2010).

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, 05/30/2024

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