

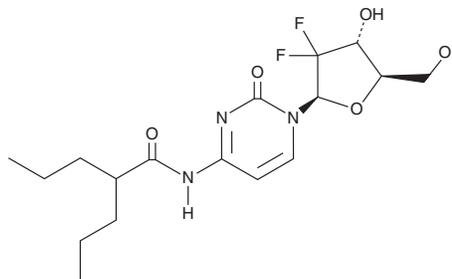
PRODUCT INFORMATION



LY2334737

Item No. 37425

CAS Registry No.: 892128-60-8
Formal Name: 2'-deoxy-2',2'-difluoro-N-(1-oxo-2-propylpentyl)-cytidine
MF: C₁₇H₂₅F₂N₃O₅
FW: 389.4
Purity: ≥95%
UV/Vis.: λ_{max}: 211, 252, 302 nm
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

LY2334737 is supplied as a solid. A stock solution may be made by dissolving the LY2334737 in the solvent of choice, which should be purged with an inert gas. LY2334737 is soluble in DMSO.

Description

LY2334737 is a prodrug form of the nucleoside analog gemcitabine (Item Nos. 11690 | 9003096).¹ *In vivo*, LY2334737 (8 mg/kg) reduces tumor volume in an MDA321-LM2-4175 breast cancer mouse xenograft model and increases survival in an SKOV3 ovarian cancer mouse xenograft model.

Reference

1. Francia, G., Shaked, Y., Hashimoto, K., *et al.* Low-dose metronomic oral dosing of a prodrug of gemcitabine (LY2334737) causes antitumor effects in the absence of inhibition of systemic vasculogenesis. *Mol. Cancer Ther.* **11**(3), 680-689 (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, 10/04/2023

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