

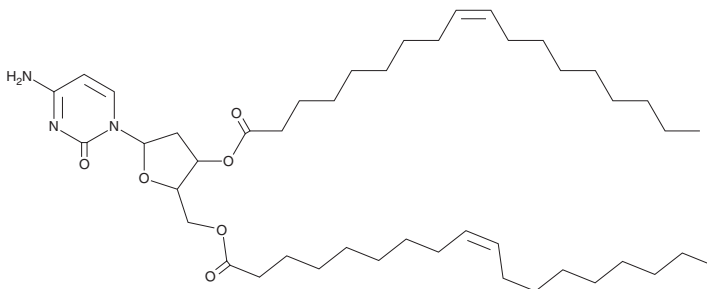
PRODUCT INFORMATION



3',5'-diOA-dC

Item No. 39087

Formal Name: 5-(4-amino-2-oxopyrimidin-1(2H)-yl)-2-((oleoyloxy)methyl)tetrahydrofuran-3-yl oleate
Synonym: 3',5'-Dioleic Acid-deoxycytidine
MF: C₄₅H₇₇N₃O₆
FW: 756.1
Purity: ≥95%
UV/Vis.: λ_{max}: 244, 273 nm
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥1 year



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

Description

3',5'-diOA-dC is a ligand for the bacterial second messenger, cyclic dinucleotide, and STING agonist cyclic di-GMP (Item No. 17144) composed of oleic acid (Item Nos. 90260 | 24659) and deoxycytidine.¹ It assembles with cyclic di-GMP to form supramolecular nanoparticles and has been used to deliver cyclic di-GMP *in vitro* and *in vivo*. Intravenous administration of nanoparticles composed of 3',5'-diOA-dC and cyclic di-GMP decrease tumor weight and volume, increase the number of CD8⁺ T cells, neutrophils, and natural killer (NK) cells in the tumor microenvironment, and increase tumoral levels of TNF-α and IFN-γ in a B16/F10 murine melanoma model.

Reference

1. Xu, L., Deng, H., Wu, L., *et al.* Supramolecular cyclic dinucleotide nanoparticles for STING-mediated cancer immunotherapy. *ACS Nano* **17**(11), 10090-10103 (2023).

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, 10/12/2023

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