

PRODUCT INFORMATION



G0-C14 analog

Item No. 40194

CAS Registry No.: 2935435-86-0

Formal Name: 3,3',3'',3'''-(ethane-1,2-diylbis(azanetriyl))tetrakis (N-(2-(bis(2-hydroxytetradecyl) amino)ethyl)propanamide)

MF: $C_{134}H_{272}N_{10}O_{12}$

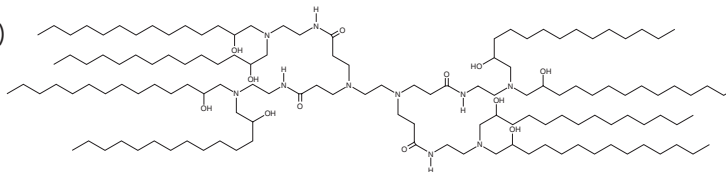
FW: 2,215.7

Purity: $\geq 95\%$ (mixture of isomers)

Supplied as: A 5 mg/ml solution in ethanol

Storage: -20°C

Stability: ≥ 3 years



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

Laboratory Procedures

G0-C14 analog is supplied as a solution in ethanol. To change the solvent, simply evaporate the ethanol under a gentle stream of nitrogen and immediately add the solvent of choice. G0-C14 analog is soluble in chloroform.

Description

G0-C14 analog is a derivative of the ionizable cationic lipid G0-C14.^{1,2}

References

1. Xiao, Y., Duda, G.D., and Shi, J. Lipid-based compositions and methods thereof. *The Brigham and Women's Hospital and The General Hospital Corporation*. **WO2023/107574A2** (2023).
2. Kong, N., Tao, W., Ling, X., et al. Synthetic mRNA nanoparticle-mediated restoration of p53 tumor suppressor sensitizes p53-deficient cancers to mTOR inhibition. *Sci. Transl. Med.* **11(523)**, eaaw1565 (2019).

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, 11/18/2025

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