

CERTIFICATE of ANALYSIS



G0-C14 analog

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

Item No. 40194 • Batch No. 0696114

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

Molecular Formula.: C₁₃₄H₂₇₂N₁₀O₁₂

CAS Number: 2935435-86-0

Formula Weight: 2215.65368

Expiry date: 08DEC2026

Overview

Tests	Results
HPLC	Purity: 100.0 %
Mass spec	[M+2H] ²⁺ : 1109.03 [M+3H] ³⁺ : 739.16
NMR	Conforms

Reviewed and approved by: Libby Perry

WARNING

THIS PRODUCT IS FOR RESEARCH USE - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE. IT IS THE RESPONSIBILITY OF THE PURCHASER TO DETERMINE SUITABILITY FOR OTHER APPLICATIONS.

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, 10/12/2018

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