

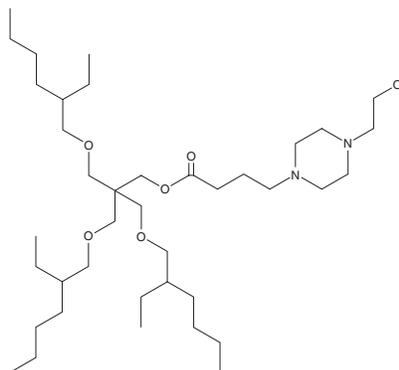
PRODUCT INFORMATION



IAJD93

Item No. 39688

CAS Registry No.: 3076338-50-3
Formal Name: 3-((2-ethylhexyl)oxy)-2,2-bis(((2-ethylhexyl)oxy)methyl)propyl 4-(4-(2-hydroxyethyl)piperazin-1-yl)butanoate
Synonym: Ionizable Amphiphilic Janus Dendrimer 93
MF: C₃₉H₇₈N₂O₆
FW: 671.0
Purity: ≥95%
Supplied as: A 25 mg/ml solution in ethanol
Storage: -20°C
Stability: ≥2 years
Special Conditions: Oxygen sensitive



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

Laboratory Procedures

IAJD93 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. Solvents such as methanol and chloroform purged with an inert gas can be used.

Description

IAJD93 is an ionizable amphiphilic Janus dendrimer (IAJD) delivery system for mRNA ($pK_a = 6.25$).¹ It co-assembles with mRNA to form dendrimer-some nanoparticles (DNPs) in acetate buffer. IAJD93 co-assembled with luciferase mRNA delivers the mRNA to the spleen, liver, lungs, and lymph nodes in mice.

Reference

1. Lu, J., Atochina-Vasserman, E.N., Maurya, D.S., *et al.* Targeted and equally distributed delivery of mRNA to organs with pentaerythritol-based one-component ionizable amphiphilic Janus dendrimers. *J. Am. Chem. Soc.* **145**(34), 18760-18766 (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, 01/27/2026

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