

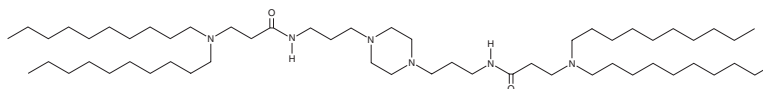
PRODUCT INFORMATION



PPZ-A10

Item No. 9004144

CAS Registry No.: 2941268-67-1
Formal Name: N,N'-(piperazine-1,4-diylbis(propane-3,1-diyl))bis(3-(didecylamino)propanamide)
MF: $C_{56}H_{114}N_6O_2$
FW: 903.6
Purity: $\geq 95\%$
Supplied as: A solution in ethanol
Storage: -20°C
Stability: ≥ 2 years



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

Description

PPZ-A10 is an ionizable cationic lipid.^{1,2} It has been used in the generation of lipid nanoparticles (LNPs) for the delivery of siRNA and mRNA *in vitro* and *in vivo*. Intraperitoneal administration of LNPs containing PPZ-A10 and encapsulating an mRNA reporter preferentially accumulates in hepatic Kupffer cells and splenic macrophages in mice.²

References

1. Subhan, M.A. and Torchilin, V.P. Biopolymer-based nanosystems for siRNA drug delivery to solid tumors including breast cancer. *Pharmaceutics* **15**(1), 153 (2023).
2. Ni, H., Hatit, M.Z.C., Zhao, K., *et al.* Piperazine-derived lipid nanoparticles deliver mRNA to immune cells *in vivo*. *Nat. Commun.* **13**(1), 4766 (2022).

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, 12/11/2023

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