

PRODUCT INFORMATION



SAINT C-18

Item No. 39709

CAS Registry No.: 187587-97-9

Formal Name: 1-methyl-4-[(10Z)-1-(9Z)-9-octadecen-1-yl-10-nonadecen-1-yl]-pyridinium, monochloride

Synonyms: SAINT-2, SAINT-7, SAINT-18, Synthetic Amphiphile INTERaction C-18

MF: $C_{43}H_{78}N \cdot Cl$

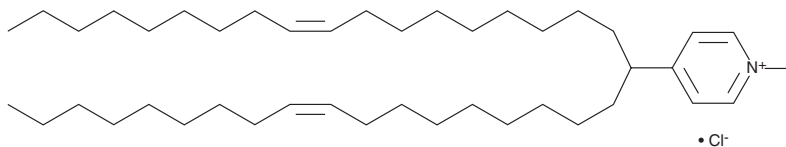
FW: 644.5

Purity: $\geq 98\%$

Supplied as: A solution in chloroform

Storage: $-20^{\circ}C$

Stability: ≥ 1 year



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

Laboratory Procedures

SAINT C-18 is supplied as a solution in chloroform. To change the solvent, simply evaporate the chloroform under a gentle stream of nitrogen and immediately add the solvent of choice, which should be purged with an inert gas. SAINT C-18 is slightly soluble in acetonitrile.

Description

SAINT C-18 is a cationic lipid that has been used in the formation of liposomes for the delivery of plasmid DNA, proteins, siRNA, and small molecules.¹⁻⁴ Complexes of SAINT C-18 with plasmid DNA encoding pigment epithelium-derived factor (PEDF) inhibit corneal neovascularization induced by complexes of SAINT C-18 with plasmid DNA encoding FGF2, also known as basic FGF, in rats.²

References

1. van der Woude, I., Wagenaar, A., Meekel, A.A.P., *et al.* Novel pyridinium surfactants for efficient, nontoxic *in vitro* gene delivery. *Proc. Natl. Acad. Sci. USA* **94**(4), 1160-1165 (1997).
2. Kuo, C.-N., Yang, L.-C., Yang, C.-T., *et al.* Inhibition of corneal neovascularization with plasmid pigment epithelium-derived factor (p-PEDF) delivered by synthetic amphiphile INTERaction-18 (SAINT-18) vector in an experimental model of rat corneal angiogenesis. *Exp. Eye Res.* **89**(5), 678-685 (2009).
3. van der Gun, B.T.F., Monami, A., Laarmann, S., *et al.* Serum insensitive, intranuclear protein delivery by the multipurpose cationic lipid SAINT-2. *J. Control Release* **123**(3), 228-238 (2007).
4. Adrian, J.E., Morselt, H.W., Süß, R., *et al.* Targeted SAINT-O-Somes for improved intracellular delivery of siRNA and cytotoxic drugs into endothelial cells. *J. Control Release* **144**(3), 341-349 (2010).

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, 03/26/2024

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