

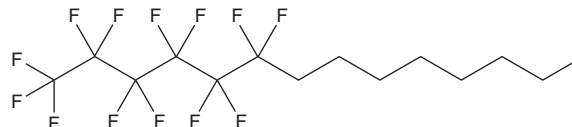
PRODUCT INFORMATION



Perfluorohexyloctane

Item No. 40490

CAS Registry No.: 133331-77-8
Formal Name: 1,1,1,2,2,3,3,4,4,5,5,6,6-tridecafluoro-tetradecane
Synonyms: F6H8, NOV03,
1-(Perfluorohexyl)octane
MF: C₁₄H₁₇F₁₃
FW: 432.3
Purity: ≥95%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Perfluorohexyloctane is supplied as a solid. A stock solution may be made by dissolving the perfluorohexyloctane in the solvent of choice, which should be purged with an inert gas. Perfluorohexyloctane is sparingly soluble (1-10 mg/ml) in ethanol.

Description

Perfluorohexyloctane is a semifluorinated alkane and perfluoroalkyl substance (PFAS).¹ It improves corneal penetration of cyclosporine A (Item No. 12088) in isolated porcine eyes when used at concentrations of 0.05 and 0.1%. Topical application of perfluorohexyloctane improves the lipid layer grade in rabbit eyes without inducing irritation.² Formulations containing perfluorohexyloctane have been used in the treatment of dry eye disease.

References

1. Agarwal, P., Scherer, D., Günther, B., *et al.* Semifluorinated alkane based systems for enhanced corneal penetration of poorly soluble drugs. *Int. J. Pharm.* **538(1-2)**, 119-129 (2018).
2. Agarwal, P., Khun, D., Krösser, S., *et al.* Preclinical studies evaluating the effect of semifluorinated alkanes on ocular surface and tear fluid dynamics. *Ocul. Surf.* **17(2)**, 241-249 (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, 12/19/2024

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