

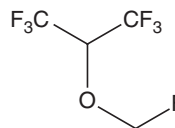
PRODUCT INFORMATION



Sevoflurane

Item No. 23996

CAS Registry No.: 28523-86-6
Formal Name: 1,1,1,3,3,3-hexafluoro-2-(fluoromethoxy)-propane
Synonyms: Fluoromethyl 1,1,1,3,3,3-Hexafluoroisopropyl ester, Fluoromethyl 1,1,1,3,3,3-Hexafluoro-2-propyl ester
MF: C₄H₃F₇O
FW: 200.1
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Sevoflurane is supplied as a neat oil. A stock solution may be made by dissolving the sevoflurane in the solvent of choice, which should be purged with an inert gas. Sevoflurane is miscible in organic solvents such as ethanol, DMSO, and dimethyl formamide.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of sevoflurane can be prepared by directly dissolving the neat oil in aqueous buffers. Sevoflurane is miscible in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Sevoflurane is a halogenated ether with anesthetic properties.¹ It enhances the activity of GABA_A and glycine receptors and inhibits the activity of nicotinic acetylcholine receptors (nAChRs) and glutamate receptors. Sevoflurane enhances the responses of α₂β₁ subunit-containing GABA_A and α₁ subunit-containing glycine receptors at submaximal agonist concentrations in HEK293 cells (EC₅₀s = 0.45 and 0.36 mM, respectively).² Sevoflurane (360 μM) also increases the amplitude of GABA_A receptor responses to GABA stimulation for receptors with an α₁β₂γ₂ subunit composition.³ It inhibits binding of the high affinity nicotinic agonist epibatidine to nAChRs in mouse brain membranes (IC₅₀ = 0.77 mM).⁴ Formulations containing sevoflurane have been used in the induction and maintenance of general anesthesia.

References

1. Campagna, J.A., Miller, K.W., Phil, D., *et al.* Mechanisms of actions of inhaled anesthetics. *N. Engl. J. Med.* **348(21)**, 2110-2124 (2003).
2. Krasowski, M.D. and Harrison, N.L. The actions of ether, alcohol and alkane general anaesthetics on GABA_A and glycine receptors and the effects of TM2 and TM3 mutations. *Br. J. Pharmacol.* **129(4)**, 731-743 (2000).
3. Nishikawa, K. and Harrison, N.L. The actions of sevoflurane and desflurane on the gamma-aminobutyric acid receptor type A: Effects of TM2 mutations in the alpha and beta subunits. *Anesthesiology* **99(3)**, 678-684 (2003).
4. Rada, E.M., Tharakan, E.C., and Flood, P. Volatile anesthetics reduce agonist affinity at nicotinic acetylcholine receptors in the brain. *Anesth. Analg.* **96(1)**, 108-111 (2003).

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, 11/08/2022

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