

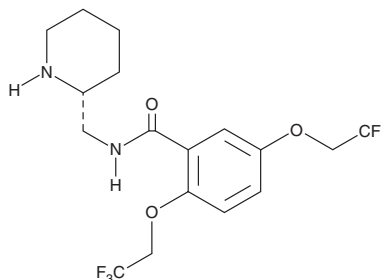
PRODUCT INFORMATION



R-(-)-Flecainide

Item No. 20682

CAS Registry No.: 99495-90-6
Formal Name: N-[(2R)-2-piperidinylmethyl]-2,5-bis(2,2,2-trifluoroethoxy)benzamide
MF: C₁₇H₂₀F₆N₂O₃
FW: 414.3
Purity: ≥98%
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

R-(-)-Flecainide is supplied as a solid. A stock solution may be made by dissolving the R-(-)-flecainide in the solvent of choice, which should be purged with an inert gas. R-(-)-Flecainide is soluble in organic solvents such as ethanol and methanol.

Description

R-(-)-Flecainide is the (R) enantiomer of the antiarrhythmic agent flecainide (Item No. 20388).¹ R-(-)-Flecainide prevents chloroform-induced ventricular fibrillation in mice (ED₅₀ = 14.5 mg/kg). It also prevents ouabain-induced ectopic ventricular tachycardia in anesthetized dogs.

Reference

1. Banitt, E.H., Schmid, J.R., and Newmark, R.A. Resolution of flecainide acetate, N-(2-piperidylmethyl)-2,5-bis(2,2,2-trifluoroethoxy)benzamide acetate, and antiarrhythmic properties of the enantiomers. *J. Med. Chem.* **29**(2), 299-302 (1986).

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, 12/15/2022

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