

PRODUCT INFORMATION



Dofetilide N-oxide

Item No. 36294

CAS Registry No.: 144449-71-8

Formal Name: N-[4-[2-[methyl[2-[4-[(methylsulfonyl)amino]phenoxy]ethyl]oxidoamino]ethyl]phenyl]-methanesulfonamide

Synonyms: CS 0019807, UK 116856

MF: $C_{19}H_{27}N_3O_6S_2$

FW: 457.6

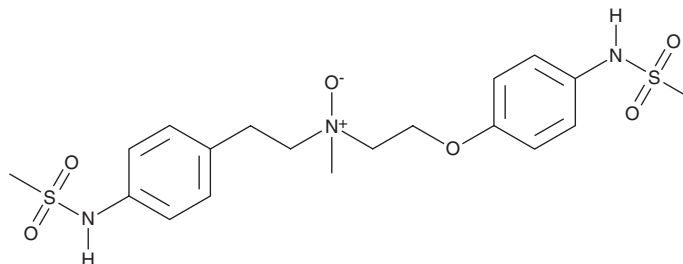
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 231 nm

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

Dofetilide N-oxide is supplied as a solid. A stock solution may be made by dissolving the dofetilide N-oxide in the solvent of choice, which should be purged with an inert gas. Dofetilide N-oxide is soluble in the organic solvent DMSO at a concentration of approximately 1 mg/ml.

Description

Dofetilide N-oxide is a metabolite of the antiarrhythmic agent dofetilide (Item No. 15045).¹ It prolongs intracellular action potential duration in isolated guinea pig papillary muscle, as a measure of antiarrhythmic activity, less potently than dofetilide.

Reference

1. Walker, D.K., Alabaster, C.T., Congrave, G.S., *et al.* Significance of metabolism in the disposition and action of the antidysrhythmic drug, dofetilide. *In vitro* studies and correlation with *in vivo* data. *Drug Metab. Dispos.* **24**(4), 447-455 (1996).

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, 11/18/2024

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