

PRODUCT INFORMATION



AM92016 (hydrochloride)

Item No. 36803

CAS Registry No.: 133229-11-5
Formal Name: N-[4-[3-[[2-(3,4-dichlorophenyl)ethyl]methylamino]-2-hydroxypropoxy]phenyl]-methanesulfonamide, monohydrochloride

MF: C₁₉H₂₄Cl₂N₂O₄S • HCl

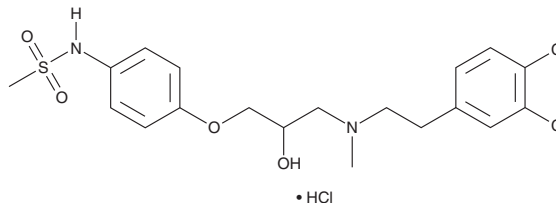
FW: 483.8

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

AM92016 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the AM92016 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. AM92016 (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of AM92016 (hydrochloride) in ethanol and DMSO is approximately 1 mg/ml and approximately 3 mg/ml in DMF.

Description

AM92016 is an inhibitor of delayed-rectifier potassium current (I_{Kr} ; $IC_{50} = 40$ nM in isolated rabbit sinoatrial node cells).¹ It is selective for I_{Kr} over L-type voltage-gated calcium (Ca_v) current and hyperpolarization-activated inward current (I_h) at 1 μ M. AM92016 (50 nM) decreases the diastolic depolarization rate and increases cycle length and action potential duration at full repolarization in isolated rabbit sinoatrial node cells. It decreases the QT interval and increases heart rate in anesthetized guinea pigs when administered at a dose of 1 mg/kg.² AM92016 (1 mg/kg) decreases the time to first arrhythmia and ventricular fibrillation in ouabain-stimulated guinea pigs.

References

1. Lei, M. and Brown, H.F. Inhibition by Compound II, a sotalol analogue, of delayed rectifier current (i_{Kr}) in rabbit isolated sino-atrial node cells. *Naunyn Schmiedebergs Arch Pharmacol.* **357(3)**, 260-267 (1998).
2. Hagerty, M.J., Wainwright, C.L., and Kane, K.A. The in-vivo cardiovascular effects of a putative class III anti-arrhythmic drug, AM 92016. *J. Pharm. Pharmacol.* **48(4)**, 417-421 (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, 12/21/2022

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