

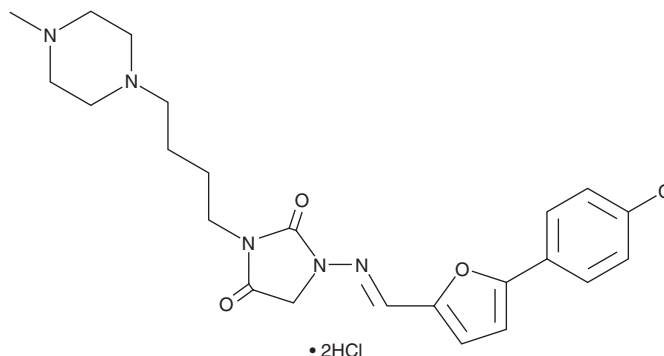
# PRODUCT INFORMATION



## Azimilide (hydrochloride)

Item No. 28300

**CAS Registry No.:** 149888-94-8  
**Formal Name:** 1-[[[5-(4-chlorophenyl)-2-furanyl]methylene]amino]-3-[4-(4-methyl-1-piperazinyl)butyl]-2,4-imidazolidinedione, dihydrochloride  
**Synonym:** NE 10064  
**MF:** C<sub>23</sub>H<sub>26</sub>ClN<sub>5</sub>O<sub>3</sub> • 2HCl  
**FW:** 530.9  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 233, 257, 343 nm  
**Supplied as:** A crystalline 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

Azimilide (hydrochloride) is supplied as a crystalline solid. Aqueous solutions of azimilide (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of azimilide (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Azimilide is a class III antiarrhythmic agent.<sup>1</sup> It blocks rapidly activating delayed rectifier (I<sub>Kr</sub>) and slowly activating delayed rectifier (I<sub>Ks</sub>) potassium currents in isolated guinea pig ventricular myocytes (IC<sub>50</sub>s = 0.3 and 3 μM, respectively). Azimilide (0.1-10 μM) prolongs action potential duration (APD) in isolated calf Purkinje fibers and ventricular trabeculae in a concentration-dependent manner. It increases the effective refractory period (ERP) in isolated perfused guinea pig hearts and ferret papillary muscles. Azimilide (0.3-30 mg/kg) increases ERP and the absolute refractory period and decreases heart rate in dogs. It reduces mean arrhythmia score in a rat model of coronary artery ligation and reperfusion-induced severe ventricular arrhythmia. Azimilide (10-30 mg/kg) also prevents sustained and non-sustained ventricular tachyarrhythmias in dogs.

### Reference

1. Salata, J.J. and Brooks, R.R. Pharmacology of azimilide dihydrochloride (NE-10064), a class III antiarrhythmic agent. *Cardiovasc. Drug Rev.* **15(2)**, 137-156 (2007).

#### 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, 11/02/2022

#### 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