

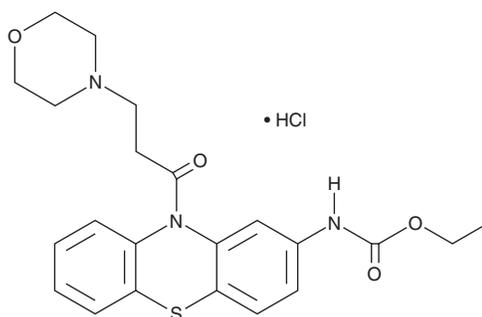
# PRODUCT INFORMATION



## Moricizine (hydrochloride)

Item No. 45203

**CAS Registry No.:** 29560-58-5  
**Formal Name:** N-[10-[3-(4-morpholinyl)-1-oxopropyl]-10H-phenothiazin-2-yl]-carbamic acid, ethyl ester, monohydrochloride  
**Synonyms:** EN-313, Moracizine  
**MF:** C<sub>22</sub>H<sub>25</sub>N<sub>3</sub>O<sub>4</sub>S • HCl  
**FW:** 464.0  
**Purity:** ≥95%  
**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

Moricizine (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the moricizine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Moricizine (hydrochloride) is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

### Description

Moricizine is a class I antiarrhythmic agent.<sup>1</sup> It induces both tonic and phasic blocks of sodium currents in isolated guinea pig atrial myocytes when used at a concentration of 30 μM and induces a tonic block of sodium currents in isolated guinea pig ventricular myocytes at 1 and 10 μM. Moricizine (1-3 mg/kg) reduces spontaneous arrhythmias occurring 24 hours after ligation of the left anterior descending (LAD) coronary artery in dogs as a model of myocardial infarction.<sup>2</sup> It also reduces arrhythmias occurring 2-5 days after LAD coronary ligation in the same model. Formulations containing moricizine have previously been used in the treatment of ventricular arrhythmias.

### References

1. Ahmmed, G.U., Hisatome, I., Kurata, Y., *et al.* Analysis of moricizine block of sodium current in isolated guinea-pig atrial myocytes. Atrioventricular difference of moricizine block. *Vascul. Pharmacol.* **38(3)**, 131-141 (2002).
2. Danilo, P., Jr., Langan, W.B., Rosen, M.R., *et al.* Effects of the phenothiazine analog, EN-313, on ventricular arrhythmias in the dog. *Eur. J. Pharmacol.* **45(2)**, 127-139 (1977).

#### 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, 02/23/2026

#### 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