

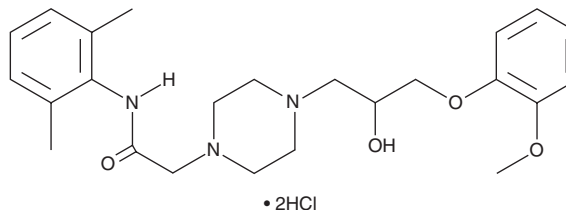
# PRODUCT INFORMATION



## Ranolazine (hydrochloride)

Item No. 15604

**CAS Registry No.:** 95635-56-6  
**Formal Name:** N-(2,6-dimethylphenyl)-4-[2-hydroxy-3-(2-methoxyphenoxy)propyl]-1-piperazineacetamide, dihydrochloride  
**Synonym:** RS 43285  
**MF:** C<sub>24</sub>H<sub>33</sub>N<sub>3</sub>O<sub>4</sub> • 2HCl  
**FW:** 500.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 272 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

Ranolazine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the ranolazine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Ranolazine (hydrochloride) is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ranolazine (hydrochloride) in these solvents is approximately 20 mg/ml.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ranolazine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ranolazine (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

Ranolazine is a piperazine derivative with cardioprotective activity.<sup>1-4</sup> It reduces the late sodium current ( $I_{Na^L}$ ) in mouse myocytes expressing the long QT syndrome 3 mutant sodium channel DKPQ, ventricular myocytes isolated from a canine model of heart failure, guinea pig ventricular myocytes exposed to hydrogen peroxide or anemone toxin-II, and HEK293 cells expressing human Na<sub>v</sub>1.5 channels ( $IC_{50}$ s = 5.9-15 μM) as well as the late potassium current ( $I_{Kr}$ ) in canine ventricular myocytes and HEK293 cells ( $IC_{50}$ s = 11.5 and 14.4 μM, respectively).<sup>1,2</sup> Ranolazine also inhibits radioligand binding to α<sub>1</sub>-, β<sub>1</sub>-, and β<sub>2</sub>-adrenergic receptors ( $K_i$ s = 8.2-19.5, 1.4-8.6, and 0.5-14.8 μM, respectively).<sup>2</sup> *In vivo*, ranolazine (480 μg/kg per min) reduces clofilium-induced prolongation of the QTc interval and Torsade de Pointes (TdP) in rabbits.<sup>3</sup> Ranolazine also reduces interstitial collagen deposition as well as atrial natriuretic peptide (ANP; Item Nos. 24539 | 24276), connective tissue growth factor (CTGF), brain natriuretic peptide (BNP; Item No. 24541), and matrix metalloproteinase-2 (MMP-2) mRNA levels, and prevents left ventricular dilation in a mouse model of cardiotoxicity induced by doxorubicin (Item No. 15007).<sup>4</sup>

### References

1. Shryock, J.C. and Belardinelli, L. *Br. J. Pharmacol.* **153**(6), 1128-1132 (2008).
2. Verrier, R.L., Kumar, K., Nieminen, T., et al. *Europace* **15**(3), 317-324 (2013).
3. Wang, W.Q., Robertson, C., Dhalla, A.K., et al. *J. Pharmacol. Exp. Ther.* **325**(3), 875-881 (2008).
4. Tocchetti, C.G., Carpi, A., Coppola, C., et al. *Eur. J. Heart Fail.* **16**(4), 358-366 (2014).

#### 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, 12/12/2022

#### 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