

PRODUCT INFORMATION

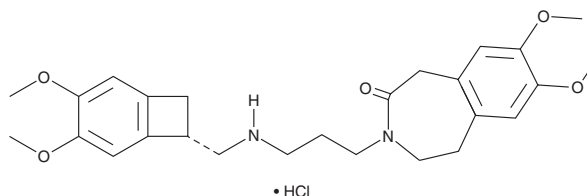


N-desmethyl Ivabradine (hydrochloride)

Item No. 27808

CAS Registry No.: 1246638-08-3
Formal Name: 3-[3-[[[(7S)-3,4-dimethoxybicyclo[4.2.0]octa-1,3,5-trien-7-yl]methyl]amino]propyl]-1,3,4,5-tetrahydro-7,8-dimethoxy-2H-3-benzazepin-2-one, monohydrochloride

Synonym: S18982
MF: C₂₆H₃₄N₂O₅ • HCl
FW: 491.0
Purity: ≥95%
UV/Vis.: λ_{max}: 287 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

N-desmethyl Ivabradine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desmethyl ivabradine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. N-desmethyl Ivabradine (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of N-desmethyl ivabradine (hydrochloride) in these solvents is approximately 1, 20, and 25 mg/ml, respectively.

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 N-desmethyl ivabradine (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N-desmethyl ivabradine (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

N-desmethyl Ivabradine is an active metabolite of ivabradine (Item No. 15868).¹ Ivabradine is metabolized by the cytochrome P450 (CYP) isoform CYP3A4.

Reference

1. Rosa, G.M., Ferrero, S., Ghione, P., et al. An evaluation of the pharmacokinetics and pharmacodynamics of ivabradine for the treatment of heart failure. *Expert Opin. Drug Metab. Toxicol.* **10(2)**, 279-291 (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, 12/21/2022

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