

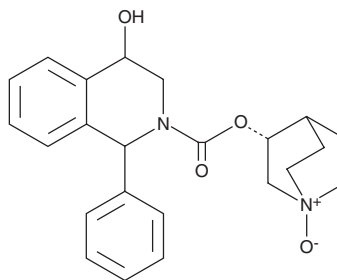
PRODUCT INFORMATION



4-hydroxy Solifenacin N-oxide

Item No. 26649

CAS Registry No.: 2747991-39-3
Formal Name: (3R)-3-((4-hydroxy-1-phenyl-1,2,3,4-tetrahydroisoquinoline-2-carbonyl)oxy)quinuclidine 1-oxide
MF: C₂₃H₂₆N₂O₄
FW: 394.5
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

4-hydroxy Solifenacin N-oxide is supplied as a solid. A stock solution may be made by dissolving the 4-hydroxy solifenacin N-oxide in the solvent of choice, which should be purged with an inert gas. 4-hydroxy Solifenacin N-oxide is soluble in the organic solvent methanol.

Description

4-hydroxy Solifenacin N-oxide is a metabolite of the muscarinic M₁, M₂, and M₃ receptor antagonist solifenacin (Item No. 17320).¹ It is formed from solifenacin by the cytochrome P450 (CYP) isoform CYP3A4 in human liver microsomes.

Reference

1. Michel, M.C., Yanagihara, T., Minematsu, T., *et al.* Disposition and metabolism of solifenacin in humans. *Brit. J. Clin. Pharmacol.* **59**(5), 647 (2005).

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, 11/18/2024

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