

PRODUCT INFORMATION



Solifenacin N-oxide

Item No. 26648

CAS Registry No.: 180272-28-0
Formal Name: (1S)-3,4-dihydro-1-phenyl-2(1H)-isoquinolinecarboxylic acid, (3R)-1-oxido-1-azabicyclo[2.2.2]oct-3-yl ester

MF: C₂₃H₂₆N₂O₃

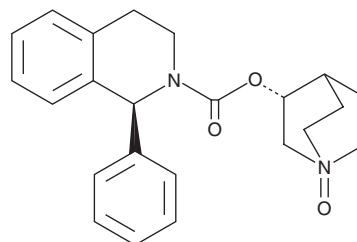
FW: 378.5

Purity: ≥98%

Supplied as: A crystalline solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Solifenacin N-oxide is supplied as a crystalline solid. A stock solution may be made by dissolving the solifenacin N-oxide in the solvent of choice. Solifenacin N-oxide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of solifenacin N-oxide in these solvents is approximately 15, 2, and 10 mg/ml, respectively.

Description

Solifenacin N-oxide is an inactive metabolite of the muscarinic receptor antagonist solifenacin (Item No. 17320).¹ It is also a potential impurity found in commercial preparations of solifenacin.² Solifenacin N-oxide is formed when solifenacin is stored under oxidative conditions.

References

1. Krauwinkel, W.J., Smulders, R.A., Mulder, H., *et al.* Effect of age on the pharmacokinetics of solifenacin in men and women. *Int. J. Clin. Pharmacol. Ther.* **43(5)**, 227-238 (2005).
2. Ganthi, H.K.R., Reddy, R.P., Park, Y.J., *et al.* Stability indicating HPLC method for quantification of solifenacin succinate & tamsulosin hydrochloride along with its impurities in tablet dosage form. *Am. J. Analyt. Chem.* **7(11)**, 840-862 (2016).

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, 07/19/2019

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