

PRODUCT INFORMATION



Solifenacin N-Glucuronide

Item No. 26650

Formal Name: (2S,3S,4S,5R,6R)-3,4,5-trihydroxy-6-(((3R)-3-(((S)-1-phenyl-1,2,3,4-tetrahydroisoquinoline-2-carbonyl)oxy)quinuclidin-1-ium-1-yl)tetrahydro-2H-pyran-2-carboxylate

MF: C₂₉H₃₄N₂O₈

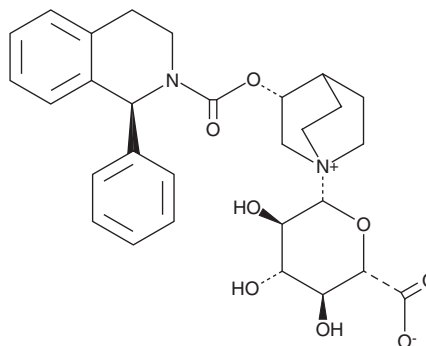
FW: 538.6

Purity: ≥85%

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

Solifenacin N-glucuronide is supplied as a solid. A stock solution may be made by dissolving the solifenacin N-glucuronide in the solvent of choice, which should be purged with an inert gas. Solifenacin N-glucuronide is soluble in organic solvents such as methanol and DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

Solifenacin N-glucuronide is a metabolite of the muscarinic receptor antagonist solifenacin (Item No. 17320) formed by direct glucuronidation.¹

Reference

1. Kuipers, M., Smulders, R., Krauwinkel, W., *et al.* Open-label study of the safety and pharmacokinetics of solifenacin in subjects with hepatic impairment. *J. Pharmacol. Sci.* **102(4)**, 405-412 (2006).

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, 11/29/2022

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