

PRODUCT INFORMATION

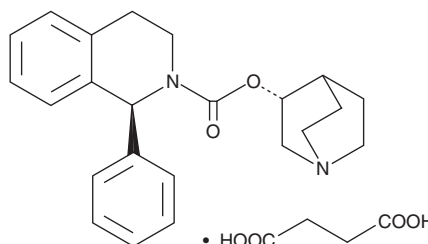


Solifenacin (succinate)

Item No. 17320

CAS Registry No.: 242478-38-2
Formal Name: butanedioic acid, compd. with (1S)-(3R)-1-azabicyclo[2.2.2]oct-3-yl 3,4-dihydro-1-phenyl-2(1H)-isoquinolinecarboxylate

Synonym: YM-67905
MF: C₂₃H₂₆N₂O₂ • C₄H₆O₄
FW: 480.6
Purity: ≥98%
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

Solifenacin (succinate) is supplied as a crystalline solid. A stock solution may be made by dissolving the solifenacin (succinate) in the solvent of choice, which should be purged with an inert gas. Solifenacin (succinate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of solifenacin (succinate) in ethanol is approximately 5 mg/ml and approximately 25 mg/ml in DMSO and DMF.

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 solifenacin (succinate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of solifenacin (succinate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Solifenacin is a competitive antagonist of M₁, M₂, and M₃ muscarinic acetylcholine receptors (K_is = 25, 125, and 10 nM, respectively, for the human receptors).¹ It inhibits calcium mobilization induced by carbachol (carbamoylcholine; Item No. 14486) in isolated guinea pig detrusor muscle cells (K_i = 4 nM).² Solifenacin inhibits carbachol-induced contraction of isolated guinea pig urinary bladder smooth muscle. *In vivo*, solifenacin (0.03-1 mg/kg) inhibits carbachol-induced increases in urinary bladder pressure in anesthetized rats. Formulations containing solifenacin have been used in the treatment of overactive bladder.

References

1. Hegde, S.S. Muscarinic receptors in the bladder: From basic research to therapeutics. *Br. J. Pharmacol.* **147(Suppl 2)**, S80-S87 (2006).
2. Ikeda, K., Kobayashi, S., Suzuki, M., *et al.* M₃ receptor antagonism by the novel antimuscarinic agent solifenacin in the urinary bladder and salivary gland. *Naunyn Schmiedebergs Arch. Pharmacol.* **366(2)**, 97-103 (2002).

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, 03/21/2024

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