

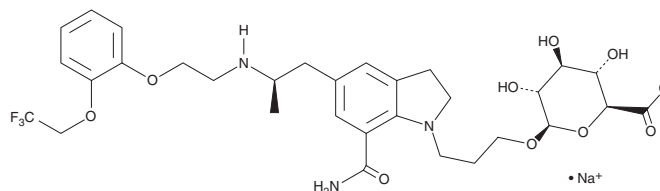
PRODUCT INFORMATION



Silodosin Glucuronide (sodium salt)

Item No. 28045

CAS Registry No.: 879292-24-7
Formal Name: 3-[7-(aminocarbonyl)-2,3-dihydro-5-[(2R)-2-[[2-[2-(2,2,2-trifluoroethoxy)phenoxy]ethyl]amino]propyl]-1H-indol-1-yl]propyl, β -D-glucopyranosiduronic acid, monosodium salt
Synonym: Silodosin β -D-glucuronide
MF: $C_{31}H_{39}F_3N_3O_{10} \cdot Na$
FW: 693.6
Purity: $\geq 90\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Silodosin glucuronide (sodium salt) is supplied as a solid. A stock solution may be made by dissolving the silodosin glucuronide (sodium salt) in the solvent of choice, which should be purged with an inert gas. Silodosin glucuronide (sodium salt) is slightly soluble in methanol.

Description

Silodosin glucuronide is an active metabolite of the α_{1A} -adrenergic receptor antagonist silodosin (Item No. 14866).^{1,2} It is formed from silodosin by the UDP-glucuronosyltransferase (UGT) isoform UGT2B7.¹ Silodosin glucuronide is toxic to rats with an LD₅₀ value of 0.347 mg/kg.²

References

1. Keating, G.M. Silodosin: A review of its use in the treatment of the signs and symptoms of benign prostatic hyperplasia. *Drugs* **75(2)**, 207-217 (2015).
2. Vishnuvardhan, C., Baikadi, S., Borkar, R.M., et al. *In vivo* metabolic investigation of silodosin using UHPLC-QTOF-MS/MS and in silico toxicological screening of its metabolites. *J. Mass. Spectrom.* **51(10)**, 867-882 (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, 12/19/2022

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