

PRODUCT INFORMATION



N-desmethyl Sildenafil (citrate)

Item No. 10010587

CAS Registry No.: 2319839-04-6
Formal Name: 5-[2-ethoxy-5-(1-piperazinylsulfonyl)phenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidin-7-one, 2-hydroxypropane-1,2,3-tricarboxylic acid

MF: C₂₁H₂₈N₆O₄S • C₆H₈O₇

FW: 652.7

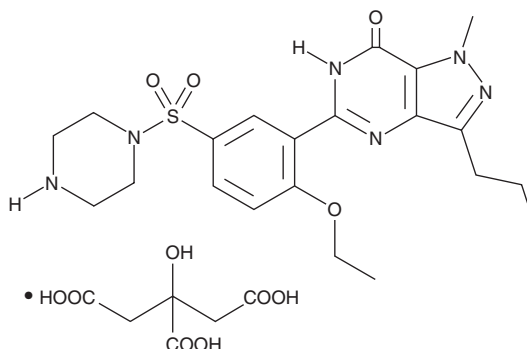
Purity: ≥98%

UV/Vis.: λ_{max}: 212, 295 nm

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

N-desmethyl Sildenafil (citrate) is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desmethyl sildenafil (citrate) in the solvent of choice. N-desmethyl Sildenafil (citrate) is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of N-desmethyl sildenafil (citrate) in these solvents is approximately 10 and 14 mg/ml, respectively.

Description

N-desmethyl Sildenafil is a major metabolite of sildenafil (Item Nos. 10008671 | 14008).¹⁻³ N-desmethyl Sildenafil is formed *via* oxidative metabolism of sildenafil by the cytochrome (CYP) P450 isoforms CYP3A4, CYP3A5, and CYP3A7.³

References

1. Walker, D.K., Ackland, M.J., James, G.C., *et al.* Pharmacokinetics and metabolism of sildenafil in mouse, rat, rabbit, dog and man. *Xenobiotica* **29(3)**, 297-310 (2015).
2. Muirhead, G.J., Wilner, K., Colburn, W., *et al.* The effects of age and renal and hepatic impairment on the pharmacokinetics of sildenafil citrate. *Br. J. Clin. Pharmacol.* **53(Suppl 1)**, 21S-30S (2015).
3. Takahiro, R., Nakamura, S., Kohno, H., *et al.* Contribution of CYP3A isoforms to dealkylation of PDE5 inhibitors: A comparison between sildenafil N-demethylation and tadalafil demethylation. *Biol. Pharm. Bull.* **38(1)**, 58-65 (2015).

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, 12/04/2024

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