

PRODUCT INFORMATION



N-Desethyl Vardenafil

Item No. 9001800

CAS Registry No.: 448184-46-1

Formal Name: 2-[2-ethoxy-5-(1-piperazinylsulfonyl)phenyl]-5-methyl-7-propyl-imidazo[5,1-f][1,2,4]triazin-4(1H)-one

MF: $C_{21}H_{28}N_6O_4S$

FW: 460.6

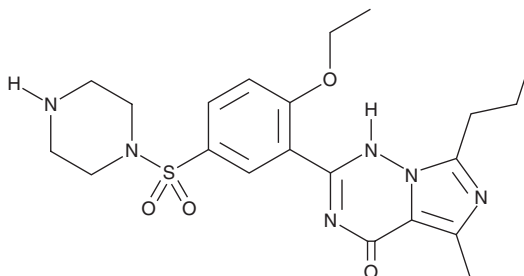
Purity: $\geq 98\%$

UV/Vis.: λ_{max} : 214 nm

Supplied as: A crystalline solid

Storage: $-20^{\circ}C$

Stability: As supplied, 2 years from the QC date provided on the Certificate of Analysis, when stored properly



Laboratory Procedures

N-Desethyl vardenafil is supplied as a crystalline solid. A stock solution may be made by dissolving the N-desethyl vardenafil in the solvent of choice. N-Desethyl vardenafil is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of N-desethyl vardenafil in these solvents is approximately 30 mg/ml.

N-Desethyl vardenafil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, N-desethyl vardenafil should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. N-Desethyl vardenafil has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

N-Desethyl vardenafil is a metabolite of vardenafil (Item No. 14930), a phosphodiesterase 5 inhibitor, that is formed by the action of cytochrome P450 (CYP)3A4 and CYP3A5.¹

Reference

1. Ku, H.-Y., Ahn, H.-J., Seo, J.-A., *et al.* The contributions of cytochromes P450 3A4 and 3A5 to the metabolism of the phosphodiesterase type 5 inhibitors sildenafil, udenafil, and vardenafil. *Drug Metab. Dispos.* **36**(6), 986-990 (2008).

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, 07/21/2016

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