

PRODUCT INFORMATION



N-desethyl-N-methyl Vardenafil

Item No. 35435

CAS Registry No.: 224785-87-9

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

MF: $C_{22}H_{30}N_6O_4S$

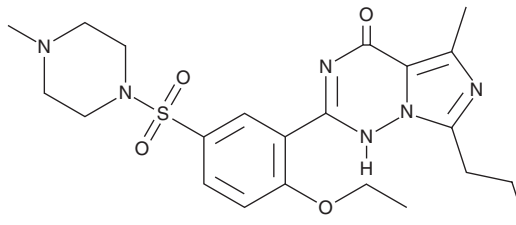
FW: 474.6

Purity: $\geq 95\%$

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

N-desethyl-N-methyl Vardenafil is supplied as a solid. A stock solution may be made by dissolving the N-desethyl-N-methyl vardenafil in the solvent of choice, which should be purged with an inert gas. N-desethyl-N-methyl Vardenafil is soluble in methanol and acetonitrile.

Description

N-desethyl-N-methyl Vardenafil is a potential impurity found in commercial preparations of the phosphodiesterase 5 (PDE5) inhibitor vardenafil (Item No. 14930).¹ It is also a PDE5 inhibitor ($IC_{50} = 2 \text{ nM}$).²

References

1. Subba Rao, D.V., Surendranath, K.V., Radhakrishnanand, P., *et al.* A stability indicating LC method for vardenafil HCl. *Chromatographia* **68**(9), 829–835 (2008).
2. Haning, H., Niewöhner, U., Schenke, T., *et al.* Comparison of different heterocyclic scaffolds as substrate analog PDE5 inhibitors. *Bioorg. Med. Chem. Lett.* **15**(17), 3900-3907 (2005).

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, 12/12/2022

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