

# PRODUCT INFORMATION



## Noracetildenafil

Item No. 9001803

CAS Registry No.: 949091-38-7

Formal Name: 5-[2-ethoxy-5-[2-(4-methyl-1-piperazinyl)acetyl]phenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidin-7-one

Synonym: Demethylhongdenafil

MF:  $C_{24}H_{32}N_6O_3$

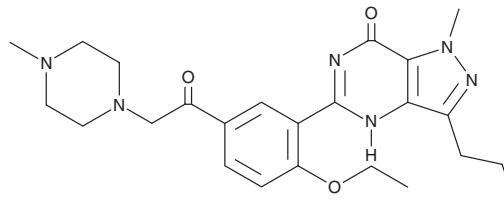
FW: 452.6

Purity:  $\geq 95\%$

Supplied as: A solid

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

Stability:  $\geq 4$  years



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

### Laboratory Procedures

Noracetildenafil is supplied as a solid. A stock solution may be made by dissolving the noracetildenafil in the solvent of choice, which should be purged with an inert gas. Noracetildenafil is soluble in the organic solvent methanol.

### Description

Noracetildenafil is a derivative of the phosphodiesterase 5 (PDE5) inhibitor sildenafil (Item No. 10008671).<sup>1</sup>

### Reference

1. Lee, E.S., Lee, J.H., Han, K.M., *et al.* Simultaneous determination of 38 phosphodiesterase-5 inhibitors in illicit erectile dysfunction products by liquid chromatography-electrospray ionization-tandem mass spectrometry. *J. Pharm. Biomed. Anal.* **83**, 171-178 (2013).

#### 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, 12/01/2022

#### 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