

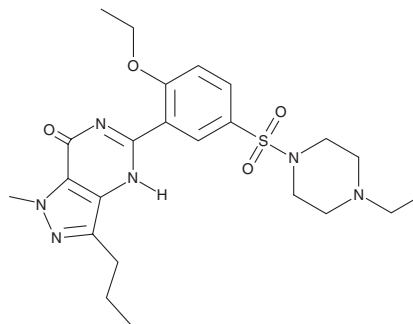
PRODUCT INFORMATION



Homosildenafil

Item No. 38891

CAS Registry No.: 642928-07-2
Formal Name: 5-[2-ethoxy-5-[(4-ethyl-1-piperazinyl)sulfonyl]phenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidin-7-one
Synonym: Methylsildenafil
MF: C₂₃H₃₂N₆O₄S
FW: 488.6
Purity: ≥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

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

Description

Homosildenafil is an inhibitor of phosphodiesterase 5 (PDE5; IC₅₀ = 8.9 nM).¹ It has been found as an adulterant in dietary supplements for erectile dysfunction.^{2,3}

References

1. Blount, M.A., Beasley, A., Zoraghi, R., *et al.* Binding of tritiated sildenafil, tadalafil, or vardenafil to the phosphodiesterase-5 catalytic site displays potency, specificity, heterogeneity, and cGMP stimulation. *Mol. Pharmacol.* **66**(1), 144-152 (2004).
2. Blok-Tip, L., Zomer, B., Bakker, F., *et al.* Structure elucidation of sildenafil analogues in herbal products. *Food Addit. Contam.* **21**(8), 737-748 (2004).
3. 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, 11/21/2023

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