

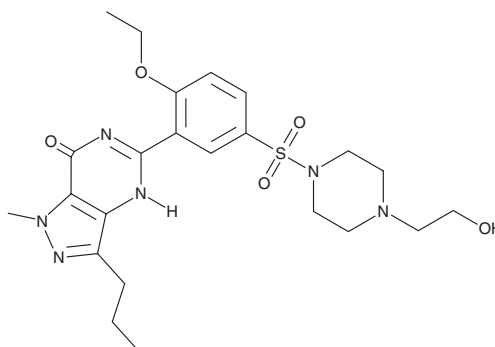
PRODUCT INFORMATION



Lodenafil

Item No. 42220

CAS Registry No.: 139755-85-4
Formal Name: 5-[2-ethoxy-5-[[4-(2-hydroxyethyl)-1-piperazinyl]sulfonyl]phenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidin-7-one
Synonym: Hydroxyhomosildenafilafil
MF: C₂₃H₃₂N₆O₅S
FW: 504.6
Purity: ≥98%
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

Lodenafil is supplied as a solid. A stock solution may be made by dissolving the lodenafil in the solvent of choice, which should be purged with an inert gas. Lodenafil is sparingly soluble (1-10 mg/ml) in DMSO.

Description

Lodenafil is an inhibitor of phosphodiesterase 5 (PDE5) and an active metabolite of lodenafil carbonate.^{1,2} It enhances acetylcholine-induced relaxation of isolated human corpus cavernosum and rabbit cavernosal strips precontracted with phenylephrine (EC₅₀s = 676 and 589 nM, respectively).²

References

1. Alshehri, Y.M., Al-Majed, A.A., Attwa, M.W., *et al.* Lodenafil. *Profiles of drug substances, excipients, and related methodology: Volume 47.* Al-Majed, A. editor, *Academic Press* 113-147 (2022).
2. Toque, H.A., Teixeira, C.E., Lorenzetti, R., *et al.* Pharmacological characterization of a novel phosphodiesterase type 5 (PDE5) inhibitor lodenafil carbonate on human and rabbit corpus cavernosum. *Eur. J. Pharmacol.* **591**(1-3), 189-195 (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, 10/30/2024

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