

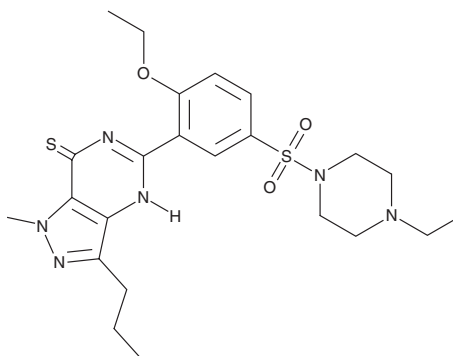
PRODUCT INFORMATION



Thiohomosildenafil

Item No. 46084

CAS Registry No.: 479073-80-8
Formal Name: 5-[2-ethoxy-5-[(4-ethyl-1-piperazinyl)sulfonyl]phenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidine-7-thione
Synonym: Sulfohomosildenafil
MF: $C_{23}H_{32}N_6O_3S_2$
FW: 504.7
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

Thiohomosildenafil is supplied as a solid. A stock solution may be made by dissolving the thiohomosildenafil in the solvent of choice, which should be purged with an inert gas. Thiohomosildenafil is slightly soluble (0.1-1 mg/ml) in DMSO.

Description

Thiohomosildenafil is a derivative of the phosphodiesterase 5 (PDE5) inhibitor sildenafil (Item No. 10008671).¹

Reference

1. Venhuis, B.J., Zomer, G., and de Kaste, D. Structure elucidation of a novel synthetic thiono analogue of sildenafil detected in an alleged herbal aphrodisiac. *J. Pharm. Biomed. Anal.* **46**(4), 814-817 (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, 07/17/2026

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