

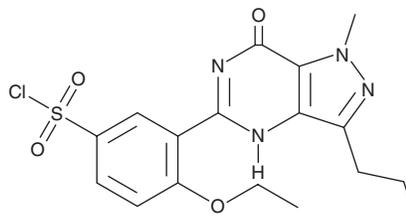
PRODUCT INFORMATION



Sildenafil Chlorosulfonyl

Item No. 34760

CAS Registry No.: 139756-22-2
Formal Name: 3-(6,7-dihydro-1-methyl-7-oxo-3-propyl-1H-pyrazolo[4,3-d]pyrimidin-5-yl)-4-ethoxy-benzenesulfonyl chloride
Synonym: CMP
MF: C₁₇H₁₉ClN₄O₄S
FW: 410.9
Purity: ≥98%
UV/Vis.: λ_{max}: 273 nm
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

Sildenafil chlorosulfonyl is supplied as a solid. A stock solution may be made by dissolving the sildenafil chlorosulfonyl in the solvent of choice, which should be purged with an inert gas. Sildenafil chlorosulfonyl is soluble in the organic solvent chloroform at a concentration of approximately 1 mg/ml. It is also slightly soluble in DMSO.

Description

Sildenafil chlorosulfonyl is a synthetic intermediate in the synthesis of the phosphodiesterase 5 (PDE5) inhibitor sildenafil (Item No. 10008671).¹ It is also a potential impurity found in commercial preparations of sildenafil.

Reference

1. Nagaraju, V., Sreenath, D., Rao, J.T., *et al.* Separation and determination of synthetic impurities of sildenafil (Viagra) by reversed-phase high-performance liquid chromatography. *Anal. Sci.* **19(7)**, 1007-1011 (2003).

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, 12/08/2022

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