

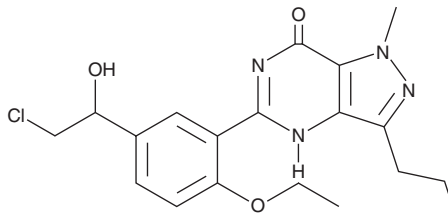
PRODUCT INFORMATION



Hydroxychlorodenafil

Item No. 29998

CAS Registry No.: 1391054-00-4
Formal Name: 5-[5-(2-chloro-1-hydroxyethyl)-2-ethoxyphenyl]-1,6-dihydro-1-methyl-3-propyl-7H-pyrazolo[4,3-d]pyrimidin-7-one
MF: C₁₉H₂₃ClN₄O₃
FW: 390.9
Purity: ≥98%
UV/Vis.: λ_{max}: 298 nm
Supplied as: A crystalline 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

Hydroxychlorodenafil is supplied as a crystalline solid. A stock solution may be made by dissolving the hydroxychlorodenafil in the solvent of choice, which should be purged with an inert gas. Hydroxychlorodenafil is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of hydroxychlorodenafil in these solvents is approximately 30 mg/ml.

Hydroxychlorodenafil is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, hydroxychlorodenafil should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Hydroxychlorodenafil has a solubility of approximately 0.16 mg/ml in a 1:5 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

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

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, 11/11/2022

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