

PRODUCT INFORMATION



Avanafil metabolite M16

Item No. 27844

Formal Name: 4-((4-((3-chloro-4-methoxybenzyl)amino)-5-((pyrimidin-2-ylmethyl)carbamoyl)pyrimidin-2-yl)amino)-5-hydroxypentanoic acid

MF: C₂₃H₂₆ClN₇O₅

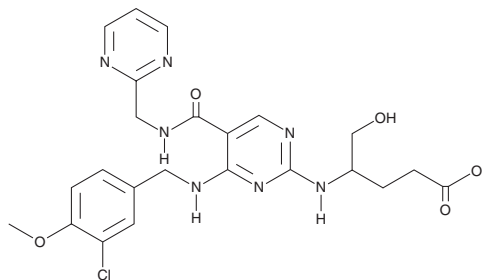
FW: 516.0

Purity: ≥95%

Supplied as: A solid

Storage: -20°C

Stability: ≥2 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Avanafil metabolite M16 is supplied as a solid. A stock solution may be made by dissolving the avanafil metabolite M16 in the solvent of choice. Avanafil metabolite M16 is soluble in the organic solvent DMSO, which should be purged with an inert gas.

Description

Avanafil metabolite M16 is a major inactive metabolite of the phosphodiesterase 5 (PDE5) inhibitor avanafil (Item No. 23024).¹ Avanafil is metabolized by the cytochrome P450 (CYP450) isoforms CYP3A4 and CYP2C to the major metabolites avanafil metabolite M16 and avanafil metabolite M4 (Item No. 27843), as well as minor metabolites.²

References

1. Kedia, G.T., Ückert, S., Assadi-Pour, F., et al. Avanafil for the treatment of erectile dysfunction: Initial data and clinical key properties. *Ther. Adv. Urol.* **5**(1), 35-41 (2013).
2. Kyle, J.A., Brown, D.A., and Hill, J.K. Avanafil for erectile dysfunction. *Ann. Pharmacother.* **47**(10), 1312-1320 (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, 04/29/2019

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