

PRODUCT INFORMATION



Gepotidacin

Item No. 44094

CAS Registry No.: 1075236-89-3
Formal Name: (2R)-2-[[4-[[[(3,4-dihydro-2H-pyrano[2,3-c]pyridin-6-yl)methyl]amino]-1-piperidinyl)methyl]-1,2-dihydro-3H,8H-2a,5,8a-triazaacenaphthylene-3,8-dione

Synonym: GSK2140944

MF: C₂₄H₂₈N₆O₃

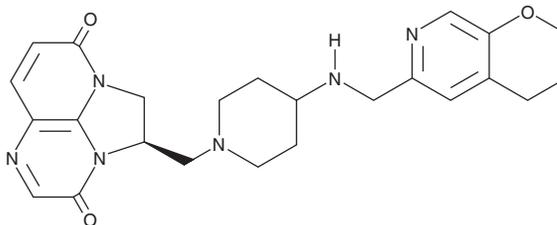
FW: 448.5

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

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

Gepotidacin is slightly soluble (0.1-1 mg/ml) in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Gepotidacin is an inhibitor of bacterial topoisomerase II.¹ It is active against 84 bacterial clinical isolates (MICs = 0.032-8 µg/ml) and vancomycin-resistant *S. aureus* (VRSA) and methicillin-resistant *S. aureus* (MRSA; MICs = 0.125 and 0.25 µg/ml, respectively).^{1,2} Gepotidacin reduces the number of lung colony-forming units (CFUs) in a mouse model of neutropenic *S. aureus* lung infection.² Formulations containing gepotidacin have been used in the treatment of uncomplicated urinary tract infections.

References

1. Waites, K.B., Crabb, D.M., Xiao, L., *et al.* *In vitro* activities of gepotidacin (GSK2140944) and other antimicrobial agents against human mycoplasmas and ureaplasmas. *Antimicrob. Agents Chemother.* **61**(10), e01064-17 (2017).
2. So, W., Crandon, J.L., and Nicolau, D.P. Pharmacodynamic profile of GSK2140944 against methicillin-resistant *Staphylococcus aureus* in a murine lung infection model. *Antimicrob. Agents Chemother.* **59**(8), 4956-4961 (2015).

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, 10/01/2025

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