

PRODUCT INFORMATION

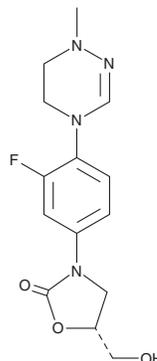


Delpazolid

Item No. 25532

CAS Registry No.: 1219707-39-7
Formal Name: (5R)-3-[4-(5,6-dihydro-1-methyl-1,2,4-triazin-4(1H)-yl)-3-fluorophenyl]-5-(hydroxymethyl)-2-oxazolidinone

Synonym: LCB01-0371
MF: C₁₄H₁₇FN₄O₃
FW: 308.3
Purity: ≥98%
UV/Vis.: λ_{max}: 238, 277 nm
Supplied as: A crystalline 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

Delpazolid is supplied as a crystalline solid. A stock solution may be made by dissolving the delpazolid in the solvent of choice. Delpazolid is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 30 mg/ml.

Description

Delpazolid is an oxazolidinone antibiotic.¹ It is active against a wide range of Gram-positive bacterial clinical isolates, including resistant strains of several species, such as methicillin-resistant *S. aureus* (MRSA), methicillin-resistant coagulase-negative staphylococci (MRCNS), and penicillin-resistant *S. pneumoniae* (MICs = 0.125-4 µg/ml). Delpazolid increases survival in mouse models of systemic MRSA, *E. faecalis*, *S. pneumoniae*, and *H. influenzae* infection (ED₅₀s = 2.96, 4.53, 2.28, and 9.96 mg/kg, respectively).

Reference

1. Jeong, J.-W., Jung, S.-J., Lee, H.-H., et al. *In vitro* and *in vivo* activities of LCB01-0371, a new oxazolidinone. *Antimicrob. Agents Chemother.* **54**(12), 5359-5362 (2010).

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, 03/20/2019

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