

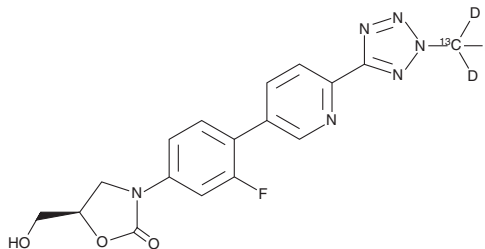
PRODUCT INFORMATION



Tedizolid-¹³C-d₃

Item No. 33290

CAS Registry No.: 2931763-72-1
Formal Name: (5R)-3-[3-fluoro-4-[6-[2-(methyl-¹³C-d₃)-2H-tetrazol-5-yl]-3-pyridinyl]phenyl]-5-(hydroxymethyl)-2-oxazolidinone
Synonym: TR-700-¹³C-d₃
MF: C₁₆[¹³C]H₁₂D₃FN₆O₃
FW: 374.4
Chemical Purity: ≥98% (Tedizolid)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₃); ≤1% d₀
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

Tedizolid-¹³C-d₃ is intended for use as an internal standard for the quantification of tedizolid (Item No. 32952) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated *versus* unlabeled).

Tedizolid-¹³C-d₃ is supplied as a solid. A stock solution may be made by dissolving the tedizolid-¹³C-d₃ in the solvent of choice, which should be purged with an inert gas. Tedizolid-¹³C-d₃ is soluble in methanol, DMSO, and acetonitrile.

Description

Tedizolid is an oxazolidinone antibiotic and active metabolite of the antibiotic tedizolid phosphate (Item No. 23729).¹ Tedizolid phosphate is converted to tedizolid by phosphatases *in vivo*. It is active against a wide variety of clinical isolates, including isolates of methicillin-resistant *S. aureus* (MRSA), *E. faecalis*, and penicillin-susceptible *S. pneumoniae* (MICs = 0.5, 0.5, and 0.25 µg/ml, respectively), as well as isolates of anaerobic bacteria (MICs = 0.25-4 µg/ml). Tedizolid protects against mortality in a mouse model of systemic methicillin-susceptible *S. aureus* (MSSA) infection (ED₅₀ = 8.8 mg/kg).²

References

1. Schaad, R., Sweeney, D.A., Shinabarger, D., *et al.* In vitro activity of TR-700, the active ingredient of the antibacterial prodrug TR-701, a novel oxazolidinone antibacterial agent. *Antimicrob. Agents Chemother.* **53**(8), 3236-3239 (2009).
2. Im, W.B., Choi, S.H., Park, J.-Y., *et al.* Discovery of torezolid as a novel 5-hydroxymethyl-oxazolidinone antibacterial agent. *Eur. J. Med. Chem.* **46**(4), 1027-1039 (2011).

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, 08/17/2023

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