

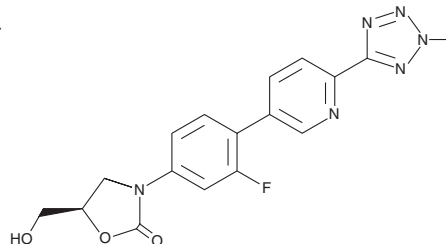
# PRODUCT INFORMATION



## Tedizolid

Item No. 32952

**CAS Registry No.:** 856866-72-3  
**Formal Name:** (5R)-3-[3-fluoro-4-[6-(2-methyl-2H-tetrazol-5-yl)-3-pyridinyl]phenyl]-5-(hydroxymethyl)-2-oxazolidinone  
**Synonyms:** DA 7157, TR-700  
**MF:** C<sub>17</sub>H<sub>15</sub>FN<sub>6</sub>O<sub>3</sub>  
**FW:** 370.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234, 302 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

Tedizolid is supplied as a crystalline solid. A stock solution may be made by dissolving the tedizolid in the solvent of choice, which should be purged with an inert gas. Tedizolid is slightly soluble in ethanol, DMSO, and dimethyl formamide.

### Description

Tedizolid is an oxazolidinone antibiotic and active metabolite of the antibiotic tedizolid phosphate (Item No. 23729).<sup>1</sup> 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 (ED<sub>50</sub> = 8.8 mg/kg).<sup>2</sup>

### References

1. Schaadt, 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, 01/04/2021

#### 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](http://WWW.CAYMANCHEM.COM)