

PRODUCT INFORMATION



PNU 100480

Item No. 28395

CAS Registry No.: 168828-58-8
Formal Name: N-[[[(5S)-3-[3-fluoro-4-(4-thiomorpholinyl)phenyl]-2-oxo-5-oxazolidinyl]methyl]-acetamide

Synonyms: Sutezolid, U 100480

MF: C₁₆H₂₀FN₃O₃S

FW: 353.4

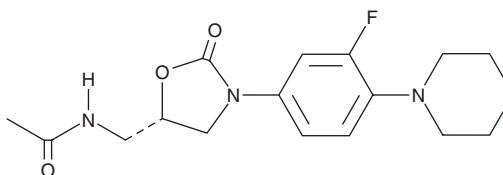
Purity: ≥98%

UV/Vis.: λ_{max}: 205, 259 nm

Supplied as: A crystalline 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

PNU 100480 is supplied as a crystalline solid. A stock solution may be made by dissolving the PNU 100480 in the solvent of choice, which should be purged with an inert gas. PNU 100480 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of PNU 100480 in these solvents is approximately 10 and 3 mg/ml, respectively.

PNU 100480 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PNU 100480 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. PNU 100480 has a solubility of approximately 0.14 mg/ml in a 1:6 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

PNU 100480 is an oxazolidinone antibiotic that is active against multidrug resistant tuberculosis.¹ It inhibits the growth of *M. tuberculosis* replicating and nonreplicating strains (MIC = ~400 ng/ml for both) and of clinical isolates that are susceptible or resistant to a combination of isoniazid (Item No. 20378), rifampin (rifampicin; Item No. 14423), ethambutol (Item No. 23713), and streptomycin (Item No. 21211; MICs = ≤0.0625-0.5 mg/L).^{1,2} It also inhibits the growth of *P. insidiosum* and *P. aphanidermatum* clinical isolates (MICs = 4-64 µg/ml).³ PNU 100480 (100 mg/kg per day) decreases the number of colony forming units (CFU) in the lung in a mouse model of systemic nonreplicating *M. tuberculosis* infection.²

References

- Alffenaar, J.W., van der Laan, T., Simons, S., et al. Susceptibility of clinical *Mycobacterium tuberculosis* isolates to a potentially less toxic derivate of linezolid, PNU-100480. *Antimicrob. Agents Chemother.* **55**(3), 1287-1289 (2011).
- Zhang, M., Sala, C., Dhar, N., et al. *In vitro* and *in vivo* activities of three oxazolidinones against nonreplicating *Mycobacterium tuberculosis*. *Antimicrob. Agents Chemother.* **58**(6), 3217-3223 (2014).
- Loreto, E.S., Tondolo, J.S.M., Oliveira, D.C., et al. *In vitro* activities of miltefosine and antibacterial agents from the macrolide, oxazolidinone, and pleuromutilin classes against *Pythium insidiosum* and *Pythium aphanidermatum*. *Antimicrob. Agents Chemother.* **62**(3), e01678-17 (2018).

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, 12/12/2022

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