

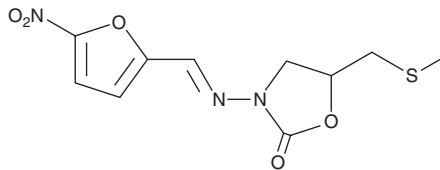
PRODUCT INFORMATION



Nifuratel

Item No. 40119

CAS Registry No.: 4936-47-4
Formal Name: 5-[(methylthio)methyl]-3-[[[(5-nitro-2-furanyl)methylene]amino]-2-oxazolidinone
Synonym: Methylmercadone
MF: C₁₀H₁₁N₃O₅S
FW: 285.3
Purity: ≥95%
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

Nifuratel is supplied as a solid. A stock solution may be made by dissolving the nifuratel in the solvent of choice, which should be purged with an inert gas. Nifuratel is soluble in methanol and DMSO.

Description

Nifuratel is a nitrofurantoin antibiotic and an antiparasitic.^{1,2} It is active against *G. vaginalis* and *A. vaginae* (MIC₉₀s = 0.25 and <0.125 µg/ml, respectively).¹ *In vivo*, nifuratel (50 mg/kg) reduces parasitic load in a mouse model of *L. donovani*-induced visceral leishmaniasis and a mouse model of *L. major*-induced cutaneous leishmaniasis.² Formulations containing nifuratel have been used in the treatment of vulvovaginal infections and urinary tract infections.

References

1. Togni, G., Battini, V., Bulgheroni, A., *et al.* *In vitro* activity of nifuratel on vaginal bacteria: Could it be a good candidate for the treatment of bacterial vaginosis? *Antimicrob. Agents Chemother.* **55**(5), 2490-2492 (2011).
2. Domínguez-Asenjo, B., Gutiérrez-Corbo, C., Álvarez-Bardón, M., *et al.* Ex vivo phenotypic screening of two small repurposing drug collections identifies nifuratel as a potential new treatment against visceral and cutaneous leishmaniasis. *ACS Infect. Dis.* **7**(8), 2390-2401 (2021).

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, 01/23/2024

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