

PRODUCT INFORMATION

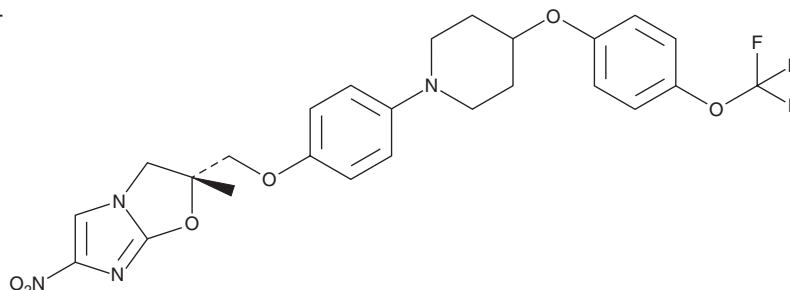


Delamanid

Item No. 29697

CAS Registry No.: 681492-22-8
Formal Name: (2R)-2,3-dihydro-2-methyl-6-nitro-2-[[4-[4-[4-(trifluoromethoxy)phenoxy]-1-piperidiny]phenoxy]methyl]-imidazo[2,1-b]oxazole

Synonym: OPC-67683
MF: C₂₅H₂₅F₃N₄O₆
FW: 534.5
Purity: ≥98%
UV/Vis.: λ_{max}: 220, 244, 316 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

Delamanid is supplied as a crystalline solid. A stock solution may be made by dissolving the delamanid in the solvent of choice, which should be purged with an inert gas. Delamanid is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of delamanid in these solvents is approximately 20 and 25 mg/ml, respectively. Delamanid is slightly soluble in ethanol.

Description

Delamanid is an antimycobacterial compound.¹ It is active against clinical isolates of *M. tuberculosis* that are susceptible (MIC_{90S} = 0.01194-0.01248 µg/ml) or resistant to rifampicin (Item No. 14423), isoniazid (Item No. 20378), ethambutol (Item No. 23713), and/or streptomycin (Item No. 21211; MIC_{90S} = 0.01221-0.01341 µg/ml). It inhibits the synthesis of methoxy- and keto-mycolic acid in the *M. bovis* strain BCG (IC_{50S} = 0.036 and 0.021 µg/ml, respectively). Delamanid (0.1, 0.3, and 1 µg/ml) inhibits the growth of the *M. tuberculosis* strain H37Rv in infected THP-1 cells. It decreases lung bacterial burden in a mouse model of *M. tuberculosis* infection in a dose-dependent manner.

Reference

1. Matsumoto, M., Hashizume, H., Tomishige, T., et al. OPC-67683, a nitro-dihydro-imidazooxazole derivative with promising action against tuberculosis in vitro and in mice. *PLoS Med.* **3**(11), e466 (2006).

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, 12/07/2022

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