

PRODUCT INFORMATION

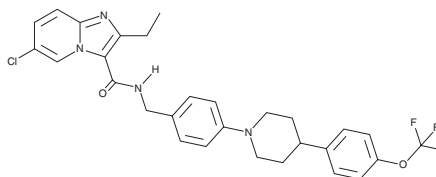


Q203

Item No. 37323

CAS Registry No.: 1334719-95-7
Formal Name: 6-chloro-2-ethyl-N-[[4-[4-(trifluoromethoxy)phenyl]-1-piperidinyl]phenyl]methyl]-imidazo[1,2-a]pyridine-3-carboxamide

Synonym: Telacebec
MF: C₂₉H₂₈ClF₃N₄O₂
FW: 557.0
Purity: ≥98%
UV/Vis.: λ_{max}: 257 nm
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

Q203 is supplied as a solid. A stock solution may be made by dissolving the Q203 in the solvent of choice, which should be purged with an inert gas. Q203 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of Q203 in DMSO and DMF is approximately 14 and 16 mg/ml, respectively. Q203 is slightly soluble in ethanol.

Description

Q203 is an imidazopyridine amide antibiotic.¹ It is active against *M. tuberculosis* and *M. ulcerans* (MIC₅₀s = 2.7 and 0.86 nM, respectively).^{1,2} Q203 binds to the oxidation site on mitochondrial complex III, also known as cytochrome bc₁ complex, and disrupts ATP synthesis in *M. tuberculosis* (IC₅₀ = 1.1 nM).¹ It reduces bacterial load, as well as the number of lung colony forming units (CFUs) and granulomatous foci in a mouse model of *M. tuberculosis* infection when administered at a dose of 10 mg/kg for three days.

References

1. Pethe, K., Bifani, P., Jang, J., *et al.* Discovery of Q203, a potent clinical candidate for the treatment of tuberculosis. *Nat. Med.* **19**(9), 1157-1160 (2013).
2. Scherr, N., Bieri, R., Thomas, S.S., *et al.* Targeting the *Mycobacterium ulcerans* cytochrome bc₁:aa₃ for the treatment of Buruli ulcer. *Nat. Commun.* **9**(1), 5370 (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, 03/01/2023

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