

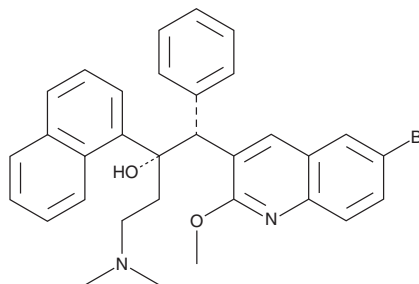
# PRODUCT INFORMATION



## TMC207

Item No. 20247

**CAS Registry No.:** 843663-66-1  
**Formal Name:** (αS,βR)-6-bromo-α-[2-(dimethylamino)ethyl]-2-methoxy-α-1-naphthalenyl-β-phenyl-3-quinolineethanol  
**Synonyms:** Bedaquiline, R207910  
**MF:** C<sub>32</sub>H<sub>31</sub>BrN<sub>2</sub>O<sub>2</sub>  
**FW:** 555.5  
**Purity:** ≥97%  
**UV/Vis.:** λ<sub>max</sub>: 226, 275, 285, 318, 333 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

TMC207 is supplied as a crystalline solid. A stock solution may be made by dissolving the TMC207 in the solvent of choice, which should be purged with an inert gas. TMC207 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of TMC207 in these solvents is approximately 0.2, 12.5, and 25 mg/ml, respectively.

TMC207 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TMC207 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TMC207 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

TMC207 is a diarylquinoline that selectively inhibits the proton pump of the mycobacterial ATP synthase.<sup>1</sup> It demonstrates potent activity against both drug-sensitive and drug-resistant *M. tuberculosis* and other mycobacterial species with MIC<sub>50</sub> values of ~0.03 µg/ml.<sup>1</sup>

### Reference

1. Matteelli, A., Carvalho, C. C., Dooley, K. E., *et al.* TMC207: The first compound of a new class of potent anti-tuberculosis drugs. *Future Microbiol.* 5(6), 849-858 (2010).

**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, 10/17/2022

### 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