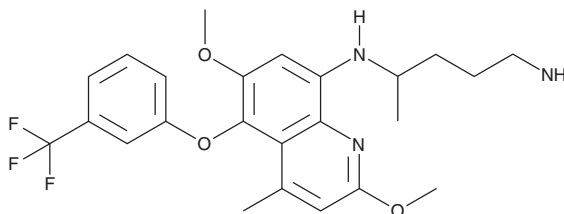


# PRODUCT INFORMATION



## Tafenoquine Item No. 30673

**CAS Registry No.:** 106635-80-7  
**Formal Name:** N<sup>4</sup>-[2,6-dimethoxy-4-methyl-5-[3-(trifluoromethyl)phenoxy]-8-quinolinyl]-1,4-pentanediamine  
**Synonym:** WR 238605  
**MF:** C<sub>24</sub>H<sub>28</sub>F<sub>3</sub>N<sub>3</sub>O<sub>3</sub>  
**FW:** 463.5  
**Purity:** ≥98%  
**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

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

### Description

Tafenoquine is an aminoquinoline antimalarial agent.<sup>1</sup> It is active against multidrug-resistant clinical isolates of *P. falciparum* (IC<sub>50</sub>s = 64.6-110.4 ng/ml). Tafenoquine (25 or 50 mg/kg) inhibits sporozoite transmission to mosquitos fed on *P. berghei*-infected mice.<sup>2</sup> It also suppresses parasitemia in a hamster model of *B. microti* infection when administered at doses ranging from 3.25 to 52 mg/kg.<sup>3</sup>

### References

1. Ohrt, C., Willingmyre, G.D., Lee, P., *et al.* Assessment of azithromycin in combination with other antimalarial drugs against *Plasmodium falciparum* in vitro. *Antimicrob. Agents Chemother.* **46**(8), 2518-2524 (2002).
2. Coleman, R.E. Sporontocidal activity of the antimalarial WR-238605 against *Plasmodium berghei* ANKA in *Anopheles stephensi*. *Am. J. Trop. Med. Hyg.* **42**(3), 196-205 (1990).
3. Marley, S.E., Eberhard, M.L., Steurer, F.J., *et al.* Evaluation of selected antiprotozoal drugs in the *Babesia microti*-hamster model. *Antimicrob. Agents Chemother.* **41**(1), 91-94 (1997).

#### 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](http://WWW.CAYMANCHEM.COM)