

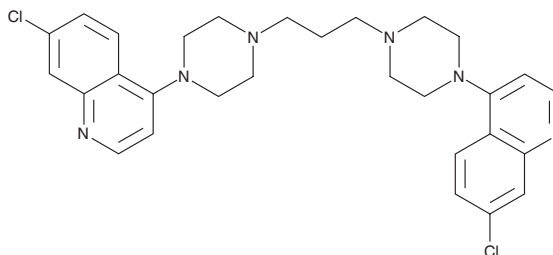
PRODUCT INFORMATION



Piperaquine

Item No. 20799

CAS Registry No.: 4085-31-8
Formal Name: 4,4'-(1,3-propanediyl-di-4,1-piperazinediyl)bis[7-chloro-quinoline]
Synonym: Piperaquinoline
MF: C₂₉H₃₂Cl₂N₆
FW: 535.5
Purity: ≥98%
UV/Vis.: λ_{max}: 225, 350 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

Piperaquine is supplied as a crystalline solid. A stock solution may be made by dissolving the piperaquine in the solvent of choice. Piperaquine is soluble in the organic solvent DMSO, which should be purged with an inert gas, at a concentration of approximately 0.2 mg/ml.

Description

Piperaquine is a *bis*-quinoline compound that inhibits the *ex vivo* growth of *P. falciparum* isolates from malaria-infected patients with IC₅₀s ranging from 11.8-217.3 nM.¹ Formulations containing piperaquine have been used in combination with artemisinin (Item No. 11816) for the treatment of malaria in HIV-1-positive populations.²⁻⁴

References

1. Akoachere, M., Buchholz, K., Fischer, E., *et al.* In vitro assessment of methylene blue on chloroquine-sensitive and -resistant *Plasmodium falciparum* strains reveals synergistic action with artemisinins. *Antimicrob. Agents Chemother.* **49(11)**, 4592-4597 (2005).
2. Askling, H.H., Bruneel, F., Burchard, G., *et al.* Management of imported malaria in Europe. *Malar. J.* **11**, (2012).
3. Henrich, P.P., O'Brien, C., Sáenz, F.E., *et al.* Evidence for pyronaridine as a highly effective partner drug for treatment of artemisinin-resistant malaria in a rodent model. *Antimicrob. Agents Chemother.* **58(1)**, 183-195 (2014).
4. Pascual, A., Parola, P., Benoit-Vical, F., *et al.* *Ex vivo* activity of the ACT new components pyronaridine and piperaquine in comparison with conventional ACT drugs against isolates of *Plasmodium falciparum*. *Malar. J.* **11**, (2012).

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, 12/19/2022

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