

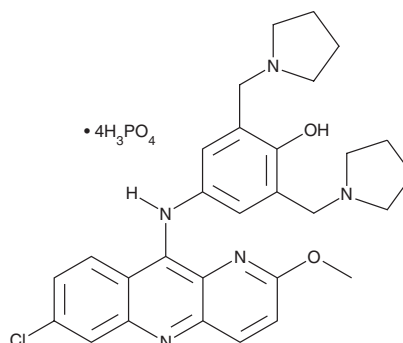
PRODUCT INFORMATION



Pyronaridine (phosphate)

Item No. 16139

CAS Registry No.: 76748-86-2
Formal Name: 4-[(7-chloro-2-methoxybenzo[b]-1,5-naphthyridin-10-yl)amino]-2,6-bis(1-pyrrolidinylmethyl)-phenol, tetraphosphate
Synonym: PND
MF: C₂₉H₃₂ClN₅O₂ • 4H₃PO₄
FW: 910.0
Purity: ≥98%
UV/Vis.: λ_{max}: 239, 278, 359, 425 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

Pyronaridine (phosphate) is supplied as a crystalline solid. Aqueous solutions of pyronaridine (phosphate) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of pyronaridine (phosphate) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Pyronaridine is an antimalarial agent.¹ It is active against six *P. falciparum* drug-sensitive or -resistant strains (IC₅₀s = 1.53-3.94 nM). Pyronaridine (3, 6, and 8 mg/kg) in combination with artesunate (Item No. 11817) reduces parasitemia in *P. chabaudi*-infected mice. It inhibits the replication of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) in infected A459 cells overexpressing angiotensin-converting enzyme 2 (ACE2; IC₅₀ = 198 nM).² Pyronaridine is also cytotoxic against a panel of human cancer cells lines, including breast, ovarian, and lung cancer cells, with 50% cytotoxic concentration values (CC₅₀s) ranging from 1.6 to 9.4 μM.³ Formulations containing pyronaridine have been used in the treatment of malaria.

References

1. Vivas, L., Rattray, L., Stewart, L., et al. Anti-malarial efficacy of pyronaridine and artesunate in combination *in vitro* and *in vivo*. *Acta Trop.* **105**(3), 222-228 (2008).
2. Puhl, A.C., Fritch, E.J., Lane, T.R., et al. Repurposing the ebola and marburg virus inhibitors tilorone, quinacrine, and pyronaridine: *In vitro* activity against SARS-CoV-2 and potential mechanisms. *ACS Omega* **6**(11), 7454-7468 (2021).
3. Villanueva, P.J., Martinez, A., Baca, S.T., et al. Pyronaridine exerts potent cytotoxicity on human breast and hematological cancer cells through induction of apoptosis. *PLoS One* **13**(11), e0206467 (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, 10/07/2022

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