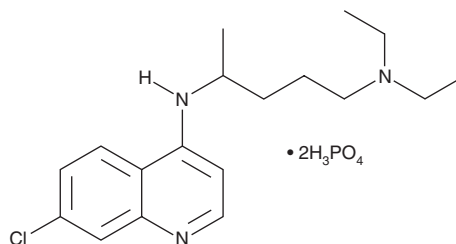


PRODUCT INFORMATION

Chloroquine (phosphate)

Item No. 30708

CAS Registry No.: 50-63-5
Formal Name: N⁴-(7-chloro-4-quinoliny)-N¹,N¹-diethyl-1,4-pentanediamine, diphosphate
Synonyms: DL-Chloroquine, NSC 14050
MF: C₁₈H₂₆ClN₃ • 2H₃PO₄
FW: 515.9
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Description

Chloroquine (phosphate) (Item No. 30708) is an analytical reference material categorized as an antimalarial, anti-inflammatory, and antiviral agent.¹⁻³ Chloroquine prevents infection by severe acute respiratory coronavirus 2 (SARS-CoV-2) *in vitro*.³ Counterfeit chloroquine has been seized by law enforcement during the COVID-19 pandemic.⁴ Formulations containing chloroquine have been associated with fatal overdoses.^{5,6} This product is intended for use in analytical research applications. This product is also available as a general research tool (Item No. 14194).

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards for reference materials.

References

1. Sidhu, A.B.S., Verdier-Pinard, D., Fidock, D.A. Chloroquine resistance in *Plasmodium falciparum* malaria parasites conferred by *pfcr* mutations. *Science*. **298**(5591), 210-213 (2002).
2. Plantone, D. and Koudriavtseva, T. Current and future use of chloroquine and hydroxychloroquine in infectious, immune, neoplastic, and neurological diseases: A mini-review. *Clin. Drug Investig.* **38**(8), 653-671 (2018).
3. Wang, M., Cao, R., Zhang, L., *et al.* Remdesivir and chloroquine effectively inhibit the recently emerged novel coronavirus (2019-nCoV) *in vitro*. *Cell Res.* **30**(3), 269-271 (2020).
4. Schneider, M. and Nam, N.H.T. Africa and counterfeit pharmaceuticals in the times of COVID-19. *J. Intellect. Prop. Law* **15**(6), 417-418 (2020).
5. Queen, H.F., Tapfumaneyi, C., Lewis, R.J. The rising incidence of serious chloroquine overdose in Harare, Zimbabwe: Emergency department surveillance in the developing world. *Trop. Doct.* **29**(3), 139-141 (1999).
6. Keller, T., Schneider, A., Lamprecht, R., *et al.* Fatal chloroquine intoxication. *Forensic Sci. Int.* **96**(1), 21-28 (1998).

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, 11/02/2020

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