

# PRODUCT INFORMATION



## Hydroxychloroquine (sulfate)

Item No. 30709

CAS Registry No.: 747-36-4

Formal Name: 2-[[4-[(7-chloro-4-quinolinyl)amino]pentyl]Cl-ethylamino]-ethanol, monosulfate

Synonyms: HCQ, NSC 4375

MF:  $C_{18}H_{26}ClN_3O \cdot H_2SO_4$

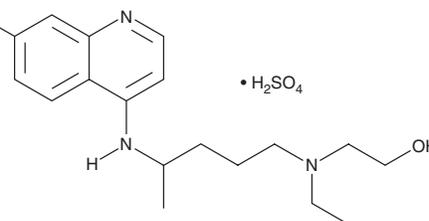
FW: 434.0

Purity:  $\geq 98\%$

Supplied as: A neat solid

Storage:  $-20^{\circ}C$

Stability:  $\geq 1$  year



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

### Description

Hydroxychloroquine (sulfate) (Item No. 30709) is an analytical reference material categorized as an antimalarial, anti-inflammatory, and antiviral agent.<sup>1-4</sup> Hydroxychloroquine reduces viral titers of severe acute respiratory syndrome coronavirus 2 (SARS-CoV-2) *in vitro*.<sup>4</sup> Formulations containing hydroxychloroquine have been associated with fatal overdoses.<sup>5</sup> This product is intended for analytical research applications. This product is also available as a general research tool (Item No. 17911).

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. Delves, M., Plouffe, D., Scheurer, C., *et al.* The activities of current antimalarial drugs on the life cycle stages of *Plasmodium*: A comparative study with human and rodent parasites. *PLoS Med.* **9(2)**, e1001169 (2012).
2. Mauthe, M., Orhon, I., Rocchi, C., *et al.* Chloroquine inhibits autophagic flux by decreasing autophagosome-lysosome fusion. *Autophagy* **14(8)**, 1435-1455 (2018).
3. da Silva, J.C., Mariz, H.A., da Rocha, L.F., Jr., *et al.* Hydroxychloroquine decreases Th17-related cytokines in systemic lupus erythematosus and rheumatoid arthritis patients. *Clinics (Sao Paulo)* **68(6)**, 766-771 (2013).
4. Maisonnasse, P., Guedj, J., Contreras, V., *et al.* Hydroxychloroquine use against SARS-CoV-2 infection in non-human primates. *Nature* **585(7826)**, 584-587 (2020).
5. Marquardt, K. and Albertson, T.E. Treatment of hydroxychloroquine overdose. *Am. J. Emerg. Med.* **19(5)**, 420-424 (2001).

#### 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, 11/03/2020

#### 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