

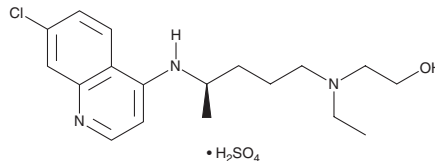
PRODUCT INFORMATION



(R)-Hydroxychloroquine (sulfate)

Item No. 33442

CAS Registry No.: 2488706-20-1
Formal Name: 2-[[[(4R)-4-[(7-chloro-4-quinolinyl)amino]pentyl]ethylamino]-ethanol, monosulfate
Synonym: (R)-HCQ
MF: C₁₈H₂₆ClN₃O • H₂SO₄
FW: 434.0
Purity: ≥95%
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

(R)-Hydroxychloroquine (sulfate) is supplied as a solid. A stock solution may be made by dissolving the (R)-hydroxychloroquine (sulfate) in the solvent of choice, which should be purged with an inert gas. (R)-Hydroxychloroquine (sulfate) is soluble in methanol and DMSO. It is also soluble in water.

Description

(R)-Hydroxychloroquine is an isomer of hydroxychloroquine (Item No. 17911).¹ (R)-Hydroxychloroquine (2.5-20 mg/ml) reduces the rate of insulin degradation in isolated liver homogenates from non-diabetic and diabetic rats. Formulations containing hydroxychloroquine have been used in the prevention or treatment of malaria, as well as in the treatment of rheumatoid arthritis and systemic lupus erythematosus.

Reference

1. Emami, J., Pasutto, F.M., Mercer, J.R., *et al.* Inhibition of insulin metabolism by hydroxychloroquine and its enantiomers in cytosolic fraction of liver homogenates from healthy and diabetic rats. *Life Sci.* **64(5)**, 325-335 (1999).

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, 11/15/2022

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