

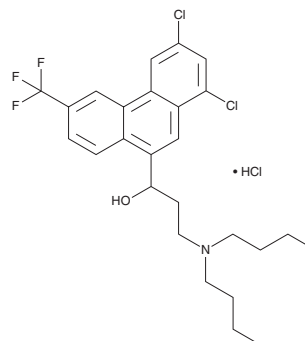
PRODUCT INFORMATION



Halofantrine (hydrochloride)

Item No. 30962

CAS Registry No.: 36167-63-2
Formal Name: 1,3-dichloro- α -[2-(dibutylamino)ethyl]-6-(trifluoromethyl)-9-phenanthrenemethanol, monohydrochloride
Synonyms: (\pm)-Halofantrine, SKF 102886, WR 171669
MF: $C_{26}H_{30}Cl_2F_3NO \cdot HCl$
FW: 536.9
Purity: $\geq 98\%$
UV/Vis.: λ_{max} : 216, 232, 259 nm
Supplied as: A crystalline solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Halofantrine (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the halofantrine (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Halofantrine (hydrochloride) is soluble in DMSO and methanol.

Description

Halofantrine is an antimalarial agent.¹ It is active against chloroquine-sensitive and chloroquine-resistant strains of *P. falciparum* (IC_{50} s = 1.5-2.5 and 1.3-3.9 $\mu g/L$, respectively). Halofantrine reduces parasitemia in a mouse model of *P. berghei* infection with a 50% curative dose (CD_{50}) value of 15 mg/kg. It also reduces parasitemia in an *Aotus* monkey model of *P. falciparum* infection (CD_{50} = 58.3 mg/kg). Formulations containing halofantrine have been used in the treatment of malaria.

Reference

1. Bryson, H. and Goa, K.L. Halofantrine. A review of its antimalarial activity, pharmacokinetic properties and therapeutic potential. *Drugs* **43**(2), 236-258 (1992).

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, 12/14/2022

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