

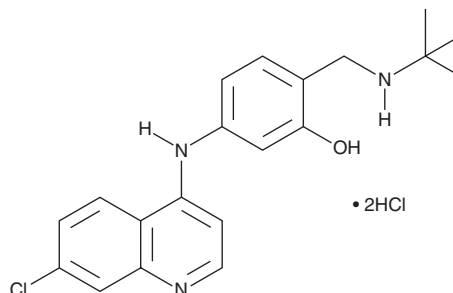
PRODUCT INFORMATION



GSK369796 (hydrochloride)

Item No. 36061

CAS Registry No.: 1010411-21-8
Formal Name: 5-[(7-chloro-4-quinolinyl)amino]-2-[[[(1,1-dimethylethyl)amino]methyl]-phenol, dihydrochloride
Synonym: N-tert-butyl Isoquine
MF: C₂₀H₂₂ClN₃O • 2HCl
FW: 428.8
Purity: ≥95%
UV/Vis.: λ_{max}: 224, 256, 344 nm
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

GSK369796 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the GSK369796 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. GSK369796 (hydrochloride) is soluble in organic solvents such as DMSO. It is also soluble in water. We do not recommend storing the aqueous solution for more than one day.

Description

GSK369796 is an antimalarial agent.¹ It is active against the chloroquine-sensitive *P. falciparum* strains 3D7 and HB3 (IC₅₀s = 11.2 and 12.6 nM, respectively) and the chloroquine-resistant strain K1 (IC₅₀ = 13.2 nM). It reduces parasitemia in mice infected with *P. berghei* (ED₅₀ = 2.8 mg/kg).

Reference

1. O'Neill, P.M., Park, B.K., Shone, A.E., *et al.* Candidate selection and preclinical evaluation of N-tert-butyl isoquine (GSK369796), an affordable and effective 4-aminoquinoline antimalarial for the 21st century. *J. Med. Chem.* **52**(5), 1408-1415 (2009).

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, 12/19/2022

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