

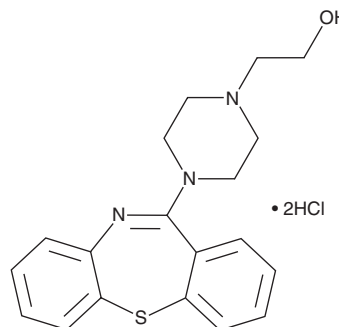
PRODUCT INFORMATION



Desethoxy Quetiapine (hydrochloride)

Item No. 35410

CAS Registry No.: 329218-14-6
Formal Name: 4-dibenzo[b,f][1,4]thiazepin-11-yl-1-piperazineethanol, dihydrochloride
Synonym: O-Desalkylquetiapine
MF: C₁₉H₂₁N₃OS • 2HCl
FW: 412.4
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

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

Description

Desethoxy quetiapine is an active metabolite of the atypical antipsychotic quetiapine (Item No. 14152).¹ It is formed from quetiapine primarily by the cytochrome P450 (CYP) isoform CYP3A5 and binds to dopamine D₂ receptors (IC₅₀ = 1,330 nM).^{1,2} Desethoxy quetiapine (20 and 40 mg/kg) reduces climbing induced by the dopamine receptor agonist apomorphine in mice. It also improves apomorphine-induced swimming deficits in mice.¹

References

1. Warawa, E.J., Migler, B.M., Ohnmacht, C.J., *et al.* Behavioral approach to nondyskinetic dopamine antagonists: Identification of seroquel. *J. Med. Chem.* **44**(3), 372-389 (2001).
2. Bakken, G.V., Rudberg, I., Christensen, H., *et al.* Metabolism of quetiapine by CYP3A4 and CYP3A5 in presence or absence of cytochrome B₅. *Drug Metab. Dispos.* **37**(2), 254-258 (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, 10/24/2022

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