

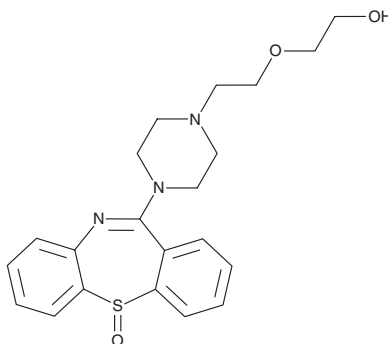
PRODUCT INFORMATION



Quetiapine Sulfoxide

Item No. 42035

CAS Registry No.: 329216-63-9
Formal Name: 11-(4-(2-(2-hydroxyethoxy)ethyl) piperazin-1-yl)dibenzo[b,f][1,4]thiazepine 5-oxide
MF: $C_{21}H_{25}N_3O_3S$
FW: 399.5
Purity: $\geq 95\%$
Supplied as: A 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

Quetiapine sulfoxide is supplied as a solid. A stock solution may be made by dissolving the quetiapine sulfoxide in the solvent of choice, which should be purged with an inert gas. Quetiapine sulfoxide is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of quetiapine sulfoxide can be prepared by directly dissolving the solid in aqueous buffers. Quetiapine sulfoxide is slightly soluble (0.1-1 mg/ml) in PBS (pH 7.2). We do not recommend storing the aqueous solution for more than one day.

Description

Quetiapine sulfoxide is an inactive metabolite of the atypical antipsychotic quetiapine (Item No. 14152).¹ Quetiapine sulfoxide is formed from quetiapine via the cytochrome P450 (CYP) isoform CYP3A.

Reference

1. 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, 10/17/2024

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