

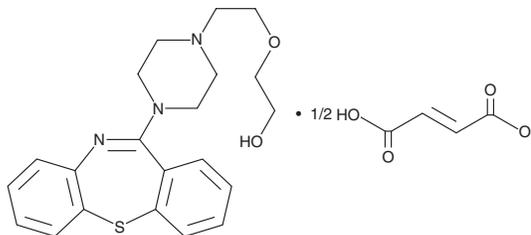
PRODUCT INFORMATION



Quetiapine (hemifumarate)

Item No. 22603

CAS Registry No.: 111974-72-2
Formal Name: 2-[2-(4-dibenzo[b,f][1,4]thiazepin-11-yl-1-piperazinyl)ethoxy]ethanol, 2E-butenedioate (2:1)
Synonym: ICI 204636
MF: C₂₁H₂₅N₃O₂S • 1/2C₄H₄O₄
FW: 441.5
Purity: ≥98%
Supplied as: A neat solid
Storage: -20°C
Stability: ≥1 year



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

Description

Quetiapine (hemifumarate) (Item No. 22603) is an analytical reference material categorized as an atypical antipsychotic.¹ It is misused and abused more commonly than other atypical antipsychotics and is associated with drug-seeking behavior.^{2,3} This product is intended for use in analytical forensic applications. This product is also available as a general research tool (Item No. 14152).

This product is qualified as a Reference Material that has been manufactured and tested to ISO/IEC 17025 and ISO 17034 international standards.

References

1. Tamminga, C.A. and Lahti, A.C. The new generation of antipsychotic drugs. *Int. Clin. Psychopharmacol.* **11(suppl 2)**, 73-76 (1996).
2. Klein, L., Bangh, S., and Cole, J.B. Intentional recreational abuse of quetiapine compared to other second-generation antipsychotics. *West. J. Emerg. Med.* **18(2)**, 243-250 (2017).
3. Kim, S., Lee, G., Kim, E., *et al.* Quetiapine misuse and abuse: Is it an atypical paradigm of drug seeking behavior? *J. Res. Pharm. Pract.* **6(1)**, 12-15 (2017).

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, 07/01/2019

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